

Operating Manual

Fluent®



제목:	Fluent 작동 설명서			부품 번호:	30256654.01
ID:	399706, ko, V2.7			번역 원본:	해당사항 없음
버전:	개정:	발행:	문서 이력:		·
1	0	2017년 8월 30 일	제1판		
1	1	2017년 9월 18 일	2.3 적용 시 위험		
1	2	2017년 11월 07일	1.3 사용 목적		
1	3	2018년 03월 05일	1.3 사용 목적		
1	4	2018년 04월 16일	3.3 치수 및 중량		
1	5	2018년 07월 20일	4.6.2 FCA 그리퍼		
1	6	2019년 02월 26일	2.2 일반 안전 정보		
1	7	2020년 10월 22일	1.8.3, 2.5, 2.8. 7.5.4, 8.2.3, 8.3 정보가 추가되었 가 섹션 4.6.3, 6 8.2.3, 8.3.14에 7.5.24가 추가되 4.2.2, 4.6, 6.3.9 8.3.13-20이 업대	1, 4.6, 4.6.3, 6.6 3.13-15 섹션에 M 습니다. 튜브 로테 .5.3, 7.4.1, 7.4.2 추가되었습니다. 었습니다. 섹션 2. 5, 7.1, 7.4, 7.5.1 베이트되었습니다.), 7.4.1, 7.4.2, lix & Pierce 관련 이터에 대한 정보 2, 7.5.5, 7.5.6, 섹션 4.6.4 및 5, 2.9, 2.10, 6, 8.2.4,
1	8	2020년 12월 01일	섹션 2.2 및 2.80	비 업데이트되었습	니다.
1	9	2021년 02월 15일	섹션 2.7.2, 7.1, 8.2.3이 업데이트 7.5.25가 업데이.	7.2, 7.4.2, 7.5.1 되었습니다. 섹션 트되었습니다.	15, 7.5.24, 2.7.5, 4.6.4,
1	10	2021년 11월 25일	섹션 3.7, 6.5.10	업데이트되었습	니다.
2	0	2022년 03월 02일	섹션 3.1이 업데() 트되었습니다.	
2	1	2022년 04월 14일	섹션 5.2.5, 6.60 업데이트되었습니	ㅣ추가되었습니다 ㅣ다.	. 섹션 6.5.1이

© 2025, Tecan Trading AG, Switzerland, all rights reserved.

Information contained in this document is subject to change without notice.

2	2	2022년 11월 16일	섹션 4.6.5가 추가되었습니다. 섹션 2.2 및 3.4가 업데이트되었습니다.
2	3	2023년 03월 15일	섹션 1.8.2, 2.4, 3.4, 7.4.2가 업데이트되었습니 다.
2	4	2023년 03월 30일	섹션 2.4, 2.8.1.1, 3.3, 4.3.2, 4.3.3, 4.3.4, 4.3.5.3, 7.4.2, 7.4.4가 업데이트되었습니다.
2	5	2024년 01월 16일	섹션 1.9, 2.2, 2.7.1, 2.8.1, 6.2, 7, 7.2.1, 7.2.2, 7.4.2, 7.5.14, 8.2.2, 8.3.6이 업데이트되었습니 다.
2	6	2024년 06월 06일	1, 1.8.4, 2.2, 2.4, 3.7, 4.6, 6.1 섹션이 업데이트 되고 4.6.6 섹션이 추가됨
2	7	2024년 12월 05일	섹션 1.8.4, 3.7, 4.6, 6.5, 7.4.1.2, 7.5.18, 10.1 이 업데이트되었습니다., 2.12가 추가되었습니다.

© 2025, Tecan Trading AG, Switzerland, all rights reserved.

Information contained in this document is subject to change without notice.

© 2025, Tecan Trading AG, Switzerland, all rights reserved.

Information contained in this document is subject to change without notice.

•TECAN•

목차

1	설명시	ㅓ 정보	9
	1.1	본 설명서의 범위	9
	1.2	제조업체	9
	1.3	용도	9
	1.4	사용 영역/적용 영역	9
	1.5	부적절한 사용	10
	1.6	보증	10
	1.7	상표	10
	1.8	참조 문서	10
	1.9	법률 및 표준 준수	11
	1.10	문서 규칙	11
n	아저		10
2	인전	아저메시지 그치	10
	2.1	안전 메시지 규칙	12
	2.2	일만 안산 경모	14
	2.3	카메라 개인정보 처리방침	15
	2.4	애플리케이션 위험	15
	2.5	운영 회사	21
	2.6	검사법 및 프로세스 검증	22
	2.7	사용자 자격 평가	22
	2.8	안전 요소	24
	2.9	제품 안전 표지판	30
	2.10	레이저 방사선	35
	2.11	오염 제거 고지	36
	2.12	사고 보고	36
3	기술	데이터	37
	3.1	타입 플레이트	37
	3.2	일련번호 레이블	38
	3.3	치수 및 중량	39



	3.4	전원 공급	40
	3.5	데이터 및 전원 연결	41
	3.6	환경 조건	41
	3.7	방출 및 내성	42
4	기능	설명	44
	4.1	개요	44
	4.2	데크	44
	4.3	로봇 암	47
	4.4	액체 시스템(Liquid FCA)	53
	4.5	세척 시스템(MCA 384)	53
	4.6	옵션 및 장치	54
5	제어	요소	65
	5.1	작동 요소	65
	5.2	사용자 인터페이스	66
	5.3	오류 신호 및 기기 상태	70
	5.4	Fluent ID 상태 LED	72
6	작동		73
	6.1	이 장의 안전 지침	73
	6.2	작동 모드	74
	6.3	작동 준비	75
	6.4	검사법 시작 전	83
	6.5	검사법 실행하기	87
	6.6	DeckCheck 작업	99
	6.7	검사법 복구	101
	6.8	기기 전원 끄기	102
7	시스텪	볰 지원	104
	7.1	오염 제거	104
	7.2	세정제	105
	7.3	시스템 관리 모드	106



	7.4	시스템 관리 표	107
	7.5	시스템 관리 활동	115
8	문제	해결	144
	8.1	이 장의 안전 지침	144
	8.2	문제 해결 표	144
	8.3	문제 해결 활동	155
9	포장,	포장 풀기, 운반, 보관 및 폐기	185
9	포장, 9.1	포장 풀기, 운반, 보관 및 폐기 포장 라벨	185 185
9	포장, 9.1 9.2	포장 풀기, 운반, 보관 및 폐기 포장 라벨 폐기	185 185 186
9	포장, 9.1 9.2	포장 풀기, 운반, 보관 및 폐기 포장 라벨 폐기	185 185 186
9 10	포장, 9.1 9.2) 고객	포장 풀기, 운반, 보관 및 폐기 포장 라벨 폐기 지원	185185186188
9 10	포장, 9.1 9.2) 고객 10.1	포장 풀기, 운반, 보관 및 폐기 포장 라벨 폐기 지원 연락처	185185186188188
9 10	포장, 9.1 9.2) 고객 10.1	포장 풀기, 운반, 보관 및 폐기 포장 라벨 폐기 지원 연락처	 185 185 186 188 188





1 설명서 정보

이 Operating Manual(작동 설명서)은 에 대해 설명하고 안전한 작동 및 양호한 작 동 순서 유지에 필요한 모든 정보를 제공합니다. 이 설명서는 에 대한 작업을 수행 하기 전과 기기를 사용하기 전에 주의 깊게 읽어야 합니다. 이 장에서는 본 설명서의 목적에 대해 설명하고 참조되는 제품을 명시합니다. 더불 어 기호와 규약 사용과 추가 일반 정보를 설명합니다. 이 설명서는 기기 자체에 적용됩니다. 하위 모듈에 대한 중요 정보는 해당 모듈의 특정 설명서를 참조하십시오.



이 Operating Manual(작동 설명서)에는 소프트웨어 설명이 포함되어 있지 않습니 다 소프트웨어에 대한 자세한 내용은 해당 소프트웨어 설명서를 참조하십시오. 참 조 문서 섹션을 참조하십시오.

1.1 본 설명서의 범위

본 설명서는 다음에 적용됩니다.

- Fluent 480(부품 번호 30042011)
- Fluent 780(부품 번호 30042021)
- Fluent 1080(부품 번호 30042031)

1.2 제조업체

제조업체 주소



Tecan Schweiz AG Seestrasse 103 CH-8708 Männedorf Switzerland

1.3 용도

Fluent은 일반적인 실험실 사용을 위한 완전 자동화된 실험실 액체 취급 플랫폼입 니다. 이 제품은 특정 검사 절차에서 실험 기구의 피펫팅, 액체 취급 및 로봇 조작 과 같은 일상적인 실험실 작업에 사용됩니다.



Fluent GX Assurance Software는 이 용도로 사용해야 합니다.

Fluent에서 사용할 수 있는 Tecan 및 타사 장치의 특정 옵션은 연구용으로만 사용 할 수 있습니다(RUO).

연구용 전용 옵션 또는 장치가 Fluent에 포함되어 있거나 Fluent GX Assurance Software가 설치되지 않은 경우 사용 목적은 다음과 같이 변경됩니다.

Fluent은(는) 연구 및 산업 분야에 사용할 수 있는 완전 자동화된 실험실 액체 취급 플랫폼입니다. 이 제품은 특정 검사 절차(예: 세포 기반 검사, 생화학 검사 및 화합 물 관리)에서 실험 기구의 피펫팅, 액체 취급 및 로봇 조작과 같은 일상적인 실험실 작업에 사용됩니다. 이 기기는 임상 진단 절차에 사용하도록 고안된 것이 아닙니 다.

1.4 사용 영역/적용 영역

Fluent은(는) 사용 목적에 따라 다양한 실험실 환경에서 사용할 수 있습니다.



각 환경에서 실험실의 적용 워크플로 또는 검사법에 사용되는 특정 액체 및 실험 기구와 함께 Fluent의 검증은 개별 실험실이 맡아 수행합니다.

1.5 부적절한 사용

부적절한 사용은 Fluent 안전 개념을 저해할 수 있습니다.

- Tecan에서 승인하지 않은 옵션 또는 구성 요소와 함께 Fluent 제품을 사용해서 는 안 됩니다.
- Fluent 제품은 방폭 장치가 아니므로 폭발의 위험이 있는 곳에 설치해서는 안 됩니다.
- Fluent은(는) 기능적 안전 장치가 없는 경우 사용해서는 안 됩니다.

1.6 보증

Tecan이 승인하지 않은 옵션 또는 구성 요소와 함께 Fluent 제품을 사용해서는 안 됩니다.

승인되지 않은 옵션을 사용할 경우 Fluent의 안전 개념이 손상될 수 있습니다.

승인되지 않은 구성 요소를 사용할 경우 NRTL 인증 및 EC 지침 등에 따라 국가 및 국제 표준에 대한 안전 및 규정 준수가 무효화됩니다.

1.7 상표

이 설명서에 언급된 제품 이름은 등록 상표 또는 미등록 상표의 여부와 관계 없이 식별 목적으로만 재생산되며 해당 소유자의 독점 재산입니다. 단순성을 위해 [®] 및 ™과 같은 상표 기호는 설명서에서 반복되지 않습니다.

1.8 참조 문서

이 섹션에서는 Fluent 사용 시 필요하거나 유용한 문서 목록을 제공합니다.

아래에 나열된 문서 ID는 루트 번호입니다. 따라서 문서의 언어, 문서 버전 또는 매 체(데이터 저장 매체, 하드카피, 다운로드 가능한 파일 등)에 대한 정보는 포함되지 않습니다.



주문 구성에 따라 옵션 장비의 작동 설명서도 해당됩니다.

해당 문서의 범위를 확인하여 올바른 버전을 소유하고 있는지 확인하십시오.

문서 ID는 주문 정보를 참조하지 않습니다. 주문 시 바인더, CD 케이스 등에 있는 번호를 참조하십시오

1.8.1 기기 설명서

- Fluent[®] 작동 설명서(문서 ID 399706)
- Fluent[®] 참조 설명서(문서 ID 399937)

1.8.2 소프트웨어 설명서

- Tecan 샘플 추적 애드온 소프트웨어 설명서(문서 ID 393933)
- FluentControl 애플리케이션 소프트웨어 설명서(문서 ID 399935)
- Introspect 애플리케이션 소프트웨어 설명서(문서 ID 400733)
- MissionControl 애플리케이션 소프트웨어 설명서(문서 ID 401940)



1.8.3 QC 키트 설명서

- QC 키트 사용 설명서(문서 ID 39709)
- QC 키트 애플리케이션 소프트웨어 설명서(문서 ID 397070)

1.8.4 기타 참조 문서

- Fluent[®] Carousel 작동 설명서(문서 ID 398350)
- HEPA 후드(문서 ID Caron 70072)
- Frida Reader[™] 사용 설명서(문서 ID 401882)
- Te-Shake™ 작동 설명서(문서 ID 391496)
- Te-VacS™ 작동 설명서(문서 ID 391236)
- Fluent[®] Stacker 작동 설명서(문서 ID 398658)
- MIO2 작동 설명서(문서 ID 394934)
- Resolvex i300 작동 설명서(문서 ID 402756)

1.9 법률 및 표준 준수

다음 선언 및 인증이 Fluent에 적용됩니다.

- 해당하는 EU 지침(CE 마크)에 대한 EC 적합성 선언
- Nationally Recognized Testing Laboratory(NRTL) 인증
- (IECEE) CB Scheme 인증(CB 마크)

표시에 대한 자세한 내용은 "타입 플레이트" [37] 섹션을 참조하십시오.

1.10 문서 규칙

상호 참조	상호 참조는 다음과 같이 표시됩니다(예:
	"안전" [▶ 12] 섹션 참조
	• "안전"은 해당 섹션 헤더를 나타냅니다
	• 페이지 번호는 대괄호로 제공됩니다
필수 조건	다음과 같은 전제 조건이 표시됩니다. 예:
	✔ "일반 안전 정보"를 읽었습니다.
팁	추가 팁은 다음과 같습니다(예:
A	안전 규칙과 기호에 대한 정보는 "안전" [▶ 12] 장을 참조하십시오.

그림에 Fluent 제품과 관련이 없는 구성 요소 버전이 표시될 수 있습니다.

그림



2 안전

이 장에서는 Fluent의 안전 개념을 설명하고, 올바른 동작에 대한 일반적인 규칙 및 Fluent 사용과 관련된 위험에 대한 경고를 제공합니다.

2.1 안전 메시지 규칙

2.1.1 신호 단어

표 1: 신호 단어

신호 단어	의미
<u>▲</u> 위험	예방하지 않을 경우 사망 또는 심각한 부상을 초래할 위험한 상황을 나타냅니다.
▲경고	예방하지 않을 경우 사망 또는 심각한 부상을 초래할 수 있는 위험한 상황을 나타냅니다.
<u> </u>	예방하지 않을 경우 경미하거나 중간 정도의 부 상을 초래할 수 있는 위험한 상황을 나타냅니 다.
공지사항	위험과 관련이 없지만, 예방하지 않을 경우 장 비의 손상 또는 오작동을 초래하거나 부정확한 프로세스 결과를 초래할 수 있는 상황을 나타냅 니다.



2.1.2 안전 기호





2.2 일반 안전 정보

▲ 경고

Fluent 제품은 현재 최첨단 기술 및 인정된 기술 안전 규정에 따라 설계 및 제작됩니다. 그럼에도 불구하고 적절한 주의와 관심 없이 Fluent 제품을 사용할 경우 사용자, 재산 및 환경에 대한 위험이 발생할 수 있습니다.

모든 사용자 및 직원의 안전은 이러한 안전 지침의 엄격한 준수와 본 설명서에 제 공된 안전 관련 경고에 대한 인식에 달려 있습니다.

- 다음 일반 안전 정보에 각별히 주의하십시오.
- 본 설명서는 여기에 설명된 작업을 수행하는 모든 사람들이 항상 사용할 수 있어야 합니다.
- 항상 장비와 함께 제공된 전원 케이블을 사용하십시오.
- 전원 케이블을 다른 제품과 함께 사용하지 마십시오.
- Fluent 제품과 관련된 위험 물질의 취급뿐 아니라 사용 또는 적용에 관한 지역, 주 및 연방 법과 같은 법규를 엄격히 준수해야 합니다.
- 운영 회사는 회사 절차 및 현지 법률 요건에 따라 지침을 정의할 책임이 있습니
 다. 운영 회사가 제공하는 지침을 엄격히 준수해야 합니다.
- 보관 및 작동을 위한 올바른 환경 조건을 준수하십시오.
- 안전 장치의 구조적 변경은 금지됩니다.
- 손상된 안전 장치는 이 설명서의 설명에 따라 즉시 교체해야 합니다.
- 이 Fluent 제품은 Tecan의 사전 자문 및 서면 승인 없이 어떠한 방식으로도 개 조해서는 안 됩니다. 승인된 시스템 개조는 Fluent의 수리 및 업그레이드에 대 해 인증을 받은 FSE만 수행할 수 있습니다. Tecan에서는 무단 개조로 인해 발 생하는 모든 클레임을 거부합니다.
- Fluent의 부적절한 사용으로 인한 화재 위험이 있습니다. 폭발 위험이 있는 곳 에 Fluent 제품을 설치해서는 안 됩니다.
- 가연성 액체 또는 시스템 액체로 인한 화재 위험이 있습니다.
- 가연성 증기가 형성되거나 축적되지 않도록 하십시오.
- 화학적, 생물학적 및 방사능 위험은 사용된 물질 또는 Fluent에서 처리되는 검 체 및 시약과 관련될 수 있습니다(예: 로딩 및 언로딩 중). 폐기물 처리에도 동 일하게 적용됩니다.
 - 이러한 물질과 관련된 잠재적 위험을 항상 인지하십시오.
 - 적절한 보호복, 보안경, 방독 마스크 및 장갑을 사용하십시오.
 - 물질 취급 및 폐기물 폐기는 보건, 환경 또는 안전에 관한 지역, 주 또는 연 방 법률의 적용을 받습니다. 해당 규정을 엄격히 준수하십시오.
- 모든 오염은 본 설명서에 설명된 대로 즉시 처리해야 합니다.
- 사용자는 Fluent 제품이 항상 적절한 조건에서 작동하고 유지 관리, 서비스 및 수리 작업이 일정에 따라 승인된 직원에 의해서만 수행되도록 할 책임이 있습 니다.
- 부정확한 측정 결과의 위험. 시스템 관리 또는 유지 관리를 수행한 후에는 올바 른 시스템 작동 상태를 확인한 후에만 작동을 재개해야 합니다.
- 시스템 성능 및 신뢰성을 보장하기 위해 항상 유효 기한 내의 권장 소모품을 사용하고 유지 관리 및 수리를 위해 정품 예비 부품을 사용하십시오.
- 피부가 기기의 시스템 액체에 접촉할 경우 부상을 입을 수 있습니다.
 - 항상 GLP에 다라 보호복을 착용하십시오.

•TECAN•

- 무거운 적재물입니다! 기기를 들어 올리지 마십시오.
- 데크 트레이 및 데크 세그먼트 없이 시스템을 작동하지 마십시오.
- 데크 트레이는 수동 데크 로딩 영역에서 발생할 수 있는 액체 유출물을 포집합 니다. 모든 액체 유출물을 수집하려면 데크 아래에 가능한 한 많은 데크 트레이 를 설치한 상태로 시스템을 작동해야 합니다. 데크 트레이 없이 시스템을 작동 하지 마십시오.
- 유출이 허용되지 않는 경우 필터와 함께 일회용 팁을 사용할 것을 강력히 권장 합니다.
- 충돌 가능성. 데크 위에 Tecan 모델 데이터가 없는 장치를 배치하지 마십시오.
- Extension 300은 최대 40kg(88파운드)의 하중을 견딜 수 있도록 설계되었으며 40kg(88lbs.) 이하의 가벼운 옵션에서만 사용할 수 있습니다.
- Fluent에는 생물학적 위험이 있는 물질을 사용할 때 사용자가 붙여야 하는 생 물학적 위험 물질 안전 표시가 포함되어 있습니다. 사용자가 볼 수 있고 적용할 때 편리하도록 전면 도어에 라벨을 부착합니다. 제품 안전 표지판 섹션을 참조 하십시오.
- 의 작업대에서 사용하는 옵션은 강한 자기장을 생성할 수 있으며, 이로 인해 심 박 조율기 또는 인슐린 펌프와 같이 운영자가 착용하거나 신체에 이식된 의료 장치의 기능을 방해할 수 있습니다. 제품에는 강력한 자기장에 관한 안전 표지 판이 함께 제공됩니다. 이 표지판은 강력한 자기장을 생성하는 옵션을 사용할 경우 사용자가 볼 수 있고 적용할 때 편리하도록 사용자가 전면 도어에 부착해 야 합니다.
- DeckCheck 카메라의 이더넷 케이블은 FSE에 의해 설치되며 항상 Fluent PC(EMC)에 설치되어야 합니다. 이더넷 인터페이스는 네트워크에 연결할 수 없습니다.
- 캘리포니아주 거주자만 해당: 이 제품은 캘리포니아주에서 암, 기형아 출산 및 생식 피해를 유발하는 물질로 규정된 납 등 화학 물질에 사용자가 노출될 수 있 습니다. 자세한 내용은 www.P65Warnings.ca.gov/product를 참조하십시오.

2.3 카메라 개인정보 처리방침

이 Fluent 시스템에는 내부 전면 프로필에 장착된 카메라가 장착되어 있습니다. 카 메라는 상부 데크와 하부 데크에 각각 초점을 맞추고 있습니다. 아크릴 유리 측면 패널을 통해 아래로 볼 수 있습니다.

- 사용자는 카메라가 작동 중인 방에 있는 사람들에게 조언할 책임이 있습니다.
- 사용자는 기기가 데스크 공간에 측면으로 인접한 경우, 후면이나 측면 패널 컷 아웃이 발생한 경우 또는 후면 벽을 교체하기 위해 아크릴 유리 패널을 사용하 는 경우와 같이 촬영한 사진에서 직원이 식별될 수 없도록 해야 합니다.

2.4 애플리케이션 위험

시스템 기능/ 모듈	가능한 고장 모드	잠재적인 장애 효과	가능한/잠재적인 원인	라벨링 또는 완화 조치
시스템	불충분한 유 지보수	사용자의 안전 또는 건강: 기기 의 잠재적 오염	사용 오류: 작동 설 명서 또는 유지보수 지침 무시	사용자는 적절한 소모품 사용 및 예방 유지보수 지침을 확인해야 합 니다(참조 시스템 지원).
				사용자는 GLP 및 일반 현지 규정에 따라 보호복, 장갑 및 보안경을 착 용해야 합니다.



시스템 기능/ 모듈	가능한 고장 모드	잠재적인 장애 효과	가능한/잠재적인 원인	라벨링 또는 완화 조치
시스템	화재	사용자의 안전 또는 건강: 실험 실 내 화재(기기 연소)	휘발성 가연성 액체 의 가스, 전자 보드 확산으로 인한 스파 크	이 기기는 방폭 장치가 아니므로 증기 농도가 높지 않은지 확인해야 합니다(참조 일반 안전 정보).
모듈 FCA 및 에어 FCA	Z축 기계 마 모(평균 사용 량 초과)	검체의 안전성 또는 임상 상태: 실험 기구 내 Z- 위치 설정이 잠 재적으로 잘못됨	기기와 일회용 팁을 함께 사용할 때의 평균 사용량이 초과 되는 경우 적용 시 피어싱 단 계의 높은 비율	Z축이 축 예상 수명의 90%에 도달 한 경우 시스템에서 이를 사용자에 게 알립니다.
모듈 FCA 및 에어 FCA	P축 기계 마 모(평균 사용 량 초과)	검체의 안전성 또는 임상 상태: 실험 기구 내 P- 위치 설정이 잠 재적으로 잘못됨	기기와 일회용 팁을 함께 사용할 때의 평균 사용량이 초과 되는 경우 적용 시 피어싱 단 계의 높은 비율	시스템은 P축이 축의 예상 수명 90%에 도달했음을 사용자에게 알 려줍니다.
모듈 FCA 및 에어 FCA	X-드라이브 톱니바퀴 마 모(평균 사용 량 초과)	검체의 안전성 또는 임상 상태: 폴리아미드 입자 에 의해 샘플이 오염될 가능성	기기 후면에 주요 실험 기구를 배치한 경우 장치의 평균 사용량 초과	실험 기구 상단에 입자 보호 도구 (예: 뚜껑)를 배치합니다. 또는 기 기 후면에 입자에 민감한 요소(예: 검체 및 시약)를 배치하지 않습니 다.
모듈 FCA 및 에어 FCA	격막 피어싱 으로 인해 발 생하는 간섭 신호	환자 검체의 안 전성 또는 임상 상태: 잘못된 cLLD로 인한 공 기 흡입 및 잠재 적으로 잘못된 결과	팁과 격막/호일 간 의 상호작용	FCA 및 에어 FCA에서 액체 레벨 감지 기능과 결합하여 피어싱 적용 에 비전도성 호일만 사용합니다. 참조 설명서를 참조하십시오. 사용자는 FCA 및 에어 FCA에서 피어싱 기능과 결합하여 액체 감지 기능을 검증해야 합니다.
모듈 FCA 및 에어 FCA	시약 용기 내 거품 또는 기 포로 인한 잘 못된 검체 처 리, 잘못된 cLLD	공정 안전: 잘못 처리된 검체	시약 용기 내 기포 또는 거품으로 인해 cLLD 기능이 잘못 되고 FCA 또는 에 어 FCA로 공기를 흡입될 수 있습니다	사용자는 적절한 검체 준비와 관련 하여 애플리케이션/프로세스를 검 증할 책임이 있습니다.
모듈 FCA 및 에어 FCA	팁 차단	환자 검체의 안 전성 또는 임상 상태: 잠재적으 로 잘못된 피펫 팅 용량	웰 하단에서 흡입 (팁 차단)	사용자 지정 실험 기구 Z-max 레 벨에 너무 가깝게 흡입되지 않도록 적용 수준을 확인해야 합니다.
모듈 FCA 및 에어 FCA	FCA 튜브 삽 입 시스템: 미생물의 성 장	환자 검체의 안 전성 또는 임상 상태: 잘못된 피 펫팅 용량 또는 검체 오염	미생물 증식(안쪽 표면에 생물막)	탈이온수를 FCA의 시스템 액체로 사용하고 일일 유지 관리를 실행하 여 일일 시스템 관리 지침(시스템 지원 참조)에 따라 시스템을 세척 합니다. 이때 이 단계에서 허용되 는 세정제도 확인합니다.



시스템 기능/ 모듈	가능한 고장 모드	잠재적인 장애 효과	가능한/잠재적인 원인	라벨링 또는 완화 조치
모듈 MCA 96 및/또는 MCA 384/96	마이크로플 레이트 내부 에서 피펫팅 과정 중에 검 체 액체가 과 도하게 넘쳐 나는 현상입 니다.	환자 검체의 안 전성 또는 임상 상태: 검체의 잠 재적 교차 오염 (물 넘침)	사용자에 의해 잘못 정의된 Z-수준(예: Z-max 위치에서 흡입)	흡입 및 분사를 위한 안전한 위치 를 정의합니다. 참조 설명서를 참 조하십시오.
모듈 MCA 96 및/또는 MCA 384/96	검체가 자유 분사 모드에 서 의도한 위 치를 완전히 또는 부분적 으로 벗어남	환자 검체의 안 전성 또는 임상 상태: 잠재적 교 차 오염	지정된 조건 이외의 기기 사용으로 인해 팁 끝의 정전기가 발생할 경우 검체가 끝에 남아 매달려 있거나 분무 기능이 제어되지 않음다	사용자는 MCA 액체 취급에 대해 지정된 작동 조건, 특히 최소 필수 습도에 대한 지침을 준수해야 합니 다("환경 조건" [▶ 41] 참조). 사용자는 항상 웰 내부에서 분배 높이를 설정해야 합니다. 참조 설 명서를 참조하십시오.
모듈 MCA 96 및/또는 MCA 384/96	혼합 피펫팅 을 위해 액체 (검체/시약) 대신 공기 혼 합	환자 검체의 안 전성 또는 임상 상태: 검체가 잠 재적으로 잘못 처리되어 잘못된 결과가 발생할 수 있습니다	팁과 마이크로플레 이트의 잘못된 조합 으로 인해 추적 매 개변수가 부적절합 니다	사용자는 가상 작업대의 실험 기구 이름을 참조하여 실제 작업대와의 일치 여부를 확인해야 합니다. 사용자는 DiTi 박스의 고유한 색상 디자인(팁 유형별) 및 라벨 표시(필 터 및 비필터의 경우)를 준수해야 합니다. 사용자는 프로세스를 시작하기 전 에 작업대 레이아웃을 확인해야 합 니다.
일회용 팁에 만 해당	DiTi 가져오 기: 잘못된 팁 유형이 장 착됨	환자 검체의 안 전성 또는 임상 상태: 흡입된 검체가 없거나 부족함 검체의 교차 오 염 가능성	사용 오류: 잘못된 데크 레이아 웃: 팁 상자를 잘못 된 위치에 놓음: 팁 이 예상보다 짧음 잘못된 데크 레이아 웃: 작업대에 여과 된 팁 대신 여과되 지 않은 팁을 포함 한 팁 상자를 놓음 잘못된 데크 레이아 웃: 팁 박스를 잘못 된 위치에 놓음: 팁 부피가 예상보다 작 음(예: 200µl 대신 100µl), 예상 수준 의 팁 길이, MCH에 서 흡입된 액체	사용자는 가상 작업대의 실험 기구 이름을 참조하여 실제 작업대와의 일치 여부를 확인해야 합니다. 사용자는 DiTi 박스의 고유한 색상 디자인(팁 유형별) 및 라벨 표시(필 터 및 비필터의 경우)를 준수해야 합니다. 사용자는 프로세스를 시작하기 전 에 작업대 레이아웃을 확인해야 합 니다. 기계식 설계로 백색 필터를 볼 수 있습니다. 참조 설명서에는 DiTi 상자의 색상 코딩과 길이 차이 및 여과된 DiTi에 대한 정보가 포함되어 있습니다. 참조 설명서를 참조하십시오.



시스템 기능/ 모듈	가능한 고장 모드	잠재적인 장애 효과	가능한/잠재적인 원인	라벨링 또는 완화 조치
일회용 팁에 만 해당	불완전한 팁 드롭: 일부 오염된 팁이 여전히 헤드 위치에 매달 려 검체 플레 이트에 떨어 짐	환자 검체의 안 전성 또는 임상 상태: 잠재적 교 차 오염	정전기 발생	사용자는 MCA 액체 취급에 대해 지정된 작동 조건, 특히 최소 필수 습도에 대한 지침을 준수해야 합니 다("환경 조건" [▶ 41] 참조). 일회용 팁은 재사용할 수 없습니 다.
모듈 RGA	잘못 정렬된 실험 기구와 의 충돌로 인 한 플레이트 손실	공정 안전성: 플 레이트 손실, 샘 플 손실	4개 이상의 마이크 로플레이트가 쌓이 면 운반 중에 정렬 오류가 발생할 수 있습니다	실제 샘플로 스크립트를 실행하기 전에 플레이트 이동을 검증해야 합 니다. 애플리케이션 소프트웨어 설 명서의 검증 체크리스트를 참조하 십시오.
모듈 FluentContro I 소프트웨어	WorktableBa se: 잘못된 DiTi 상태가 보고되었습 니다	공정 안전성: 교 차 오염/잘못된 결과	팁의 사용 상태에 대한 잘못된 정보로 인한 교차 오염	결함 모드로 인해 위험 심각도가 높아지면 "팁 뒤로 설정"을 사용하 지 마십시오.
모듈 FluentContro I 소프트웨어	코어 스크립 팅 프로그래 밍 실행 중 변수 설정: 잘못된 값	공정 안전: 잘못 된 결과	소프트웨어 오류: 변수가 잘못된 값으 로 설정되었습니다	특정 변수 소스, 대상 및 범위에 대 한 애플리케이션의 유효성을 검사 합니다. 애플리케이션 소프트웨어 설명서의 검증 체크리스트를 참조 하십시오.
모듈 FluentContro I 소프트웨어	런타임 또는 스크립트 실 행 시 Core.Scripti ng.Program ming QueryVariabl e에서 UI 값 이 잘못 표시 되거나 수락 되는 문제가 발생합니다.	공정 안전: 잘못 된 결과	숫자 값이 UI에서 잘못 포맷되거나 변 환되었습니다	특정 변수 소스, 대상 및 범위에 대 한 애플리케이션의 유효성을 검사 합니다. 애플리케이션 소프트웨어 설명서의 검증 체크리스트를 참조 하십시오.
모듈 FluentContro I 소프트웨어	코어 스크립 팅 프로그래 밍에서는 런 타임에 ImportVariab le를 사용하 여 잘못된 값 을 가져옵니 다.	공정 안전: 잘못 된 결과	가져오기 소스에서 잘못된 값이 검색되 었습니다	특정 변수 소스, 대상 및 범위에 대 한 애플리케이션의 유효성을 검사 합니다. 애플리케이션 소프트웨어 설명서의 검증 체크리스트를 참조 하십시오.



시스템 기능/ 모듈	가능한 고장 모드	잠재적인 장애 효과	가능한/잠재적인 원인	라벨링 또는 완화 조치
모듈 FluentContro I 소프트웨어	런타임 시 Core.Scripti ng.Program ming ExportVariab le: 잘못된 값 을 파일로 내 보냄	공정 안전: 잘못 된 결과	파일을 내보내기 위 해 잘못된 값이 작 성되었습니다	특정 변수 소스, 대상 및 범위에 대 한 애플리케이션의 유효성을 검사 합니다. 애플리케이션 소프트웨어 설명서의 검증 체크리스트를 참조 하세요.
모듈 FluentContro I 소프트웨어	API: 변수 가 져오기/설정 또는 표현식 확인 실패	공정 안전: 잘못 된 결과	잘못된 변수 값이 검색되었거나 잘못 된 값이 할당되었거 나 잘못된 표현식 결과가 반환되었습 니다	특정 변수 소스, 대상 및 범위에 대 한 애플리케이션의 유효성을 검사 합니다. 애플리케이션 소프트웨어 설명서의 검증 체크리스트를 참조 하십시오.
UVC 조명	애플리케이 션에서 잘못 된 사용	효과 부족	애플리케이션에서 잘못된 사용	"광학 방사선(UVC)" [▶ 29] 섹션의 구체적인 지침을 참조하십시오.
튜브 로테이 터/피어싱 팁 (Mix & Pierce)	애플리케이 션에서 잘못 된 사용	효과 부족	애플리케이션에서 잘못된 사용	"Mix & Pierce" [▶ 58] 섹션의 구체 적인 지침을 참조하십시오.
Frida Reader	애플리케이 션에서 잘못 된 사용	효과 부족	애플리케이션에서 잘못된 사용	"Frida Reader" [▶ 61] 섹션의 구체 적인 지침을 참조하십시오.
모두	애플리케이 션에서 비효 율적인 사용	적용에 효과가 부족합니다	시스템 관리 부족	시스템 지원 장의 부품별 지침을 참조하십시오.
잠재적 위험 물질 처리	오염 물질	사용자, 자산 및 환경에 대한 잠 재적 위험.	일반 안전 정보 준 수 부족	일반 안전 정보 섹션의 부품별 지 침을 참조하십시오.
MCA 96, cLLD	측정 결과가 올바르지 않 습니다	잘못된 측정: 감 지된 액체 레벨 은 마이크로플레 이트의 모든 웰 에 적용되지 않 습니다. 검사 결 과가 잘못되었거 나 검체 손실이 발생했습니다.	부적합한/비호환 실험실 제품 사용	그러면 cLLD 기능은 수조에서만 사용할 수 있습니다.



시스템 기능/ 모듈	가능한 고장 모드	잠재적인 장애 효과	가능한/잠재적인 원인	라벨링 또는 완화 조치
MCA 96, 일 반	소모품/구성 요소/모듈의 비호환성	처리 과정에서 DiTi 손실이 발 생했습니다. 잠 재적으로 복구 불가능한 검체 손실. 잠재적 교 차 오염.	부적합/비호환 소 모품 사용. DiTi를 장착할 상자 또는 트레이와 호환되지 않는 오프셋 픽업 명령이 지시되었습 니다(예: DiTi 트레 이의 잘못된 버전). DiTi가 부적절하게 장착되었고 공정 중 에 분실되었습니다	Tecan 콤비 트레이와 함께 DiTi 상 자를 사용하십시오.
MCA 96, cLLD	측정 결과가 올바르지 않 습니다	잘못된 측정: cLLD 하위 시스 템의 위양성 감 지.	사용자가 소프트웨 어에 반영되지 않은 다른 DiTi를 작업대 에 놓습니다. 사용 자의 조치(예를 들 어 사용자가 일부 DiTi를 삭제하는 등)으로 인해 소프 트웨어가 예상하는 DiTi의 수와 실제로 설치된 DiTi의 수가 다릅니다. 잘못된 cLLD 임계치로 인 해 위양성 감지.	중요한 점입니다. cLLD 검출에 사 용되는 DiTi의 수는 소프트웨어에 반영된 수와 일치해야 합니다.
MCA 96, 일 반	표본 오염	충돌 후 누출로 인한 검체 오염. 잘못된 검사 결 과 또는 검체 손 실.	잘못된 입력 매개 변수(픽업 또는 드 롭 위치를 잘못 설 정하거나 도구 유형 을 잘못 선택한 경 우) 벡터 이동 중 MCA 96 헤드 충 돌. 예를 들어 주변 네스트에 높은 실험 기구가 있는 경우.	벡터 이동 기능의 잘못된 사용과 관련된 위험이 있습니다.
MCA 96, 일 반	표본 오염	환자 검체의 안 전성 또는 임상 상태: 표본 오염: 그리퍼 벨트의 마모로 인한 교 차 오염 가능성	수명 기간 동안 벨 트 재질의 마모: G 축 벨트의 마모 입 자가 작업대의 환자 검체에 떨어져서 최 종적으로 검체가 화 학적으로 오염될 수 있습니다.	벨트의 마모 입자 및 먼지가 작업 대에 떨어져 검체/화학 물질을 오 염시킬 수 있습니다. 이를 방지하 려면 뚜껑을 닫아 민감한 액체가 들어 있는 실험 기구를 보호해야 합니다.



시스템 기능/ 모듈	가능한 고장 모드	잠재적인 장애 효과	가능한/잠재적인 원인	라벨링 또는 완화 조치
MCA 96, 일 반	표본 오염	환자 검체의 안 전성 또는 임상 상태: 잠재적으 로 잘못된 결과: 인접한 플레이트 캐비티로 누출되 어 교차 오염 가 능성이 있습니 다.	검체 액체가 튀거나 누출됨 그리핑된 실 험 기구의 액체에 의한 DiTi 콘 오염 (예: 전체 실험 기구 가 폐기함에 들어가 고 액체가 유출된 경우).	폐기하기 전에 먼저 실험 기구를 비우십시오.
MCA 96, 일 반	검체 또는 시 약 유출(FC)	DiTi 콘형 및 실 린더 블록 오염 으로 인해 잠재 적으로 발생하는 검체의 교차 오 염.	장치가 DiTi를 픽업 하려고 시도합니다. 소프트웨어에서 누 락된 DiTi를 감지할 수 없으므로 오류 없이 계속 진행합니 다. DiTi 콘은 잠재 적으로 수조와 접촉 한 상태로 액체를 흡입합니다.	팁의 가장자리가 상자의 상단 표면 과 같은 높이인 팁 상자 유형을 사 용할 경우 MCA 96에서 빈 팁 상자 를 감지할 수 없습니다.
MCA 96, 일 반	검체 액체가 튀거나 누출 되는 경우	충분히 튼튼한 실험 기구가 부 족하여 발생하는 누출로 인한 검 체 오염.	충분히 튼튼하지 않 은 실험 기구 사용 으로 인한 검체 액 체가 튀기거나 누출 되는 문제.	실험 기구의 변형을 방지하려면 액 체를 운반할 수 있을 만큼 충분히 단단한 실험 기구만 사용하십시오.
MCA 96 일 반	충돌 후 그리 퍼 핑거가 손 상될 수 있습 니다	그리퍼 핑거가 손상됨		그리퍼 핑거를 확인하고 손상된 경 우 교체합니다.
세척 스테이 션 Mix & Pierce	표본 오염	폐기물이 올바르 게 배출되지 않 았습니다. 잘못된 폐기물 배출 및 잘못된 팁 세척으로 인 한 검체 오염.	세척 스테이션 연결 부가 막혀 세척 스 테이션에서 물이 넘 쳤기 때문에 팁이 잘못 세척되었습니 다.	시스템을 정기적으로 유지 관리해 야 합니다. 세척 스테이션 커넥터는 막힘을 방 지하기 위해 전혈 작업 시 2~3개 월 주기로 교환해야 합니다. 세척 스테이션, 커넥터 및 튜브를 통해 부식성 액체(예: 2%의 표백 제)를 배출할 경우 부식을 방지하 려면 물과 같은 중성 액체를 사용 해 추가적으로 헹구어야 합니다.

2.5 운영 회사

운영 회사는 Fluent 및 특히 안전 기능, 기능이 제대로 작동하는지 여부 그리고 기 기와 접촉하는 모든 직원이 적절한 교육을 받았는지 확인해야 합니다.

책임

- 검사법 및 프로세스 검증.
 - 표준 운영 절차에 따라 프로세스 정의.
 - 설치 및 운영 적합성(IQ OQ)이 완료되었는지 확인.

핵심 운영자의

책임



- Fluent 제품과 접촉하는 모든 직원이 적절한 교육을 받는지 확인.
- 적절한 보호복과 장비를 사용할 수 있는지 확인.
- Fluent 제품의 유지 관리 및 안전한 작동 확인.
- 실험실 안전 규정 및 지침 준수.

2.6 검사법 및 프로세스 검증

검사법 및 프로세스 검증을 수행하는 동안 다음 사항에 유의하십시오.

- MCA 384 또는 FCA에서 고정 팁을 사용할 경우 세척 절차가 예상 검체 농도 범위 및 검사 민감도에 대해 유효한지 확인합니다.
- 피펫팅된 용량이 자동화 중인 공정의 정밀도와 정확도 요건을 충족하는지 확인 하십시오.
- Tecan 이외 제품 또는 사용자 지정 실험 기구를 사용하고 추적 기능을 켜고 흡 입할 경우 공기 흡입을 방지하기 위해 용기 관련 정의가 올바른지(예: 추적에 적절한 속도가 사용됨) 확인합니다.
- 위상 분리기 기능은 표준 Tecan 1ml 일회용 팁 및 1ml Tecan 와이드 보어 일 회용 팁과 함께 사용하도록 검증되었습니다. 지원되는 Tecan 소모품에 대한 자세한 내용은 참조 설명서(참조 "참조 문서" [▶ 10])를 참조하십시오.
- Fluent Stacker 전송 스테이션에서 액체 감지를 확인합니다.
 - 애플리케이션에서 MCA 세척 스테이션을 올바르게 사용하고 있는지 검증합니다.
 - 피펫팅 용량이 정확하고 추적이 가능한지 확인하여 애플리케이션을 검증합니다.
 - 사용자 지정 실험 기구 Z-max 레벨에 너무 가깝게 흡입되지 않도록 적용 수준 을 검증합니다.
 - 필요한 다운홀더(활성 또는 수동)에 따라 피어싱 작업을 검증합니다.
 - 화학물질 및 실험 기구를 제거하지 않은 경우 데크에 있는 화학물질과 실험 기 구에 대한 UVC 조명의 영향을 평가하고 검사를 검증해야 합니다.
 - 피펫팅 용량이 올바른지 확인하는 수동 실행 후 점검을 포함합니다.
 - 직원들에게 카메라 개인정보 취급 방침에 대한 정보를 제공해야 합니다("카메 라 개인정보 처리방침" [▶ 15] 참조).

2.7 사용자 자격 평가

실험실 직원은 Fluent 작동을 위해 완전한 자격을 갖추고 교육을 받아야 합니다. 이 Operating Manual(작동 설명서)에 설명된 작업은 아래에 명시된 자격을 갖춘 허가 받은 직원만 수행해야 합니다.

실험실 직원은 다음을 수행해야 합니다.

- 적절한 기술 교육 이수.
- 실험실 안전 규정 및 지침 숙지.
- 기기의 안전 요소 지침 숙지.
- 보호복과 장비 사용.
- 모범 실험실 관행의 숙지 및 준수,
- 작동 설명서의 지침을 읽고 이해.



Tecan은 운영자가 운영자 교육 과정을 수강할 것을 권장합니다. Tecan 고객 서비 스 부서에 수강 가능한 강좌에 대해 문의해 주십시오. "고객 지원" [▶ 188] 섹션을 참조하십시오.

2.7.1 운영자

운영자(실험실 기술자)는 운영 회사에서 근무합니다.

• 특정 애플리케이션 또는 시스템에 대한 지식이 필요하지 않음

필수 기술

필수 기술

책임

- 현지 언어 구사
- 영어 사용이 선호됨

운영자는 애플리케이션 소프트웨어 액세스 권한을 가지고 있어 검사법을 실행하고 시스템 관리를 수행할 수 있으며 핵심 운영자로부터 필요한 교육을 받을 수 있습니 다.

2.7.2 핵심 운영자

핵심 운영자(애플리케이션 전문가)는 운영 회사를 지원하거나 동일한 회사에서 근 무합니다.

- 광범위한 애플리케이션 지식
 - 제한된 시스템 지식
 - 현지 언어 구사
 - 영어 구사
 - 해당 소프트웨어 설명서에 대한 심층적인 지식
- 운영자에게 지시
 - 검사법 작성, 실행 및 유효성 검사
 - 운영자가 기기의 문제를 해결하도록 지원



2.8 안전 요소

▲ 주의

구동 부품

작업 중에 Fluent에 설치된 보호 및 안전 요소를 제거하거나, 비활성화하거나, 무 시해서는 안 됩니다.

장치를 제거한 경우(예: 유지 관리 작업), 작업을 다시 시작하기 전에 모든 보호
 및 안전 장치를 다시 설치하고 다시 활성화하고 점검해야 합니다.

안전 패널 및 안전 센서는 Fluent의 핵심 부품인 반면, 장비 도어 잠금 장치 및 캐 비닛 도어 잠금 장치는 특정 시스템 구성에만 포함될 수 있습니다.

2.8.1 안전 패널

Fluent 안전 패널로 보호됩니다.

전면 안전 패널은 열 수 있으며 활성 정지를 트리거하는 도어 센서가 장착되어 있 습니다. 전면 안전 패널은 도어 잠금 장치(옵션)로 잠글 수 있습니다.

MCA 96 또는 MCA 384 암이 설치된 Fluent는 전면 안전 패널에서만 사용할 수 있 습니다.

도어 개구부는 가스 충전 스프링으로 지지됩니다. 사용자는 장비에 안전하고 완전 하게 접근할 수 있도록 도어를 완전히 열어야 합니다.

희석기 패널은 Fluent 작동에 영향을 주지 않고 열 수 있습니다(UVC 조명 옵션이 설치된 Fluents 제외 - 희석기 패널 센서는 패널이 열릴 때 빠른 중지를 트리거함).

상단 및 측면 안전 패널은 고정되어 있습니다.

2.8.1.1 전면 안전 패널

전면 안전 패널은 작동 중에 로봇 암과 기기 데크의 요소에 직접 접근하는 것을 방 지합니다. 이는 개인 안전을 위한 것으로 검사법 안전을 강화합니다. 또한 전면 안 전 패널은 검체 또는 시약 누출로부터 사용자를 보호합니다. 다양한 유형의 전면 안전 패널이 있습니다.



2 – 안전 안전 요소

전체 전면 안전 패널



• 구동 부품(구동 부품, 기계적 위험)에 접근할 수 없음

- 외부 영향으로부터 검체 보호(검사법 안전)
- 검체 또는 시약 누출로부터 보호



전체 전면 안전 패널의 경우 배치 방식 로딩만 가능합니다.



전체 전면 안전 패널(UVC)



그림 2: 전체 전면 안전 패널(UVC)

전체 전면 안전 패널(UVC)에는 다음과 같은 기능이 있습니다.

- 구동 부품(구동 부품, 기계적 위험)에 접근할 수 없음
- 외부 영향으로부터 검체 보호(검사법 안전)
- 검체 또는 시약 누출로부터 보호
- 광학 방사선(UVC)으로부터 보호



전체 전면 안전 패널의 경우 배치 방식 로딩만 가능합니다.

▲ 주의

구동 부품!

MCA, FCA 및 에어 FCA를 이동할 때 실행 중 기기 절반 전면 안전 패널을 지나거 나 연장부가 포함된 전면 안전 패널에 닿아 기기 안에 손이 들어와 다칠 수 있습니 다.

• 실행 중에 기기 안으로 손을 넣지 마십시오.



2 – 안전 안전 요소

절반 전면 안전 패널



그림 3: 절반 전면 안전 패널

절반 전면 안전 패널에는 다음과 같은 기능이 있습니다.

- 구동 부품(구동 부품, 기계적 위험)에 접근할 수 없음
- 검체 또는 시약 누출로부터 보호



절반 전면 안전 패널을 사용하는 경우 운영자는 기기 데크에 접근하지 않도록 합니 다. 패널을 열지 않고도 러너를 로드 및 언로드할 수 있으므로 운영자가 검사법 실 행 중에 검체 또는 시약을 다시 로드할 수 있습니다.



전면 안전 패널 (확장 포함)



그림 4: 전면 안전 패널(확장 포함)

전면 안전 패널에는 다음과 같은 기능이 있습니다.

• 구동 부품(구동 부품, 기계적 위험)에 접근할 수 없음

- 검체 또는 시약 누출로부터 보호
- 전면 DiTi 폐기물 스테이션을 사용할 수 있습니다. 이 스테이션은 데크에서 돌 출되어 있고 전면 안전 패널에서 입구가 아래쪽을 항해야 합니다.



확장부를 포함한 전면 안전 패널의 경우 배치 방식 로딩만 가능합니다.

2.8.1.2 옵션 장치용 안전 패널

옵션 장치를 Fluent 측면에서 추가하거나 제거하는 경우 적절한 측면 안전 패널을 설치해야 합니다. "고객 지원" [▶ 188]을 참조하십시오.

2.8.2 장비 도어 잠금 장치(옵션)

2개의 선택사양 도어 잠금 장치는 전면 안전 패널이 열리지 못하도록 방지하며, 진 행 중인 프로세스를 보호하는 데 도움이 됩니다. 이렇게 하면 공정 실행이 부당하 게 중단되는 문제를 방지할 수 있습니다. 프로세스를 중지하려면 터치스크린을 통 해 일시 중지 요청을 입력할 수 있습니다.



2.8.3 캐비닛 도어 잠금 장치

데크 아래에서 RGA의 장축에 접근할 수 있는 경우, 접근 지점과 가장 가까운 캐비 닛 도어에 도어 잠금 센서 옵션을 장착해야 합니다. 갑판 아래에 액세스 포인트가 두 개 이상 구현되거나 기기 수명 동안 액세스 포인트가 변경되는 경우 액세스 포 인트 근처의 각 도어에는 도어 잠금 센서가 장착되어 있어야 합니다.

기기에 HEPA 후드가 있는 경우 모든 캐비닛 도어에 도어 잠금 센서를 장착해야 합니다.

2.8.4 광학 방사선(UVC)

Fluent에는 HEPA 후드(옵션)가 장착될 수 있습니다. 이 후드에는 UVC 조명 또는 별도의 UVC 조명 옵션이 포함됩니다.

UVC 조명 방사선 노출은 부상을 초래할 수 있으므로 피해야 합니다. 전면 안전 패 널이 열리면 UVC 조명이 자동으로 꺼집니다. UVC 조명 옵션의 경우 희석기 커버 가 열려도 자동으로 꺼집니다. 특수 UVC 내성 안전 패널은 UVC 조명과 함께 Fluent에 설치됩니다.

UVC 조명은 오염 제거 절차에 사용할 수 있습니다. 개별 프로세스에 대한 UVC 사용의 적합성 및 효과성은 사용자가 검증해야 합니다.



또한 HEPA 후드 제조업체에서 제공하는 설명서를 참조하십시오.

2.8.5 외부 도어 잠금 장치

외부 도어 잠금 장치는 외부 인클로저에 설치하는 Fluent에 구현됩니다. 외부 인클 로저의 도어 패널은 Fluent 안전 패널과 캐비닛 도어의 기계적 안전 기능을 대체하 고, 센서가 내장된 외부 도어 도크는 Fluent 전면 안전 패널과 캐비닛 도어의 도어 센서 및 도어 잠금 기능을 대체합니다.



외부 도어 잠금장치는 ActiveStop을 허용하지 않습니다. 프로세스를 중지 또는 일 시 중지하려면 터치스크린을 통해 일시 중지 요청을 입력할 수 있습니다.



2.9 제품 안전 표지판

안전 표지판은 안전을 위해 Fluent에 부착되어 있습니다. 손상, 분실 또는 판독이 불가능한 안전 표지판은 그림과 같이 즉시 교체해야 합니다. 안전 기호의 의미는 "안전 메시지 규칙" [▶ 12] 섹션을 참조하십시오.

표준 기기



그림 5: 표준 기기

UVC





생물학적 위험



그림 7: 내부도



그림 8: 생물학적 위험





반 전면 안전 패 널이 있는 기기 Fluent 에는 생물학적 위험 물질 사용 시 사용자가 적용해야 하는 생물학적 위험 물 질 안전 표지가 포함되어 있습니다. 사용자가 볼 수 있고 적용할 때 편리하도록 전면 도어에 라벨을 부착합니다.



Fluent® Operating Manual, 399706, ko, V2.7



전면 안전 패널 이 있는 기기(확 장 포함)



그림 10: 전면 안전 패널이 있는 기기(확장 포함)

MCA





2 – 안전 제품 안전 표지판



데크 확장



그림 12: 데크 확장

2.9.1 Mix & Pierce 워크스테이션

FCA 안전 실드



그림 13: 안전 실드



2.10 레이저 방사선

Fluent에는 레이저 바코드 스캐너를 장착할 수 있습니다. 이 바코드 스캐너의 레이 저 방사선은 가시 스펙트럼에서 저전력, 평행 빔입니다. 각 바코드 스캐너 및 전체 Fluent 시스템의 레이저 클래스는 해당 하드웨어에 부착된 레이저 안전 라벨에 표 시되어 있습니다.

레이저가 있는 모든 모듈에는 적절한 레이저 안전 라벨이 표시되어 있습니다.

Fluent 기기는 IEC 60825-1:2007 및 IEC 60825-1:2014에 따라 테스트를 거쳐 인증되었습니다.



▲ 주의

Fluent 제품은 레이저 방사선을 방출하는 IEC 60825-1:2014에 따른 클래 스 1 레이저 제품입니다.

레이저 빔으로 인해 눈부심, 섬광맹, 잔상이 발생할 수 있습니다. • 레이저 빔이나 반사된 빔을 응시하지 마십시오.

2.10.1 레이저 방사선 장치

독립형 바코드 스캐너를 장치에 장착할 수 있습니다.

항상 안전 라벨이 바코드 스캐너에 올바르게 부착되어 있는지 확인하십시오:

- 설명 레이저 방사선 라벨(A): 내장형 저전력 레이저 바코드 스캐너가 포함된 IEC 60825-1에 따라 클래스 2 레이저 제품을 식별합니다. 사용자에게 레이저 빔이나 반사되는 광선을 응시하지 않도록 지시하십시오.
- 클래스 2 레이저는 시스템이 실행 중일 때만 작동하며 운영자와의 인터페이스 가 없습니다.

라벨 배치	설명
A LASER RADIATION Do NOT STARE INTO BEAM Dubus duration Stape Pulse duration Stape CLASS 2 LASER PRODUCT IEC/EN 60825-1 : 2007	호텔에 설치된 독립형 바코드 스캐너: 라벨이 스캐너의 아래쪽에 있습니다.
LASER RADIATION DO Utput 85,00 Wavelength 650nm Pulse duration 112,05 CLASS 2 LASER PRODUCT IEC/EN 60825-1: 2007	로봇 그리퍼 암에 장착된 독립형 바코 드 스캐너: 스캐너에 있는 라벨.



라벨 배치	설명
	로딩 ID : 스캐너 하우징의 후면에 있는 라벨입 니다.
	로딩 ID : 스캐너 하우징의 측면에 있는 라벨입 니다.

2.11 오염 제거 고지

일반적인 시스템 관리 및 표준 검사실 규정에 따라 다음 상황에서 Fluent 및 관련 부품과 액세서리의 오염을 완전히 제거해야 합니다.

- Fluent에 대한 유지 관리 또는 서비스 작업을 수행하기 전, 특히 Fluent에 대한 FSE의 개입 전
- 사고 발생 시(예: 충돌, 누출 등)
- Fluent 또는 관련 부품이나 액세서리를 Tecan에 반납하기 전(예: 수리를 위해)
- 보관 전
- 폐기 전
- 일반적으로 Fluent 또는 관련 부품을 해당 위치에서 옮기기 전

기기 소유자는 모든 장비의 효과적인 오염 제거에 대한 전적인 책임을 집니다.

FSE가 Fluent에 개입하기 전에 Fluent 또는 관련 부품이나 액세서리를 Tecan에 반 납하기 전에, 기기 소유자는 반드시 오염 제거 고지 양식을 작성하고 이에 서명하 여 적절한 실험실 실무 지침에 따라 오염 제거가 수행되었음을 확인해야 합니다. 이 양식을 얻으려면 현지 서비스 조직에 문의하고 오염 제거 섹션을 참조하십시오.



Tecan에서는 오염 제거 고지 양식과 함께 제공되지 않는 Fluent 또는 관련 부품이 나 액세서리의 처리를 거부할 권한을 보유합니다.

2.12 사고 보고

장치와 관련하여 발생한 모든 심각한 사고는 제조사와 사용자 및/또는 환자가 거주 중인 회원국의 관할 기관에 보고해야 합니다. 제조업체 주소는 제조업체 섹션을 참 조하십시오. (부작용 보고 관련 문의처 (한국의료기기안전정보원, 080-080-4183).
TECAN

3 기술 데이터

3.1 타입 플레이트



- Vor Service oder Wartungsarbeiten Netzstecker ziehen
- Prior to any repair or maintenance job disconnect mains power cord
- Avant tout type d'intervention, retirer la prise de raccordement au secteur et lire attentivement le manuel
- Prima di esegurie qualsiasi lavoro di manutenzione o servizio, disconnettere il cavo di alimentazione dalla presa di corrente Antes de cualquier intervención de servicio o mantenimiento
- apagar y desconectar el instrumento

https://www.tecan.com/manuals



(01)07640137481124(11)991231(21)9912123456



ACLASS 1. LASER RFODUCT. THIS PRODUCT COMPLIES WITH 21 CFR 1040.10 AND 1040.11 EXCEPT FOR CONFORMANCE WITH IEC 60825-1 Ed. 3., AS DESCRIBED IN LASER NOTICE No. 56, dated May 8, 2019. IEC 60825-1:2014*



그림 14: 타입 플레이트

타입 플레이트는 Fluent 뒷면에 있으며 다음 정보를 포함합니다.

식별 데이터	모델
	REF: 주문 정보(자재 번호 및 개정 수준)
	제조일(YYYYMMDD)
	SN: 일련번호
기술 데이터	U, f: 공급 전압(볼트), 주파수(헤르츠)
	P: 전력 소비량(W)
주소 데이터	제조업체 이름 및 주소
적합성 데이터	적합성 표시





3.2 일련번호 레이블

Model Instrument Fluent 780



그림 15: 일련번호 레이블

일련번호 레이블은 기기 뒷면의 오른쪽에 있는 하우징 안에 부착되며 다음 데이터 가 포함되어 있습니다.

식별 데이터	모델
	REF: 주문 정보(자재 번호 및 개정 수준)
	SN: 일련번호
주소 데이터	제조업체 이름 및 주소



3.3 치수 및 중량



	치수	Fluent 480	Fluent 780	Fluent 1080
а	전체 길이	1150mm(45. 28인치)	1650mm(64. 96인치)	2150mm(84. 65인치)
b	풋프린트 깊이	780mm(30.71인치)		
С	전체 깊이	923mm(36.34인치)		
d	캐비닛 전체 높이	19	977mm(77.8인치	:)

구성 요소	Fluent 480 Fluent 780		Fluent 1080
베이스 유닛	120kg(264.5파운 드)	140kg(308.6파운 드)	190kg(418.9파운 드)
포장	61kg(135파운드)	83kg(183파운드)	106kg(234파운 드)
FCA	10.4kg(22.9파운드)		
MCA 384	12.6kg(27.8파운드)		
384채널 헤드	7.2kg(15.9파운드)		
RGA	10.2kg(22.4파운드)		



구성 요소	Fluent 480	Fluent 780	Fluent 1080
RGA-Z	10.6kg(23.4파운드)		
СХР		1.2kg(2.6파운드)	
MCA 96(옵션 그 리퍼 포함)		19.7kg(43.43파운드))

3.4 전원 공급

공지사항

전원 공급 장치 과열

전원 공급 장치가 손상되거나 파괴될 수 있습니다.

- 전원 공급 장치를 덮지 마십시오.
- 전원 공급 장치의 열 방출이 보장되어야 합니다.



외부 장치를 전원 공급 장치에 연결해서는 안 됩니다. 그러면 이 경우에는 Fluent 가 다시 초기화되거나 일시 중단될 수 있습니다.

표 2: Fluent 전원 입력

공급	등급
라인 전압(단상)	100~240VAC
입력 전류	9.8A(100V에서) - 4A(240V에서)
주파수	50~60Hz

표 3: Fluent 전원 출력

공급	등급
출력 전압	24~28V 기본 설정: 25.2V
지속적인 전력	500W
피크 전력(시간 제한)	3초 동안 15,000W
중량	3.8kg(8.5파운드)

최대 주 전원 공급 전압 변동폭: 공칭 전압 기준 ±10%.

EN/IEC 표준에 따른 전기 안전 관련 분류:

표 4: 전기 사양(안전)

과전압 범주	II	IEC 60664-1
오염 정도	2	(KO) IEC 61010-1





3.5 데이터 및 전원 연결



그림 16: 데이터 및 전원 연결

- A Fluent 7|7|
- C 전원 공급 장치
- I1 USB 인터페이스
- B 제어 PC
- D 벽면 콘센트
 - l2 전원 코드

이 그림은 데이터 및 전원 연결이 포함된 샘플 시스템의 구성 요소를 보여줍니다. Fluent 기기의 각 부분은 사각형 안에 표시됩니다. 기기 전원 스위치는 전원 공급 장치의 일부입니다. 전원 코드는 주 전원 공급 장치의 벽면 콘센트에 연결되어 있 습니다.

USB 인터페이스를 통해 Fluent에서 송수신되는 모든 데이터 트래픽. USB 케이블 은 기기를 제어하는 PC에 연결됩니다.

3.6 환경 조건



잘못된 피펫팅 용량

피펫팅 결과는 작동 조건의 영향을 받을 수 있습니다.

응결이 전자 부품에 영향을 미칠 수 있습니다.

 Fluent를 실온보다 낮은 온도에서 보관하거나 운송하는 경우 설치 후 몇 시간 동안 적응 기간을 거쳐야 합니다.





Fluent 실내 작동 및 보관 전용으로만 사용해야 합니다.

작동 조건

작동 온도	15~32°C(59~90°F)
작동 습도	30°C(86°F)에서 상대(비응축) 30~80%
작동 고도	해발 최대 0~2000m

액체 취급 및 피펫팅을 위한 작동 조건:

실내 온도	20~25°C(68~77°F)
작동 습도	상대(비응축) 30~60%
작동 고도	해발 약 0~500m
증발	층류, 에어컨 또는 환기 등으로 인해 공 기 흐름이 증가하는 환경에서는 특히 소량 또는 휘발성 물질의 경우 증발 위 험이 증가하여 피펫팅 정밀도가 떨어질 수 있습니다. 공지사항! 검증 조건이 실행 조건과 일 치하는지 확인하십시오.

운송 조건

보관 조건

운송 온도	-20~60°C(-4~140°F)
운송 습도	상대(비응축) 20~80%
보관 온도	1~60°C(34~140°F)

3.7 방출 및 내성

보관 습도

소음 방출 기기로부터 1m 거리에서 측정한 < 60 dBA(음압). 프로세스가 실행되는 동안 짧은 시간 동안 소음 수준이 78dB를 초과할 수 있습니 다.

EMC Fluent 제품은 IEC 61326-1 및 IEC 61326-2-6에 설명된 방출 및 내성 요구사항 을 준수합니다. 그러나 전자기 환경은 Fluent의 작동 전에 평가되어야 합니다. Fluent가 의도한 대로 작동하기 위해 Fluent와 호환되는 전자기 환경을 유지할 수 있도록 하는 것은 운영자의 책임입니다. Fluent는 그룹 1 클래스 B 장비(CISPR 11)로 분류됩니다.

> 이 장비는 기본 전자기 환경(IEC 61326-1) 및 전문 의료 시설 환경(IEC 61326-2-6)에서 사용하도록 설계되었습니다. 산업용 전자기 환경(IEC 61326-1)과 가정 의료 설비 환경(IEC 61326-2-6)에서 사용할 경우 제대로 작동하지 않을 수 있습니다. 성능이 전자기 간섭의 영향을 받는 것으로 의심되는 경우 장비와 간섭원 사이의 거 리를 늘려 올바른 작동을 복원할 수 있습니다.

> 강한 전자파 방사원(예: 차폐되지 않은 의도적인 RF 소스) 가까이에서 Fluent 제품 을 사용하지 마십시오. 올바른 작동을 방해할 수 있습니다.

30°C(86°F)에서 상대(비응축) 5~80%



FCC15 본 장비는 FCC 규정 파트 15에 의거, 클래스 B 디지털 장치 허용 한계에 따라 시 험한 결과 이를 준수하는 것으로 판명되었습니다. 이러한 허용 한계는 주거 환경 내 설치 시 유해한 간섭에 대한 합리적 보호를 제공하기 위해 마련되었습니다. 이 장비는 무선 주파수 에너지를 생성, 사용 및 방출할 수 있으며 지침에 따라 설치 및 사용하지 않을 경우 무선 통신에 유해한 간섭을 일으킬 수 있습니다. 그러나 특정 설치에서 간섭이 발생하지 않는다는 보장은 없습니다. 이 장비가 라디오 또는 텔레 비전 수신에 유해한 간섭을 일으키는 경우(이는 장비를 껐다가 켜면 알 수 있음), 사용자는 다음 중 한 가지 이상의 조치를 시도하여 간섭을 해결하는 것이 좋습니 다.

- 안테나 방향을 재조정하거나 재배치합니다.
- 장비와 수신기 사이의 거리를 멀리합니다.
- 수신기가 연결된 회로와 다른 회로의 콘센트에 장비를 연결합니다.
- 대리점 또는 숙련된 라디오/TV 기술자에게 도움을 요청합니다.



4 기능 설명

이 장에서는 Fluent의 기본 기능을 설명하고 구조를 보여 주며 어셈블리의 기능 설 명을 제공합니다.

4.1 개요

Fluent 제품은 로봇 암을 이용한 피펫팅 작업에 사용됩니다. 로봇 암은 검체 튜브 또는 마이크로플레이트와 같은 다양한 용기에서 흡입 및 분사 작업을 수행할 수 있 습니다.

Fluent 제품은 다음의 두 가지 크기로 제공됩니다.

- Fluent 480
- Fluent 780
- Fluent 1080



그림 17: 기기 개요(기기가 그림과 다를 수 있음)

- A 데크
- C 옵션 및 장치



지진이 발생하기 쉬운 지역을 위한 지진 보호 키트도 사용할 수 있습니다. 자세한 내용은 "고객 지원" [▶ 188] 섹션을 참조하십시오.

В

로봇 암

4.2 데크

- **세그먼트** Fluent 기기의 샘플 구역인 데크는 세그먼트로 구성됩니다. 데크 세그먼트는 치수 및 기능이 다양한 교환식 데크 구성품입니다. 주의! 데크 세그먼트 없이 시스템을 작동하지 마십시오.
- **그리드 번호** 세그먼트 너비는 그리드 번호로 표시됩니다. 그리드는 폭이 25mm이며 세그먼트 의 위치 지정 핀 사이 간격에 상응합니다.

그리드 번호는 데크에서 세그먼트 또는 러너의 위치를 나타내는 데도 사용됩니다.



4.2.1 캐리어



그림 18: Fluent 데크

A 러너

B 세그먼트

캐리어는 데크에 실험 기구 또는 소모품을 고정할 수 있도록 고안된 데크 구성품입 니다.

러너는 그리드 세그먼트 슬라이드 기능을 켜고 끄는 캐리어로 일반적으로 검체 튜 브 또는 시약 수조를 고정합니다.

세그먼트는 데크에 고정된 정적 구성요소입니다. 일부 세그먼트에는 마이크로플레 이트나 딥 웰 플레이트와 같은 실험 기구나 DiTi 상자 등의 소모품을 보관하는 네스 트(네스트 세그먼트)가 있습니다. 일부 세그먼트에는 러너를 로드 및 언로드하기 위한 그리드 핀(그리드 세그먼트)이 있습니다.

4.2.2 데크 트레이



그림 19: 데크 트레이



동적 데크 세그먼트 아래 배치되는 데크 트레이는 수동 데크 로딩 영역에서 발생할 수 있는 액체 유출물을 포집합니다. 모든 액체 유출물을 수집하려면 데크 아래에 가능한 한 많은 데크 트레이를 설치한 상태로 시스템을 작동해야 합니다. **주의! 데** 크 트레이 및 데크 세그먼트 없이 시스템을 작동하지 마십시오.

도구 및 기기에서 데크 트레이의 컷아웃은 캐비닛 버전에서만 허용됩니다.

각 데크 트레이의 융기된 부분은 집수 부피에 영향을 주지 않고 컷아웃할 수 있도 록 설계되었습니다. 이러한 컷아웃은 캐비닛 장착 버전을 사용할 때 데크를 통과하 는 폐기물 슈트 또는 기타 장치 통합을 수용합니다. 데크를 통과하는 접근 지점의 경우 트레이를 통합 장치 아래 캐비닛의 선반에 배치할 수 있습니다.



그림 20: 데크 세그먼트 아래 데크 트레이

데크 아래 장치에 RGA가 접근해야 경우 데크 트레이가 없습니다. 데크 트레이 한 세트가 기기에 포함되어 있습니다. 데크 트레이는 필요에 따라 세척하거나 교체할 수 있습니다. "일과 종료" [▶ 109] 섹션을 참조하십시오.

4.2.3 세그먼트에 배치



그림 21: 잠금핀과 위치 지정 핀

잠금핀

А

B 위치 지정 핀

Fluent는 핀을 사용하여 러너, 어댑터 또는 옵션을 세그먼트에 올바르게 배치합니다. 러너는 핀에 밀어 넣을 수 있도록 설계되었습니다. 기기 전면에 있는 그리드 번 호를 읽어서 위치를 확인할 수 있습니다. 잠금 핀은 러너를 제자리에 고정합니다.



4.2.4 세그먼트 위치



그림 22: 측면 위치 및 그리드 위치

A 측면 위치

B 격자 좌표

측면 위치(AB, YZ)를 사용하여 RGA에서 처리하는 실험 기구를 배치할 수 있습니 다.



FCA 또는 MCA를 사용하는 경우에는 측면 위치에서 피펫팅이 불가능합니다.

번호가 매겨진 그리드 위치(1~n)에서 피펫팅 암에 접근할 수 있습니다. 그러나 다 중 암 구성에서 일부 숫자 그리드는 일부 피펫팅 암에 액세스할 수 없습니다. 장치 의 암 구성에 따라 제한이 적용할 수 있습니다.

4.3 로봇 암

Fluent에는 다양한 로봇 암을 장착할 수 있습니다.

- 유연한 채널 암(FCA)
- 다중 채널 암(MCA 96 및 MCA 384)
- 로봇 그리퍼 암(RGA)

로봇 암에는 다양한 암 부속품을 장착할 수 있습니다.





4.3.1 유연한 채널 암(FCA)



그림 23: 유연한 채널 암



유출이 허용되지 않는 경우 필터와 함께 일회용 팁을 사용할 것을 강력히 권장합니 다.

FCA(A)에는 피펫팅 팁이 장착되어 있으며 최대 8개의 개별 채널에 대한 액체 취급 을 제어할 수 있습니다.

DiTi 어댑터로 구성된 FCA에는 특정 실험 기구 이동을 활성화하는 옵션 FCA 그리 퍼가 있습니다("FCA 그리퍼" [▶ 56]).

4.3.1.1 액체 시스템을 갖춘 FCA(Liquid FCA)

액체 변위 시스템을 갖춘 FCA는 주사기 펌프에서 공급되는 시스템 액체로 충전됩 니다. 여러 가지 용량 범위의 액체(사용하는 팁과 주사기 크기에 따라 달라짐)를 피 펫팅하는 데 사용됩니다. 액체 FCA 는 세척 가능한 고정 팁 또는 일회용 팁 어댑터 를 선택하여 구성할 수 있습니다.



Tecan은 시스템 액체로 탈이온수를 사용할 것을 권장합니다.

4.3.1.2 에어 시스템 포함 FCA(에어 FCA)

에어 배치 시스템이 있는 FCA는 피펫팅 채널 내부에서 플런저를 움직여 액체를 피 펫팅하는 데 사용됩니다. Air FCA는 일회용 팁 어댑터로 구성됩니다.



4.3.2 다중 채널 암 384(MCA 384)



그림 24: 다중 채널 암



유출이 허용되지 않는 경우 필터와 함께 일회용 팁을 사용할 것을 강력히 권장합니 다.

MCA(A)는 다채널 피펫팅 헤드를 갖추고 있는 로봇 암입니다. 모든 펌핑 헤드 채널 은 동시에 흡입하고 분사합니다. 피펫팅 헤드는 헤드 어댑터를 교환할 수 있습니 다. 다양한 유형의 헤드 어댑터로 다양한 피펫팅 형식을 사용할 수 있습니다.

- MCA DiTi 포트폴리오와 호환
- 최대 125ul 피펫팅
- 피펫팅 범위 최대 500ul(옵션/확장 용량 어댑터 포함)
- 액체 레벨 감지 없음
- MCA384(일회용 팁 384개 포함)
- MCA384(96개의 일회용 팁 포함)(어댑터 플레이트)
- MCA 384(384개의 고정, 세척 가능한 팁 포함)
- MCA 384(96개의 고정, 세척 가능한 팁 포함)

4.3.3 다중 채널 암 96(MCA 96)

MCA 96는 다채널 피펫팅 헤드를 갖춘 로봇 암입니다. 96개의 피펫팅 헤드 채널 모두 동시 흡입 및 분사가 가능합니다. MCA 96 암은 다음과 같은 핵심 기능을 지 원합니다.





그림 25: MCA 96 헤드/그리퍼

- 최대 1000µL를 포함하는 전도성 FCA-DiTi 포트폴리오와 호환
- 피펫팅 범위: 1ul~1000µl
- (정전식) 액체 수준 감지 시스템(cLLD) 수행
- Tecan 콤비 트레이가 있는 팁 제품을 사용하여 부분 DiTi 픽업 및 오프셋 피펫 팅(예: 희석) 수행
- 간단한 실험 기구의 취급 작업을 위해 옵션으로 제공되는 그리퍼입니다.

MCA 96 헤드는 cLLD와의 호환성을 위해 MCA 96의 전도성 일회용 팁과 함께 작 동합니다.



6

MCA 96의 cLLD는 수조에서 사용하는 경우만 검증되었습니다.

액체로 들어가는 첫 번째 팁은 cLLD 신호를 트리거합니다. 보충 높이가 동일하지 않으면(예: 마이크로플레이트의 웰을 수동으로 보충함) 의도하지 않은 방식으로 처 리되거나 잘못된 결과가 나올 수 있습니다.

6

MCA 96에는 cLLD 호환 캐리어 및 실험 기구만 사용하십시오. 부분 DiTi 픽업을 위해 호환 가능한 트레이 사용: Tecan SLAS 트레이 자세한 내용은 참조 설명서(참 조 [4])를 참조하십시오.

MCA 96의 필드 등급을 구분할 수 없는 그리퍼(옵션)는 간단한 실험 기구 처리 작 업을 위해 사용할 수 있습니다.

그리퍼는 다음과 같은 작업에 사용할 수 있습니다.



- 작업대의 목표 위치에서 마이크로플레이트 및 DWP를 가로 모드로 이동
- 폐기물에 빈 레이어를 넣고 DiTi 트레이를 겹쳐서 작업합니다.
- 마이크로플레이트 뚜껑 취급

MCA 그리퍼는 MCA 헤드에서 하위 모듈로 제작되므로 전용 로봇 그리퍼 암(RGA) 을 대체할 수 있는 비용 효율적인 컴팩트형 로봇 그리퍼입니다. MCA 96 그리퍼는 다음과 같은 복잡한 실험 기구 처리 작업을 수행할 수 없습니다.

- 호텔 액세스
- 세로 방향으로 실험 기구 처리
- 회전을 포함한 작업

이러한 복잡한 작업을 수행하려면 여전히 전용 로봇 그리퍼 암(RGA)을 시스템에 추가해야 합니다. MCA 96 그리퍼는 RGA 그리퍼와 동일한 핑거 교환 스테이션으 로 구성됩니다.

4.3.4 로봇 그리퍼 암(RGA)



그림 26: 로봇 그리퍼 암

RGA(A)는 그리퍼 헤드 및 그리퍼 핑거를 갖춘 로봇 암입니다. RGA는 데크 위치, 주변 장치 및 실험 기구 보관 사이에서 마이크로플레이트 및 기타 실험 기구를 전 송합니다.

표준 높이 로봇 그리퍼 암(RGA 표준 Z)은 데크나 하단 데크에 있는 물체에 접근할 수 있습니다.

높은 로봇 그리퍼 암(RGA 롱 Z)은 데크나 하단 데크 및 그 아래에 있는 물체에 접 근할 수 있습니다.

4.3.4.1 로봇 그리퍼 헤드

RGA 에는 두 가지 로봇 그리퍼 헤드 옵션이 장착될 수 있습니다.

일반 그리퍼 헤드는 수동으로 교환할 수 있는 그리퍼 핑거 옵션을 제공합니다.

핑거 교환 시스템(FES)은 그리퍼 핑거 세트 옵션과 함께 자동 핑거 교환 기능을 제 공합니다. 핑거 세트는 표준 네스트 세그먼트에 장착된 도킹 스테이션에 장착됩니 다. 핑거 세트는 자동으로 픽업되어 로봇 암에 의해 배치됩니다. 핑거 유형 및 핑거 교환은 모니터링됩니다. 한 가지 검사법 내에서 모든 핑거를 사용할 수 있습니다.

TECAN

4.3.5 암액세서리



그림 27: 암 액세서리

- A 고정 팁
- C 그리퍼 핑거

4.3.5.1 고정 팁

Ť

유출이 허용되지 않는 경우 필터와 함께 일회용 팁을 사용할 것을 강력히 권장합니 다.

В

D

일회용 팁

RGA용 바코드 스캐너

흡인 및 분사를 위한 세척 및 재사용 가능 팁을 FCA 및 MCA에 사용할 수 있습니 다.

4.3.5.2 일회용 팁

팁은 유형에 따라 트레이 또는 상자(단일 또는 중첩)에 제공됩니다. 팁이 폐기되거 나 흡인 후 다시 랙에 배치됩니다. 팁은 일회용 팁 배출 시스템을 통해 데크 세그먼 트에 장착된 폐기물 슈트로 폐기됩니다.

4.3.5.3 그리퍼 핑거

핑거 교환 시스템(FES)를 사용하는 RGA 일반 그리퍼 헤드 및 MCA 96 그리퍼에서 다양한 유형의 그리퍼 핑거를 사용할 수 있습니다.

편심 그리퍼 핑 거는 피펫팅 영역 내외의 마이크로플레이트 기반 물체를 이송합니거다. 플레이트 물체를 측면에서 잡습니다. 두 가지 변형 모델을 사용할 수 있습니다.

- 마이크로플레이트를 호텔 및 장치에 로드하는 표준 길이 핑거.
- 더 깊은 장치(예: 4슬롯, 셀 플레이트, 모니터링되는 인큐베이터)를 로딩하기 위한 편심 롱 핑거.
- MCA 96 편심 그
 MCA 96 편심 그리퍼 핑거는 피펫팅 영역 내에서 마이크로플레이트 기반 물체를 이송합니다. 이 핑거 유형은 MCA 96 그리퍼 헤드 아래 플레이트 물체를 잡을 수 있습니다. (이 핑거 유형은 RGA 그리퍼와 호환되지 않습니다.)
- 중심 그리퍼 핑중심 그리퍼 핑중역 내부 및 아래의 마이크로플레이트 기반 물체를 이
송합니다. 플레이트 물체를 위에서 잡습니다. RGA 일반 그리퍼 헤드에만 적용됩니
다.
 - **튜브 핑거** 튜브 핑거는 피펫팅 영역 내부 및 아래의 튜브 기반 물체를 이송합니다. RGA 일반 그리퍼 헤드에만 적용됩니다.



4.3.5.4 바코드 스캐너

RGA는 마이크로플레이트와 DiTi 박스의 바코드를 읽기 위한 수평 스캐너를 장착 할 수 있습니다.



레이저 등급의 안전 지침을 주의하여 읽어야 하며 반드시 따라야 합니다. 바코드 스캐너 제조업체가 제공한 설명서도 참조하십시오.

4.4 액체 시스템(Liquid FCA)



그림 28: 액체 시스템(Liquid FCA)

액체 시스템은 고정 피펫 팁의 내부 및 외부를 효율적으로 세척할 수 있도록 설계 되었습니다.

4.5 세척 시스템(MCA 384)



그림 29: 세척 시스템(MCA)

MCA 384 세그먼트에 설치된 세척 블록(A)은 각 피펫팅 사이클 이후 고정 팁 어댑 터의 팁을 세척합니다.



4.6 옵션 및 장치

1	
	이 섹션에서는 여구 전용 옵션 및 장치에 별표(*)가 표시되어 있습니다
	자세한 내용은 "용도" [》 9] 섹션을 참조하십시오.
패시브 옵션	• 호텔(플레이트 저장 장치)
	• 캐비닛
	• 머지 던개
	• FCA 그리머
화서 오셔	
	• Te-vacs
	• 세적 및 모중 센터(WRC) 타워(예: MCA 세적 스테이션)
	Resolvex i300
	• 피어싱 팁 및 튜브 로테이터
	FRIDA Reader
	자세한 내용은 "참조 문서" [▶ 10] 섹션을 참조하십시오.
	• HydroControl 기반 워셔
	• MT-SICS 레벨 1 표준에 따른 저울
	• SiLA 호환 장치 *
	• Agilent 씰러 *
	Inheco ODTC
	• Inheco의 MTC/STC 컨트롤러를 사용하여 가열 및 냉각 수행
	• Cytomat 10*, 20*, 200*, 6000*
바코드 판독기	• Fluent ID 튜브 바코드 스캐너
	• Keynence BL-1300 시리즈의 바코드 판독기
판독기	• Magellan으로 제어되는 Tecan 판독기
	• Spark 및 SparkControl Magellan*
	• Ziath 2D 플랫베드 판독기*
U	주의 깊게 읽고 따라야 합니다.



4.6.1 Fluent ID 튜브 바코드 스캐너



그림 30: Fluent ID

- A 스캐너 하우징
- C 로딩 영역
- E 리플렉터

- B 레이저 바코드 스캐너
- D LED

Fluent ID는 튜브 러너가 데크에 장착될 때 튜브 바코드 라벨을 스캔하는 옵션 모듈 입니다. 각 Fluent ID 모듈에는 최대 6개 러너의 바코드 라벨을 로드하고 스캔하는 데 사용할 수 있는 최대 6개의 전용 그리드 위치가 포함되어 있습니다. 리플렉터는 러너에서 빈 튜브 위치를 감지하는 데 사용됩니다. 터치스크린 모니터의 그래픽 인 터페이스는 Fluent ID 작동에 대한 지침을 제공합니다.

바코드 스캐너의 레이저 방사선은 다음 속성을 가진 가시 스펙트럼에서 저전력, 평 행 빔입니다.

- 파장: 655nm
- 펄스 지속 시간: 150µs
- 최대 에너지 출력 전력: 1.0mW

4.6.1.1 Fluent ID 호환 튜브 러너

Fluent ID 튜브 러너는 각각 한 가지 유형의 튜브에 맞게 설계되었습니다.

- 10mm 직경의 튜브용 32개 위치가 있는 러너
- 13 mm 직경의 튜브용 32개 위치가 있는 러너
- 16mm 직경의 튜브용 26개 위치가 있는 러너
- 2mL의 Eppendorf Safe-Lock 튜브용 32개 위치가 있는 러너



선택사양 플러그를 사용하여 26위치 러너의 두 위치를 차단하여 24위치 러너로 사 용할 수 있으며, 8의 배수로 튜브에서 병렬 피펫팅을 수행할 수 있습니다.



4.6.2 FCA 그리퍼

개요



그림 31: FCA 그리퍼

FCA 그리퍼는 DiTi 어댑터에 구성된 FCA의 옵션으로, 피펫팅 외에도 FCA에서 일 부 실험 기구를 이동할 수 있도록 지원합니다. FCA는 실행 중에 FCA 그리퍼 핑거 를 자동으로 가져와 드롭할 수 있습니다.

FCA 그리퍼 핑 거



그림 32: FCA 그리퍼 핑거

FCA 그리퍼 핑거는 2년 또는 20000회 사이클 사용 이후에 교체해야 합니다(1회 사이클은 픽업, 사용 및 파킹으로 정의됨). 주기는 Fluent Control 소프트웨어에서 정의된 카운터를 통해 모니터링됩니다.



FCA 그리퍼 도 킹 스테이션 네 스트



그림 33: FCA 그리퍼 도킹 스테이션 네스트

FCA 그리퍼 도킹 스테이션 네스트는 FCA 그리퍼 핑거를 보관하는 데 사용됩니다. 데크 세그먼트에 표준 마이크로플레이트 둥지처럼 장착할 수 있습니다.



4.6.3 Mix & Pierce

Fluent Mix & Pierce 워크스테이션은 고무 캡을 통과하여 피어싱함으로써 캡을 제 거하지 않고도 고무 캡이 장착된 검체 튜브로 또는 해당 검체 튜브에서 액체를 이 송하는 용도로 설계되었습니다.



튜브 격막은 한 번만 뚫을 수 있습니다. 동일한 튜브의 반복적인 피어싱이 지원되지 않습니다.

매끄러운 풀 캡 및 안전 트위스트 캡을 사용하는 Greiner Vacuette® 튜브와 Hemogard 마개를 사용하는 BD Vacutainer® 튜브는 피어싱 테스트를 거쳤습니 다._____

고무 스토퍼/플러그가 있는 튜브의 피어싱이 지원되지 않습니다.



피어싱 팁은 정기적으로 교환해야 합니다. 20000개의 피어싱/팁에 대한 사용이 확인되었습니다.

Fluent Mix & Pierce 워크스테이션은 최대 2개의 액체 FCA, 딥 세척 스테이션 및 최대 4개의 튜브 로테이터로 구성됩니다(Fluent 기본 유닛 크기에 따라 달라짐). 튜브 로테이터는 모든 Fluent 기본 유닛 크기에 통합될 수 있으며 튜브 바코드 판 독, 검체 혼합, 피어싱 및 분주와 같은 작업을 지원합니다. 지원되는 튜브 유형에 대한 자세한 내용은 "튜브 로테이터 러너" [▶ 59] 섹션을 참조하십시오.

워크플로우는 다음 단계로 나눌 수 있습니다.

- 1. 튜브 로딩 중 바코드 스캔
- 2. 튜브 내용물 혼합
- 3. 피어싱 팁을 포함한 튜브 로테이터에서 FCA로 피어싱 및 액체 처리
- 4. 딥 세척 스테이션 및 오염 제거 수조에서 피어싱 팁의 세척 및 오염 제거
- 5. 2단계 반복 및 계속
- 6. 스크립트 끝에서 중앙 폐기물, 전면 및 후면 클리너를 포함한 세척 스테이션의 연장 세척이 권장됩니다.

4.6.3.1 튜브 로테이터

튜브 로테이터 모듈의 기본 용도는 튜브의 액체 내용물을 혼합하고 피어싱 및 피펫 팅 작업을 위한 캐리어 역할을 하는 것입니다. 단일 튜브 로테이터는 5개의 튜브 로테이터 러너를 포함하여 각자 24개의 튜브를 수용할 수 있는 총 120개의 튜브 용량을 가집니다.

장치에는 다음과 같은 하위 구성 요소가 포함되어 있습니다.

- 장착 중에 검체 바코드를 스캔하기 위한 통합형 튜브 바코드 스캐너
- 피어싱 팁과 오류 튜브 홀더의 오염 제거를 위해 깊은 수조를 포함하는 딥 워시 스테이션(옵션). 오류 튜브 홀더는 피어싱 오류가 발생할 경우 검체를 보관하는 데 사용할 수 있습니다. 세척 스테이션이 회전하는 드럼 옆에 배치됩니다.
- 최대 5개의 튜브 로테이터 러너를 수용할 수 있는 튜브 다운 홀더가 있는 회전 식 드럼. 드럼은 다양한 각도와 속도에서 360도 회전 또는 진동을 통해 시료 혼 합을 수행합니다. 다운 홀더(커버)는 피어싱 공정을 지지합니다.
- 튜브 로테이터 러너가 있는 튜브 로테이터는 흡입 전후 및 폐쇄형 튜브(옵션 설 정)를 통한 액체 분사(액체 도착 확인) 후에도 정전식 액체 레벨 감지를 지원합 니다.

⁶



- 튜브 로테이터는 FSE에 의해 설치되며, 핵심 운영자나 사용자가 이동해서는 안 됩니다.

4.6.3.2 튜브 로테이터 러너

튜브 로테이터 러너는 튜브 로테이터에서 사용하도록 설계되었으며 피어싱 기능을 지원합니다. 피어싱을 위해 지원되는 튜브 유형을 수용할 수 있는 다양한 튜브 러 너가 있습니다.

- 13x75mm BD 튜브 로테이터 러너, 24개의 튜브 위치
- 13x100mm BD 튜브 로테이터 러너, 24개의 튜브 위치
- 13x75mm Greiner 튜브 로테이터 러너, 24개 튜브 위치
- 13x100mm Greiner 튜브 로테이터 러너, 24개 튜브 위치
- 16x100mm 튜브 로테이터 러너, 24개 튜브 위치



다중 분사는 매끄러운 풀 캡 및 안전 트위스트 캡을 사용하는 Greiner Vacuette® 튜브에서만 지원됩니다.

표 5: 튜브 및 러너 호환성

제품 라인	튜브			러너	
	직경 [mm]	길이 [mm]	실험 기구 정의	호환 러너	브리지 색상
Greiner Vacuette	13	100	13x100mm Greiner Vacuette(격막 포함)	1x24 13x100mm Greiner 튜 브 로테이터 러너	회색
	13	75	13x75mm Greiner Vacuette(격막 포함)	1x24 13x75mm Greiner 튜브 로테이터 러너	
	16	100	16x100mm Greiner Vacuette(격막 포함)	16x100mm 튜브 로테이터 러 너 24개	검은색
BD Vacutainer	13	100	13x100mm BD Vacutainer(격막 포함)	1x24 13x100mm BD 튜브 로 테이터 러너	흰색
	13	75	13x75mm BD Vacutainer(격막 포함)	1x24 13x75mm BD 튜브 로 테이터 러너	
	16	100	16x100mm BD Vacutainer(격막 포함)	16x100mm 튜브 로테이터 러 너 24개	검은색



4.6.3.3 피어싱 팁 보호



그림 34: 피어싱 팁 보호

피어싱 팁 보호캡은 팁 교체 및 문제 해결 과정에서 피어싱 팁의 날카로운 끝부분 을 보호하는 데 사용됩니다. 사용자가 부상을 입거나 팁이 손상되지 않도록 보호합 니다.



피어싱 팁 보호는 일회용입니다. 사용 후에는 모든 피어싱 팁 보호 장치를 생물학 적 폐기물 용기에 폐기해야 합니다.



4.6.3.4 피어싱 팁 제거 도구



그림 35: 피어싱 팁 제거 도구

피어싱 팁 제거 도구는 소프트웨어 명령으로 수축될 수 없는 튜브에 끼인 피어싱 팁 수축에 사용됩니다.

4.6.4 Frida Reader



그림 36: Frida Reader



Frida Reader는 핵산의 자동 정량화 및 정규화를 위해 고안되었습니다. Frida 리더 로 측정할 검체는 검체 증발로 인해 측정 결과가 손상되지 않도록 4°C에서 냉각해 야 합니다.

▲ 주의

진동으로 인해 잘못된 결과가 발생할 수 있습니다!

검체가 떨어질 때 진동이 발생하면 측정 결과가 부정확해지고 환자 검체의 안전이 나 임상 상태가 손상될 수 있습니다.

- 적절한 설치 위치를 위해서는 바닥이 안정되어 있어야 합니다.
- Frida Reader 측정 중에는 내부 또는 외부 진동원이 근처에 있어서는 안 됩니다.
- 저주파수와 관련된 출처를 피하십시오. 특히, 36Hz(2160rpm) 및 약 42Hz(2520rpm)의 진동은 매달려 있는 드롭의 공명 주파수이기 때문에 피해야 합니다.

▲ 주의

실내 조명으로 인해 잘못된 결과가 발생할 수 있습니다!

모듈 위의 실내 조명은 측정을 방해하고 잘못된 측정 결과를 초래하며 환자 시료의 안전 또는 임상 상태를 손상시킬 수 있습니다.

 로봇 시스템은 Frida 리더의 측정 위치에서 주변광을 차단하기 위해 전면 및 후 면 패널 상단에 불투명 커버가 있어야 합니다.

4.6.5 위상 분리기

위상 분리기는 서로 다른 점도의 액체 사이의 분리 위상을 감지하도록 설계되었습 니다. 따라서 실험 기구 외부에서 볼 수 있는 분리 위상과 독립됩니다. 위상 분리기 는 소스에서 대상 실험 기구로 액체 위상을 깨끗하게 이송해야 하는 작업에 사용할 수 있습니다.

다음은 일반적인 워크플로를 나타냅니다.

- ✓ 액체 혼합물을 담은 소스 실험 기구의 원심 분리로 액체 사이의 개별 위상을 생 성합니다.
- ✔ 원심 분리 중 위상이 형성되도록 액체의 점도가 달라야 합니다.
- 마개를 제거한 튜브를 Fluent 데크에 로드하고 완전한 추적을 위해 바코드 스 캐너(예: Fluent ID)를 사용합니다. 로드하는 동안 위상/액체 부분 사이의 층을 흩뜨리지 않도록 주의하십시오.
- 분수 분리를 위해 정의된 프로토콜을 시작합니다. Air FCA의 위상 분리기 기능 은 액체 사이의 위상을 감지하고 대상 실험 기구로 관심 부분의 이송을 시작합 니다.

소스 실험 기구에서 둘 이상의 위상을 추출할 수 있습니다. 위상 감지 중에 일회용 팁 위 피펫팅 채널의 잠재적 오염을 방지하기 위해 데크에 액체가 유출되지 않도록 (실험 기구에서 넘침) 상부 위상의 일부를 제거해야 합니다.

기술 사양은 참조 설명서를 참조하십시오. 소프트웨어에 대한 자세한 내용은 설명 서를 참조하십시오. "참조 문서" [▶ 10]를 참조하십시오.

액체 단계들은 명확하게 구분되어야 합니다. 전혈을 혈장 및 혈액 세포를 분리하는 경우 검체의 품질과 전처리가 필수적입니다. 혈액 검체 내 위상 감지에 영향을 미 칠 수 있는 매개 변수는 검체 품질(지방혈증, 용혈), 보관 시간, 보관 온도, 이송 조 건, 원심 분리 조건(시간, rcf, 온도, 램프, 로터 유형), 원심 분리 후 위상 왜곡 등



입니다.

위상 분리 결과를 최적화하려면 혈액 검체는 추출 후 가능한 한 빨리 처리해야 합니다. 검체 처리 및 보관 조건은 특정 튜브 제조업체의 권장 사항을 따라야 합니다.

속도 저하 램프를 사용해 실온에서 10분 동안 2500rcf로 원심 분리할 경우 혈장 검 체가 깨끗하게 위상 분리됩니다(Tecan 내부 검사 조건).

흡인 및 위상 감지 중에 튜브에서 유출을 방지하려면 튜브를 림까지 채우지 마십시 오. 위상 감지는 일반적으로 흡입 속도가 느린 상태에서 튜브에서 아래쪽으로 빠르 게 이동해야 합니다. 그러면 감지 중에 액체 레벨이 증가하게 됩니다. 보충 용량이 많은 튜브의 경우 위상 감지를 시작하기 전에 위쪽에서 액체를 덜어내 는 것이 좋습니다.

4.6.6 Resolvex i300



그림 37: Resolvex i300 작업대 모듈

Resolvex i300은 플랫폼에 통합할 수 있는 모듈입니다. 플랫폼은 동일한 소프트웨 어 프레임워크 내 Resolvex i300 모듈에서 작업대로부터 직접 액체를 처리하고 작 업대를 이송할 수 있습니다.

Resolvex i300은 다음과 같은 주요 기능을 사용하여 다양한 실험실 프로세스를 자 동화할 수 있는 실험실 기기입니다.

 공기 또는 질소를 사용하여 전체적으로 또는 부분적으로 로드된 필터 컬럼 가 압



- 통합된 실험 도구 리프트를 사용하여 수십 플레이트에 필터 플레이트 스택 및 언스택
- 통합 드립 보호 장치를 사용하여 필터 및 수집 실험 기구의 스택 및 언스택 시 검체를 교차 오염으로부터 보호
- 필터 또는 수집 실험 기구에 액체 분주(선택 사항)
- 가열 가스(예: 공기 또는 질소 등)에 의해 실험실 용구에서 액체 증발을 일으키 는 방법입니다.

Resolvex i300의 일반적인 애플리케이션 워크플로에는 다음과 같은 실험실 프로세 스가 포함됩니다.

- 질량 분석 검체 전처리 워크플로를 위한 고체상 추출 프로세스(SPE)
- 액체 지원 추출(SLE)
- 양압을 사용하는 기타 여과, 정제 또는 농축 워크플로

Resolvex i300을 사용하기 전에 자세한 내용을 알아보려면 Resolvex i300 작동 설명서를 참조하십시오. "참조 문서" [▶ 10]를 참조하십시오.

•TECAN•

5 제어 요소

5.1 작동 요소



A 클램프 레버 B 터치스크린 클램프 레버로 세그먼트를 잠그고 잠금 해제합니다.

운영자가 기기를 제어할 수 있도록 터치스크린에 검사법 및 설명이 표시됩니다.



5.2 사용자 인터페이스



А	탐색 경로	В	작업 영역
С	표시/옵션/작업 버튼		

FluentControl 의 사용자 인터페이스를 통해 사용자는 작동 및 시스템 관리를 위한 검사법 실행에 액세스할 수 있습니다.

5.2.1 탐색 경로

탐색 경로를 사용하면 FluentControl의 계층 구조를 이해하고 탐색할 수 있습니다.

표 6: 탐색 경로 버튼

버튼	이름	기능
$ \Delta\rangle$	· 空 日	버튼을 누르면 홈페이지 로 돌아갑니다.
System Care 🛞	탐색 창	현재 및 이전 선택 항목을 표시합니다.
<	메뉴 확장	메뉴 확장 버튼을 누르면 조명 컨트롤과 같은 옵션 이 표시되고 운영자를 전 환할 수 있습니다.

5.2.2 작업 영역

사용자 인터페이스의 작업 영역을 통해 검사법 및 설명에 액세스합니다. 검사법 실행 상태에 대한 세부 정보도 여기에 표시됩니다.



표 7: 작업 영역 버튼

버튼	이름	기능
	실행	누르면 선택한 검사법을 시작합니다.
+	추가	누르면 빠른 시작 목록에 검사법을 추가합니다.
Assay 1 unknown	선택된 검사법	'계속(Continue)'을 누를 때 실행되는 현재 선택된 검사법입니다.
Assay 3 unknown	사용 가능한 검사법	클릭하여 선택할 수 있는 검사법입니다.
Start of Day Due: is not executed	빠른 시작 버튼	누르면 선택한 검사법을 즉시 시작합니다.

표 8: 작업 영역 디스플레이

디스플레이	디스플레이 기능
00:16:52 10:09 10:43 Vot Time Tot Time	검사법 실행의 상태 및 남은 시간을 표 시합니다.
Assay 1 is ready to be started.	현재 선택된 검사법에 대한 설명 또는 현재 작업에 대한 추가 정보입니다.

5.2.3 디스플레이, 옵션 및 동작 버튼

표 9: 디스플레이, 옵션 및 동작 버튼

버튼	이름	기능
✓ Ok	확인	누르면 확인합니다.
× Cancel	취소	누르면 취소합니다.
Continue	계속	누르면 계속 진행합니다.



버튼	이름	기능
II Pause	일시 중지	누르면 현재 작업이 끝날 때 실행 일시 중지를 요청합니다.
Stop	정지	누르면 현재 작업 중이어도 실행이 즉 시 중단됩니다. 가능한 경우 시스템은 실행을 복원하거 나 계속할 수 있습니다.
THE Remove	제거	누르면 빠른 시작 보기에서 검사법을 제거합니다.
View Mode	보기 모드	누르면 목록 보기와 빠른 시작 보기 간 을 전환합니다.
Sort by	정렬 기준	누르면 검사법 실행 표시를 알파벳 순 정렬과 최근 항목 정렬 사이에서 전환 합니다.

5.2.4 검사법 복구 버튼

표 10: 디스플레이, 옵션 및 동작 버튼

버튼	이름	기능
Discard	폐기	누르면 복구된 검사법 상태를 폐기합니 다.
\rightarrow	_	누르면 다음 화면으로 이동합니다.
Recovery Point	복구 지점	누르면 이전 화면("복구 지점")으로 돌 아갑니다.
Run Recovery	복구 실행	누르면 실행을 계속합니다.



5.2.5 DeckCheck 버튼

표 11: DeckCheck 버튼

버튼	이름	기능
Left	왼쪽 카메라	왼쪽 카메라에서 촬영한 카메라 사진을 표시합니다(Fluent 780/1080 전용). 이 카메라와 관련하여 레이아웃 불일치가 발견되면 아이콘에 느낌표가 표시됩니 다.
Center	중앙 카메라	중앙 전망 카메라에서 촬영한 사진을 표시합니다. 이 카메라와 관련하여 레 이아웃 불일치가 발견되면 아이콘에 느 낌표가 표시됩니다.
Right	오른쪽 카메라	오른쪽 카메라에서 촬영한 카메라 사진 을 표시합니다(Fluent 780/1080 전용). 이 카메라와 관련하여 레이아웃 불일치 가 발견되면 아이콘에 느낌표가 표시됩 니다.
Pause Alternate	일시 중지 전환	화면에서 참조 사진과 실시간 사진을 번갈아 표시합니다. 참조 사진 또는 실시간 사진을 표시할 때 해당 사진을 정적 상태로 유지하려 면 이 버튼을 누릅니다.
Resume Alternate	재개 전환	정적 사진: 버튼을 누르면 참조 사진과 실시간 사 진을 번갈아 재개합니다.
Check	검토	시스템 재검토(예: 일부 수정이 이루어 진 경우)를 활성화합니다. 도어가 닫힙 니다. 3개 암 시스템의 경우 가운데 암이 이 동해야 합니다. 도어가 닫히지 않으면 점검이 수행되지 만 가운데 암이 카메라 시야를 차단합 니다.
Ignore & Continue _{al}	무시하고 계속하기	검사법에서 해당 명령에 대해 구성된 경우에만 나타납니다. 강조 표시된 불 일치를 무시할 수 있으며 스크립트 실 행이 계속됩니다.
Continue	계속	모든 불일치가 해결되거나 시스템에서 불일치를 찾지 않고 명령에 show always(항상 표시) 옵션이 선택된 경우 표시됩니다. 이렇게 하면 시스템이 인 식하지 못한 미묘한 색상 변화를 육안 으로 확인할 수 있습니다.



표 12: 디스플레이

디스플레이	설명	기능
	참조 사진	참조 사진은 원하는 데크 레이아웃을 표시하는 스크립트 명령에 저장됩니다.
LIVE	실시간 사진	스크립트를 실행하는 동안 카메라에서 촬영한 라이브 사진입니다.
	불일치(참조 사진과의 차이)	빨간색 사각형은 참조와 실시간 그림 사이에 불일치가 발견된 영역을 표시합 니다. 표시된 영역에는 두 가지 이상의 오류가 포함될 수 있습니다.

5.3 오류 신호 및 기기 상태



그림 40: 상태 램프

A 전원 상태 램프

B 상단 상태 램프

6

상태 램프는 다양한 색상, 지속 또는 깜박이는 표시등을 통해 기기 상태를 나타냅 니다. 상단 상태 램프는 소프트웨어가 실행 중일 때만 활성화됩니다.



표 13: 상태 표시등의 광 신호

신호	색상	모드	기기 상태
	_	꺼짐	기기의 스위치가 꺼져 있습니다(전원 공급 장치에서 연결 해제됨).
	흰색	"하트비트"	기기의 스위치가 켜져 있습니다(제어 소프트웨어 연결됨. 모듈이 아직 초기 화되지 않음).
	흰색(전원 램 프만 해당)	연속	기기 "전원 켜짐" 상태(제어 소프트웨 어가 연결되지 않음).
	FluentControl 사용자 인터페 이스의 색 구 성표	"하 <u>트비트</u> "	유휴 모드 모든 모듈이 초기화되었으며 기기가 검 사법을 실행할 준비가 되었습니다. 유휴 모드에서 약 1시간 후, 기기는 대 기 모드로 전환됩니다. 보류 모드 모든 축이 제동됩니다. 암이 ZeroG에 있지 않으며 수동으로 움직일 수 없습 니다. ZeroG 모드를 사용하려면 Move Tool(이동 도구)를 선택하도록 핵심 운 영자에게 요청하여 기기를 활성화할 수 있습니다.
	노란색	연속	티치 모드 기기가 위치를 "학습"합니다. 이 모드에서 사용자는 로봇 암을 수동 으로 움직일 수 있습니다.
	녹색	연속	검사법(스크립트 또는 프로세스)이 실 행 중입니다. 이것은 일반 "프로덕션" 모드입니다.
	빨간색	깜박임	오류 상태 제어 컴퓨터 화면 또는 터치스크린에 오류 메시지가 표시됩니다.
	색상은 사용자 가 구성할 수 있음	깜박임	사용자 프롬프트 시스템이 사용자 상호 작용을 기다리고 있습니다.
	녹색	깜박임	활성 정지 런타임 컨트롤러에 의해 또는 안전 패 널을 열어서 트리거된 의도적인 일시 정지입니다. 사용자가 데크와 상호 작용할 수 있도 록 기기가 일시 정지합니다. 운영자는 검사법을 재개할 수 있습니다.



5.4 Fluent ID 상태 LED



그림 41: Fluent ID LED Fluent ID LED는 다음 상태를 나타냅니다.

표 14: Fluent ID LED

신호	색상	모드	기기 상태
	_	꺼짐	Fluent ID가 유휴 상태입니다.
	흰색	연속	Fluent ID 전원이 켜졌습니다(아직 초기 화되지 않음).
	파란색 또는 사용자 정의 색상	깜박임	러너 로딩 또는 언로딩 준비가 완료되 었습니다.
	녹색	연속	바코드가 성공적으로 스캔되었습니다. 러너를 감독합니다. 실행이 중단되므로 언로드하지 마십시오.
	빨간색	깜박임	오류 상태 오류 메시지 및 필요한 조치가 터치스 크린에 표시됩니다.
•TECAN•

6 작동

6.1 이 장의 안전 지침

▲ 주의

잘못된 결과 또는 기기 오염!

설치 적합성 평가 및 작동 적격성 평가를 수행하지 않았거나 본 설명서에 제공된 작동 절차를 따르지 않을 경우 잘못된 결과 또는 장비 오염이 발생할 수 있습니다.

- 설치 자격 평가와 운영 자격 평가의 기록은 모두 존재하며 알려져 있습니다.
- 피펫팅 매개변수를 포함한 검사법 및 프로세스는 핵심 운영자가 검증해야 합니다.
- FCA 및 에어 FCA용 피어싱 작업에서 액체 레벨 감지는 핵심 작업자가 검증해 야 합니다.
- 피어싱 팁과 튜브 로테이터 및 농축 부식성 오염 제거 용액을 사용하는 전혈 작 업의 경우에는 커넥터의 막힘과 마모를 방지하기 위해 3개월마다 세척 스테이 션 커넥터(세척 스테이션과 시스템 액체 용기에 있음)를 모두 교체할 것을 권장 합니다. 배액 또는 세척 스테이션을 반드시 물과 같은 중성 액체로 헹구기 전에 부식성 액체를 중화해야 합니다.
- 운영자는 작동 절차, 검사법 및 프로세스에 대한 교육을 받아야 합니다.



▲ 주의

사용자의 생물학적 및 화학적 오염!

FCA 그리퍼 핑거가 손상되면 플레이트가 손실될 수 있습니다. 떨어뜨린 플레이트 는 위험 물질에 의한 오염을 유발할 수 있습니다.

• 충돌 후 FCA 그리퍼 핑거를 점검합니다.

▲ 주의

날카로운 모서리 및 지점!

Fluent Mix & Pierce 워크스테이션의 피어싱 팁에는 부상을 초래할 수 있는 뾰족한 팁과 날카로운 모서리가 있습니다.

- 기기를 로드할 때 소프트웨어 명령을 사용해 FCA를 저장 위치로 이동합니다.
- 오류가 발생한 후에는 피어싱 팁을 피어싱 팁 보호 장치로 덮고 FCA를 수동으로 저장 위치로 이동합니다. "피어싱 팁 보호" [▶ 60] 섹션을 참조하십시오.

▲ 주의

시스템의 생물학적 오염!

Fluent Mix & Pierce 워크스테이션에서 혈액이 튜브 캡을 오염시킬 수 있습니다.

- 튜브를 주의하여 취급하십시오.
- · 보호 장비를 착용하십시오.

공지사항

부식성 액체로 인한 오작동!

세척 스테이션 및 튜브를 통해 2% 표백제와 같은 부식성 농축 액체가 배출되면 세 척 스테이션 커넥터가 오작동할 수 있습니다.

 배액 또는 세척 스테이션을 반드시 물과 같은 중성 액체로 헹구기 전에 부식성 액체를 중화해야 합니다.

6.2 작동 모드

Fluent는 다음 세 가지 작동 모드로 실행할 수 있습니다.

운영자 일반 작동 모드

- 애플리케이션 또는 일상적인 시스템 관리 작업이 실행되는 정상 작동 모드입니 다.
- Fluent FluentControl 소프트웨어 런타임 컨트롤러가 모니터링합니다.

핵심 운영자 검사법 정의 모드

• 이 작동 모드는 검사법을 설정하는 조정과 같은 특수 작업을 수행하는 데 사용 됩니다.



FSE

서비스 모드

- 이 작동 모드는 작동 준비 상태를 확인하기 위해 테스트와 같은 특수 작업을 수 행하는 데 사용됩니다.
- 이 모드를 실행하려면 서비스 인증서가 필요합니다.



▲ 주의

암이 작업대에 있는 물체와 충돌합니다.

유연한 암은 수동으로 움직일 수 있습니다. 암이 고정된 상태에서 암이 수동으로 원활하게 움직이는지 확인합니다. 암 범위 기계적 정지 위치를 포함하여 암을 단단 한 물체에 충돌하지 않도록 합니다.

6.3 작동 준비

6.3.1 기기 켜기

기기를 켜려면 다음과 같은 절차를 따르십시오.

1. 외부 전원 공급 장치 뒷면에 있는 전원 스위치(A)의 전원을 켭니다.



기기의 전원이 켜지면, 전원 램프가 파란색으로 켜집니다. 섹션을 참조하십시 오.

상태 표시등이 켜지지 않으면 PC를 시작하거나 핵심 운영자에게 문의하십시 오.

2. FluentControl 소프트웨어를 시작합니다. "FluentControl 시작" [▶ 76] 섹션을 참조하십시오.



6.3.2 FluentControl 시작

- ✔ 운영 절차가 있어야 하며 알 수 있어야 합니다.
- ✔ 설치 자격 평가와 운영 자격 평가의 기록은 모두 존재하며 알려져 있습니다.
- ✔ 시스템 관리가 수행되었습니다.
- ✔ 기기의 전원이 켜져 있습니다.
- Start(시작) > All programs(프로그램) > Tecan > FluentControl을 통하여 소 프트웨어를 실행합니다.
 몇 초 후 시작 화면이 나타납니다.

6.3.3 사용자 로그인



사용자 관리는 Fluent GX Assurance Software 옵션에서만 사용할 수 있습니다.

FluentControl에 로그인하려면 다음과 같은 절차를 따르십시오.

- ✔ Fluent GX Assurance 소프트웨어를 설치합니다.
- ✔ FluentControl을 시작합니다.
- ✔ 사용자 관리는 FluentControl에서 활성화하고 프로세스를 정의합니다.
- 1. 할당된 사용자 프로필(A)을 선택합니다.



2. 키보드(B)에 암호를 입력합니다.



3. **OK**(C)를 누릅니다.



로그인 후 기기가 자동으로 초기화됩니다.

6.3.4 세그먼트 배치

세그먼트를 배치하려면 다음과 같은 절차를 따르십시오.

- ✓ 선택한 검사법에 따라 모든 세그먼트, 캐리어, 옵션 및 기기를 배치해야 합니 다.
- ✔ 세그먼트가 깨끗하고 완벽한 상태입니다.
- ✔ 세그먼트는 해당 그리드 위치에 배치됩니다.
- 1. MCA 384 활성 캐리어를 설치할 때 케이블을 연결합니다.



2. 데크의 후면으로 세그먼트를 내려놓으십시오.



- 3. 후면 가장자리는 후면 채널 덮개 또는 기기 익스텐션과 일치시키십시오.
- 4. 세그먼트(A)의 전면 부분을 조심스럽게 내립니다.



5. 클램프 레버를 왼쪽에서 오른쪽으로, 즉 닫힘 위치로 돌립니다. "세그먼트 확 인" [▶ 181] 섹션을 참조하십시오.

6.3.5 세그먼트 제거



Fluent ID 세그먼트는 제거할 수 없습니다! 기기 전자 장치에 직접 연결됩니다. 전 자 장치에 대한 연결은 자격을 갖춘 FSE에서만 수행할 수 있습니다. 고객 지원 섹션을 참조하십시오.

세그먼트를 제거하려면 다음과 같은 절차를 따르십시오.

- ✔ 모든 시약, 검체, 랙, 러너 및 플레이트가 세그먼트에서 제거되었습니다.
- ✔ 세그먼트에 아무것도 배치되지 않습니다.
- 1. 클램프 레버(A)를 오른쪽에서 왼쪽으로 돌려서 엽니다.





세그먼트가 잠금 해제되고 클램프 레버에 노란색 표시가 나타납니다.

2. 세그먼트를 약 4 mm 앞으로 밉니다.





3. 전면의 세그먼트(B)를 들어 올립니다.





4. 손상을 방지하기 위해 세그먼트를 깨끗하고 건조한 장소에 보관합니다.

6.3.6 표준 러너 로드



- 데크에 수평으로 러너를 맞춥니다.
- 한 손으로 러너의 앞쪽 끝을 지지합니다.
- 언로드 중에 러너를 들기 전에 러너가 모든 핀을 해제하는지 확인합니다.

러너를 로드하려면 다음과 같은 절차를 따르십시오.



1. 러너를 해당 그리드 위치(A)에 정렬하십시오.



- 2. 러너를 정지 위치로 밉니다.
- 러너가 세그먼트가 단단히 잠겼는지 확인합니다.
 러너가 정지 위치에 닿기 전에 마지막 몇 밀리미터에서 느껴질 수 있습니다.



러너를 언로드하려면 다음과 같은 절차를 따르십시오.

- 로딩 영역에서 완전히 제거될 때까지 평평한 데크에서 러너를 수평으로 당깁니 다.
- 2. 한 손으로 러너의 앞쪽 끝을 지지합니다.
- 3. 러너를 들기 전에 러너가 모든 핀을 해제하는지 확인합니다.



6.3.7 데크 레이아웃 확인 중

데크에 설치된 캐리어, 실험 기구, 장치가 검사법에 대해 정의된 데크 레이아웃과 일치하는지 확인합니다.

공지사항

장비 손상!

작업대에서 잘못된 세그먼트 및 실험 기구 위치로 인해 암이 충돌할 할 수 있습니다.

- 물리적 데크 구성 및 로드된 실험 기구가 FluentControl 작업대 구성과 일치하 는지 항상 확인하십시오.
- 항상 실험 기구가 네스트에 올바르게 장착되었는지 확인하십시오. "실험 기구 배치" [▶ 156]를 참조하십시오.

공지사항

자기장은 간섭을 유발합니다!

흡입 위치의 강한 자기장(위쪽: 북극)이 팁 감지 센서를 방해할 수 있으며 예상치 못한 오류(예: DiTi 손실)로 이어질 수 있습니다.

흡입 위치 근처의 SBS 위치에 강한 자석을 두지 마십시오.



Fluent ID 세그먼트는 데크 아래의 전자 보드에 연결되어 있기 때문에 FSE에 의해 서만 제거할 수 있습니다.

- ✔ 검사법은 핵심 운영자가 준비해야 합니다.
- ✔ 소모품은 이 검사법에서 정의된 소모품과 일치합니다.
- ✓ Fluent ID 튜브 러너는 터치스크린에 메시지가 표시될 때 검사법이 시작된 후에 만 로드되어야 합니다.
- 1. 터치스크린에 표시되는 지침을 따르십시오.





그림에는 터치스크린에 표시되는 지침(A)의 예가 나와 있습니다.

6.4 검사법 시작 전

검사법을 시작하기 전에 다음 체크리스트를 작성해야 합니다.

표 15: 검사법 시작 전 확인 사항

기기/구성요소	작업	참고/활동
공정 밸리데이션	프로덕션 실행을 시작하 기 전에 선택하려는 검사 법이 검증되었는지 확인 합니다.	자세한 내용은 핵심 운영 자에게 문의하십시오.
터치스크린	터치스크린의 지침을 따 르십시오. 공지사항! 핵심 운영자가 제공하는 지침을 엄격히 준수해야 합니다. 지침이 표시되지 않는 경 우 아래 작업 목록을 따릅 니다.	_
세그먼트, 캐리어, 옵션 및 장치	모든 세그먼트, 캐리어, 옵션 및 장치가 설치되고 고정되었는지 확인합니 다. 검사법에서 사용할 물체 만 데크에 있는지 확인하 십시오. 테스트 실행이 성공적으 로 완료되었는지 확인합 니다.	검사 실행이 실패할 경우, 핵심 운영자에게 연락하 여 검사 실행을 다시 수행 합니다.



기기/구성요소	작업	참고/활동
검체 및 시약	모든 검체, 시약 및 실험 기구가 올바르게 장착되 었는지 확인합니다. 공지사항! 바코드 스캐닝 은 검사법이 시작된 후에 만 발생합니다. 이 방법을 시작하기 전에 Fluent ID 데크에 러너가 없는지 확 인하십시오. 러너는 터치 스크린에 메시지가 표시 될 때만 로드되어야 합니 다.	_
폐기물 튜브(액체 시스템 만 해당)	폐기물 튜브가 올바로 배 치되었는지 확인합니다.	폐기물 튜브가 꼬이거나 구겨지지 않았는지 육안 으로 검사합니다. 결함이 있는 폐기물 배관 을 교체합니다. "세척 스 테이션 연결(MCA 384)" [▶ 155] 섹션을 참조하십 시오.
세척 시스템(액체 시스템 만 해당)	시스템 액체 및 폐기물 용 기가 올바르게 연결되어 있는지 확인합니다.	"시스템 액체 용기 및 폐 기물 용기의 튜브 점검" [▶ 85] 섹션을 참조하십 시오.
세척 시스템(액체 시스템 만 해당)	시스템 액체 용기가 올바 른 레벨로 채워져 있는지 확인합니다. 폐기물 용기가 비어 있는 지 확인합니다.	"시스템 액체 용기 및 폐 기물 용기 연결" [▶ 134] 섹션을 참조하십시오.
세척 시스템(액체 시스템 만 해당)	검사법에 정의된 대로 올 바른 시스템 액체를 사용 해야 합니다.	-
세척 시스템(MCA 세척 센터만 해당)	세척 블록의 액체 수위를 확인합니다.	-
일회용 팁 폐기물 처리장 치 및 세척 기구	일회용 팁 폐기물과 세척 스테이션 장치가 깨끗한 지 확인합니다.	"일회용 팁 폐기물 및 세 척 스테이션 유닛 세척" [▶ 128] 섹션을 참조하십 시오.
	MCA 384 팁 유도 또는 에어로졸 격리를 위해 폐 기물 슈트 커버가 장착되 었는지 확인합니다.	_



기기/구성요소	작업	참고/활동
일회용 팁	올바른 팁이 장착되어 있 는지 확인하십시오. 팁 폐기물이 비어 있는지 확인합니다.	_
고정 팁	고정 팁이 깨끗하고 손상 이 없는지 확인합니다.	고정 팁을 육안으로 검사 하여 깨끗한지 확인합니 다. 치과 거울로 고정된 끝부 분을 시각적으로 검사하 여 코팅이 완전히 유지되 는지 확인하십시오.
데크	데크에 설치된 캐리어, 실 험 기구, 장치가 검사법에 대해 정의된 데크 레이아 웃과 일치하는지 확인합 니다.	"데크 레이아웃 확인 중" [▶ 82] 섹션을 참조하십시 오.
실험 기구	모든 실험 기구가 단단히 배치되었는지 확인합니 다. 마이크로플레이트가 옆으로 이동하는 경우 실 험 기구 포지셔너 포지셔 너가 올바른지 확인합니 다.	"실험 기구 배치" [▶ 156] 섹션을 참조하십시오.
튜브 로테이터	튜브 로테이터에 위치 지 정 핀 또는 잠금 핀이 없 는지 확인합니다.	"잠금 핀 및 위치 지정 핀 교체" [▶ 183] 섹션을 참조 하십시오.

6.4.1 시스템 액체 용기 및 폐기물 용기의 튜브 점검

⚠ 주의

표본 오염!

쌍액 FCA 구성인 경우 각 암에 사용할 시스템 액체의 유형이 다를 수 있습니다. 잘 못된 시스템 액체 용기를 암에 연결하면 검체가 오염될 수 있습니다. • 각 시스템 액체 용기에 해당 시스템 액체 이름을 표시합니다.

- ✔ 세척 시스템이 올바르게 설치되어 있어야 합니다.
- 1. 튜브(B)가 시스템 액체 용기(D)에 올바르게 연결되어 있는지 확인합니다.





2. 액체 감지 센서(C)가 시스템 액체 용기(D)에 올바르게 연결되어 있는지 확인하 십시오.

6.4.2 폐기물 용기 튜브 점검

1. 튜브(A)가 폐기물 용기(D)에 연결되어 있는지 확인합니다.



- 2. 튜브(C)와 (D)가 폐기물 용기에 올바르게 연결되어 있는지 확인합니다.
- 액체 감지 센서(E)가 있는 경우 폐기물 용기에 올바르게 연결되어 있는지 확인 합니다.



4. 뚜껑을 돌려 닫습니다(F).



6.5 검사법 실행하기

방법은 FluentControl 소프트웨어에서 정의된 스크립트 또는 프로세스의 모음입니 다. 이 방법은 즉시 실행할 수 있습니다.

핵심 운영자는 다음과 같이 실행할 수 있는 검사법을 작성합니다.

공지사항

기구 손상!

데크가 제대로 설정되지 않거나 소프트웨어가 잘못 작동하거나 잘못 사용될 경우 기기 손상이 발생할 수 있습니다.

- 모든 안전 장치가 설치되어 작동하고 있는지 확인합니다.
- 데크에 설치된 캐리어, 실험 기구, 장치가 검사법에 대해 정의된 데크 레이아웃 과 일치하는지 확인합니다.
- 검사법에서 사용할 물체만 데크에 있는지 확인하십시오.

6.5.1 검사법 시작

- ✓ FluentControl이 시작되었습니다.
- ✔ "사용자 로그인" [▶ 76] 섹션이 수행되었습니다.
- ✔ "검사법 시작 전" [▶ 83] 섹션이 수행되었습니다.
- 1. **Method Starter**(검사법 스타터)(A)를 선택합니다.



버튼을 누르면 바로 버튼이 켜집니다.



- 2. 실행할 검사법(B)을 선택합니다. *선택한 검사법이 강조 표시됩니다.*
- 3. **OK**(확인)를 누릅니다.





4. **Run**(실행)(C)을 누릅니다.



- 5. 터치스크린의 지침을 따르십시오.
- 6. 스크립트에 DeckCheck가 포함된 경우 예상 참조 데크 레이아웃과 실제 실시 간 데크 레이아웃의 차이에 주의하십시오. "DeckCheck 작업" [▶ 99]을 참조하 십시오.
- 검사법 실행이 끝날 때까지 기다립니다.
 화면에는 실행 중인 검사법이 종료하는 시점의 예상 시간(D)이 표시됩니다.







실행을 완료했으며 오류가 발생했습니다! 메시지가 나타나면 Log(로그)를 눌러 오류 및 경고를 검토합니다.

9. Next(다음)를 눌러 홈 화면으로 돌아갑니다.

11.5	Run finished with	erro	or(s)!
Start Tir	a Ne		End Tim
6	omment Process	Log	
1	E: Script execution terminated by End Script statement. Run paused.		12:00:48 (2017-08-18) 11:59:35 (2017-08-18)
=			

10 Exit(종료)를 누릅니다.

.



검사법 실행이 종료되면 아래 화면(E)이 표시됩니다.



6.5.2 Fluent ID 러너 로드 및 언로드



Fluent 제품은 레이저 방사선을 방출하는 IEC 60825-1:2014에 따른 클래 스 1 레이저 제품입니다.

레이저 빔으로 인해 눈부심, 섬광맹, 잔상이 발생할 수 있습니다.

- 레이저 빔이나 반사된 빔을 응시하지 마십시오.
- 6.5.2.1 Fluent ID 러너 로드

⚠ 주의

공지사항

부적절한 로드 또는 언로드로 인한 손상

러너 및 핀 손상

- 데크에 수평으로 러너를 맞춥니다.
- 한 손으로 러너의 앞쪽 끝을 지지합니다.
- 언로드 중에 러너를 들기 전에 러너가 모든 핀을 해제하는지 확인합니다.
- ✔ Fluent 에는 Fluent ID 튜브 바코드 스캐너가 장착되어 있습니다.
- ✔ 튜브는 바코드 라벨이 왼쪽을 향하게 하여 러너에 로드됩니다.
- ✓ 러너의 모든 튜브는 크기와 모양이 동일합니다. 튜브 러너의 경우에는 "Fluent ID 호환 튜브 러너" [▶ 55] 섹션을 참조하십시오.
- 터치스크린을 사용하여 검사법을 선택하고 시작합니다.
 LED가 깜박이기 시작하며 Please load Tubes(튜브를 로드하십시오) 메시지가 터치스크린에 표시됩니다.
 서로 다른 튜브 유형을 사용하는 경우 각 그리드에 올바른 유형의 러너를 사용 해야 합니다.
- 2. 한 손으로 러너의 앞쪽 끝을 지지합니다.



- 3. 데크의 높이에서 러너를 수평으로 잡습니다.
- 4. 러너를 정지 위치로 밉니다.



5. 튜브 러너를 차례로 Fluent ID 로딩 영역의 전용 그리드로 밉니다.



 모든 바코드가 성공적으로 스캔되었는지 확인합니다.
 튜브 러너가 로딩 위치에 있고 모든 바코드 라벨이 성공적으로 스캔되면 LED 가 녹색으로 바뀝니다.



6 - 작동 검사법 실행하기



Fluent ID LED 상태 설명은 "Fluent ID 상태 LED" [▶ 72] 섹션을 참조하십시오.

- 바코드 스캔 오류가 발생한 경우, 러너를 언로드하고 문제를 해결한 후 러너를 다시 로드하십시오.
- 8. 러너가 완전히 제거될 때까지 데크를 따라 수평으로 당깁니다.



Fluent ID는 스캐너를 통과할 때 모든 코드를 여러 번 읽습니다. 작은 튜브 및 좁은 튜브(즉, 직경 ≤10mm)의 경우 수동 로드 속도를 줄여 모든 판독 작업을 가능하게 하고 오류 보고를 줄입니다.

	0000000000000000		000000000000000000000000000000000000000	000000000000000000000000000000000000000	000000000000000000000000000000000000000	000000000000000000000000000000000000000							
	1	2	3	4	5	6							
Er	rors:		Missi	ng Ru	nner		Wrong Runne	r 🔸	Unreadable	•	Duplicates	0	Missing

그림 42: 바코드 판독 확인 메시지가 터치스크린에 표시됨

표 16: GUI 의미(러너)

사각형(러너)	의미
녹색	러너의 모든 튜브 바코드가 성공적으로 판독되었습니다.
빨간색 윤곽선이 있는 흰색	이 그리드 위치에 대한 러너 유형이 잘 못되었습니다.



사각형(러너)	의미
빨간색 윤곽선이 있는 회색	러너가 없습니다. 러너가 이 그리드 위 치에 로드되어야 합니다.

표 17: GUI 의미(튜브 위치)

원(튜브 위치)	의미
녹색	바코드를 성공적으로 읽었습니다.
빨간색	읽을 수 없는 바코드
주황색	중복 바코드
빨간색 윤곽선이 있는 흰색	튜브가 없습니다. 튜브는 이 위치에 로 드되어야 합니다.



2ml 안전 잠금 튜브 러너를 사용하는 경우 읽을 수 없는 바코드에서 누락된 튜브를 구분할 수 없습니다. 누락된 튜브는 읽을 수 없는 바코드로 보고됩니다.

6.5.2.2 Fluent ID 러너 언로드

- ✓ 실행이 완료되었거나 실행이 진행 중이며, LED가 깜박이고 동시에 터치스크린 에는 Please unload tubes(튜브를 언로드하십시오) 메시지가 표시됩니다.
- 1. 러너가 완전히 제거될 때까지 데크를 따라 수평으로 당깁니다.
- 6.5.3 튜브 로테이터 러너 로드 및 언로드
- 6.5.3.1 튜브 로테이터 러너 로드

▲ 주의

시스템 및/또는 사용자의 생물학적 오염!

손상된 검체 튜브가 파열되어 튜브 로테이터에서 검체가 유출될 수 있습니다.

- 손상된 튜브가 튜브 로테이터에 장착되어 있지 않은지 확인합니다.
- ✔ Fluent에는 튜브 로테이터가 장착되어 있습니다.
- ✔ 튜브는 바코드 라벨이 왼쪽을 향하게 하여 튜브 로테이터 러너에 로드됩니다.
- ✓ 러너의 모든 튜브는 크기와 모양이 동일합니다. 튜브 러너의 경우에는 "튜브 로 테이터 러너" [▶ 59] 섹션을 참조하십시오.
- 1. 터치스크린을 사용하여 검사법을 선택하고 시작합니다.

LED가 깜박이기 시작하며 Please load Tube(튜브를 로드하십시오) 메시지가 터치스크린에 표시됩니다. 다른 튜브 유형을 사용하는 경우 튜브 유형(BD 또는 Greiner)당 올바른 러너를 선택했는지 확인하십시오. 또한, 다양한 높이의 튜브를 해당 러너에 넣으십시 오. 튜브는 항상 러너 브리지에 의해 캡 높이의 위치로 고정됩니다. 튜브 하단 은 항상 러너의 튜브 인서트에 단단히 고정되어 있어야 합니다.



2. 러너 잠금 레버를 엽니다.



3. 한 손으로 러너의 앞쪽 끝을 지지합니다.



- 4. 데크의 높이에서 러너를 수평으로 잡습니다.
- 5. 러너를 정지 위치로 밉니다.





6. 튜브 로테이터 러너를 차례로 튜브 로테이터의 전용 그리드로 밉니다.

7. 모든 바코드가 성공적으로 스캔되었는지 확인합니다.

튜브 러너가 로딩 위치에 있고 모든 바코드 라벨이 성공적으로 스캔되면 LED 가 녹색으로 바뀝니다. 튜브 로테이터 LED 상태 설명은 "Fluent ID 상태 LED" [▶ 72] 섹션을 참조하십 시오.

 바코드 스캔 오류가 발생한 경우, 러너를 언로드하고 문제를 해결한 후 러너를 다시 로드하십시오.



9. 러너 잠금 레버를 닫습니다.





그림 43: 바코드 판독 확인 메시지가 터치스크린에 표시됨

표 18: GUI 의미(러너)

사각형(러너)	의미
녹색	러너의 모든 튜브 바코드가 성공적으로 판독되었습니다.
빨간색 윤곽선이 있는 흰색	이 그리드 위치에 대한 러너 유형이 잘 못되었습니다.
빨간색 윤곽선이 있는 회색	러너가 없습니다. 러너가 이 그리드 위 치에 로드되어야 합니다.

표 19: GUI 의미(튜브 위치)

원(튜브 위치)	의미
녹색	바코드를 성공적으로 읽었습니다.
빨간색	읽을 수 없는 바코드
주황색	중복 바코드
빨간색 윤곽선이 있는 흰색	튜브가 없습니다. 튜브는 이 위치에 로 드되어야 합니다.

6.5.3.2 튜브 로테이터 러너 언로드



튜브가 장착된 튜브 로테이터 러너를 액체 취급 작동 조건 외부에 보관하지 마십시 오. "환경 조건" [▶ 41] 섹션을 참조하십시오.

- ✓ 실행이 완료되었거나 실행이 진행 중이며, LED가 깜박이고 동시에 터치스크린 에는 Please unload tubes(튜브를 언로드하십시오) 메시지가 표시됩니다.
- ✔ 튜브 로테이터가 가로 홈 위치에 있습니다.





1. 러너 잠금 레버를 엽니다.



2. 러너가 완전히 제거될 때까지 데크를 따라 수평으로 당깁니다.



튜브 로테이터 세그먼트는 데크 아래의 전자 보드에 연결되어 있기 때문에 FSE에 의해서만 제거할 수 있습니다.

- 검사법은 핵심 운영자가 준비해야 합니다.
- 소모품은 이 검사법에서 정의된 소모품과 일치합니다.
- 튜브 로테이터 러너는 터치스크린에 메시지가 표시될 때 검사법이 시작된 후에 만 로드해야 합니다.

6.5.4 오류 재설정

메시지가 표시되면 다음과 같은 절차를 따르십시오.

- 메시지 1. 디스플레이 기능, 버튼 기능 또는 오류 메시지를 확인합니다. "작업 영역" [▶ 66] 및 "검사법 복구 버튼" [▶ 68] 섹션을 참조하십시오.
 - 2. 이 설명서 및 터치스크린의 지침에 따라 오류를 수정합니다.
 - 3. 검사법 실행을 계속합니다. "디스플레이, 옵션 및 동작 버튼" [▶ 67] 섹션을 참조 하십시오.

상태 표시등이 켜지거나 색상이 변경되면 다음과 같은 절차를 따르십시오.

- 상태 램프 1. 기기 상태를 확인하십시오. 섹션을 참조하십시오.
 - 2. Fluent에 Fluent ID 튜브 바코드 스캐너가 장착된 경우, Fluent ID 튜브 바코드 스캐너 LED 상태를 확인하십시오. "Fluent ID 상태 LED" [▶ 72] 섹션을 참조하 십시오.
 - 3. 디스플레이 기능, 버튼 기능 또는 오류 메시지를 확인합니다. "작업 영역"
 [▶ 66] 및 "검사법 복구 버튼" [▶ 68] 섹션을 참조하십시오.
 - 4. 문제 해결 표를 확인하십시오. "문제 해결 표" [> 144] 섹션을 참조하십시오.
 - 5. 문제를 해결할 수 없는 경우 "고객 지원" [▶ 188]을 참조하십시오.



6.6 DeckCheck 작업

스크립트에 DeckCheck 사용이 포함된 경우 DeckCheck 카메라 시스템은 로드 후 작업대 사진을 찍고 실제 실시간 레이아웃을 참조 레이아웃과 비교합니다.

DeckCheck에서 데크의 사진을 찍고 실시간 및 참조 레이아웃의 비교를 표시하는 데 3개 암/3개 카메라 시스템의 경우 약 20초, 1개 또는 2개 암/단일 카메라 시스 템의 경우 약 12초가 소요됩니다(이때 PC 구성이 적합하다고 가정 -FluentControl 애플리케이션 소프트웨어 설명서 참조).

기기 전원만 사용한 후 처음 사용할 경우 DeckCheck 명령이 첫 번째 결과를 표시 하는 데 시간이 오래 걸립니다. 이 작업에 몇 분 정도 걸릴 수 있습니다.

이 시간 동안 후면 LED가 켜집니다.

3개 암을 사용하는 Fluent 시스템에서 가운데 암은 왼쪽 및 오른쪽 위치 사이로 이 동해야 합니다(1개 또는 2개 암을 사용하는 시스템에서는 왼쪽 및 오른쪽 암이 각 각 왼쪽 및 오른쪽 끝에 위치함). 이 암을 이동하려면 전면 도어를 닫아야 합니다. 도어가 열려 있는 상태로 이미지를 촬영할 경우 일반적으로 가운데 암 또는 수동으 로 이동한 암에 의해 하나의 카메라 시야가 차단됩니다.

DeckCheck 프로세스 중 터치스크린에 화면을 가로질러 이동하는 형상이 표시되고 Taking Images(이미지 촬영 중) 후 Checking(확인 중)이 표시됩니다. 데크 이미지들은 장치 크기와 구성에 따라 12-20초 후에 교대로 표시됩니다. 표시된 이미지는 첫 번째 카메라에서 촬영한 사진으로, 표시된 차이는 왼쪽부터 시작됩니다.

참조 사진과 관련된 불일치를 표시하는 DeckCheck 화면. 여기서 참조 사진은 플 레이트가 있어야 하고 가운데 카메라가 불일치를 감지했음을 보여줍니다.



하단에 Ignore & Continue(무시 및 계속)가 표시되면 현재 작업 테이블을 계속 진 행합니다. 필요한 작업대와 도어를 닫기 전에 차이가 없음이 확실하면 Ignore & Continue(무시 및 계속) 버튼을 선택합니다. 데크의 새 이미지를 촬영하려면 Check(확인)를 선택합니다. 3개 암 시스템에서 도어가 닫히지 않으면 이미지가 촬 영되지만 가운데 암이 카메라를 가로막습니다. 그렇지 않으면 도어가 닫힐 때 점검 이 자동으로 실행됩니다.





스크립트에 show always(항상 표시) 옵션이 포함된 경우 위의 화면이 표시되고 불 일치가 강조 표시되지 않습니다. 그러나 참조 및 실시간 사진이 번갈아 표시될 수 있으며, 시스템에 의해 캡처되지 않지만 육안으로 쉽게 확인 가능한 작은 차이가 있을 수 있습니다(예: 일부 색상 차이, 단일 누락 튜브/팁 또는 작은 측면 이동). 아 래 나열된 한도를 참조하십시오.

불일치가 발견될 경우 해당 사항이 강조 표시됩니다.

차이를 수정하려면:

- 1. 도어를 열고 강조 표시된 항목의 위치를 교체하거나 교정합니다.
- 2. DeckCheck는 교정된 실시간 상황과 참조 레이아웃을 비교하기 위해 계속 작 동합니다.
- DeckCheck 버튼을 사용하여 각 카메라에 의해 캡처된 차이를 확인하거나 보 기를 일시 중지하고 필요한 경우 참조 사진을 유지합니다. 차이가 더 이상 감지되지 않으면 녹색 continue(계속) 버튼이 나타납니다.
- 4. Continue(계속)를 선택하여 검사법을 계속 진행합니다.



나머지 차이가 현실적으로 허용 가능한 수준이라고 판단되면(예: 검사 시작 시 총 팁 수가 다를 수 있으며 실행 시작 시 액체 레벨이 크게 다를 수 있음) 핵심 운영자 가 스크립트에 제공한 경우 Ignore & Continue(무시 및 계속)를 선택할 수 있습니 다.

DeckCheck에서 일부 레이아웃 차이를 강조 표시할 수 없습니다(예: 다음과 같은 색상의 FCA 팁 트레이).

MCA 헤드 어댑터 유형 간 차이:

- 노란색/주황색
- 흰색/주황색



• 회색/모든 색상

MCA 384 다양한 팁 유형

MCA 96 다양한 팁 유형

일부 로드된 튜브 러너에 튜브 누락

수조 300 SBS

마이크로플레이트가 180도 회전됨

마이크로플레이트 웰 형상(예: 둥근 바닥 또는 PCR 웰)

주변 호텔 내 플레이트 10ml/25ml 수저(인서트용)

투명 뚜껑

하지만, 이러한 차이점 중 상당수는 라이브 레이아웃과 참조 레이아웃을 전환할 때 명확하게 확인할 수 있습니다.

6.7 검사법 복구

FluentControl은 다음과 같은 오류 복구 옵션을 제공합니다. 예:

이전 검사법 실행이 중단되었거나 심각한 오류가 있었습니다. 검사법 복구 옵션에 서는 이전 실행에서 실패가 발생한 지점부터 다시 시작할 수 있는 가능성을 제공합 니다.



검사법이 중단되거나 치명적인 오류가 발생한 후, 일일 유지관리를 실행해야 합니 다. "일일 시스템 관리" [▶ 108]를 참조하십시오.

6.7.1 검사법 복구 모드로 전환

- ✔ 핵심 운영자가 FluentControl 에서 검사법 복구 옵션을 활성화했습니다.
- ✔ 이전 검사법 실행이 취소되었습니다.
- 1. Method Recovery(검사법 복구)(A)를 선택하십시오.





6.7.2 검사법 실행 복구

- ✔ "검사법 복구 모드로 전환" [▶ 101] 섹션이 수행되었습니다.
- 다음 화면(A)에 계속 가려면 을 선택하십시오.
 화면에 마지막으로 실행된 스크립트 줄(C) 및 오류가 발생한 스크립트 줄(복구 지점(B))이 표시됩니다.

₫〉	Method Recovery 🛞	
	Reagent Distribution	
	Sample Transfer	— C
	Get Tips	FCA 1 FCA, 200ul SBSB
	Aspirate	FCA 1 100 [µl] Water Video Single 1x24 13x75mm Tube Runner[001] Q1 - X1
	Jispense	FCA 1 100 [µ] Water Video Single 96 Well Flat[001] Al - H1
	Drop Tips	FCA.1 Wash Station Thru Deck DiTi Waste_1
	Incubate & Wash	
*=	Read & Export Data	
Ŵ		A → ×
Discard		Worktable Cancel

- 2. "검사법 복구 버튼" [▶ 68] 섹션에서 설명한 필수 버튼을 선택합니다.
- Fluent의 물리적 데크 레이아웃이 터치스크린에 표시된 작업대 레이아웃(D)과 일치하는지 확인합니다.
- Run Recovery(복구 실행)(E)를 선택합니다.
 시스템이 시작됩니다.



6.8 기기 전원 끄기

실행 중인 검사법이 없을 경우 기기는 STANDBY(대기) 모드로 전환됩니다. 주 전 원에서 기기를 끌 필요는 없습니다.



기기를 끄려면 다음과 같은 절차를 따르십시오.

- 1. 어떤 검사법이든 중지하고 터치스크린에서 STANDBY(대기) 모드를 선택합니다.
- 2. 기기가 대기 모드에 있는지 확인합니다. 섹션을 참조하십시오.
- 3. 로봇 암을 자유롭게 움직일 수 있는 공간에 배치합니다.
- 4. 외부 전원 공급 장치 뒷면에 있는 전원 스위치(A)를 사용하여 전원을 끕니다.





7 시스템 지원

이 장에서는 Fluent의 작동 상태를 양호하게 유지하기 위해 수행해야 하는 모든 시 스템 관리 작업에 대한 지침을 제공합니다.



작동 상태가 양호한 경우에만 Fluent을(를) 작동하십시오. 본 설명서에 기술된 시 스템 관리 지침을 엄격히 준수하십시오. 최적의 성능 및 신뢰성을 보장하기 위해 정기적으로 유지 관리 및 청소 작업을 수행하십시오.

문제가 있거나 질문이 있는 경우 "고객 지원" [> 188] 섹션을 참조하십시오.



암이 작업대에 있는 물체와 충돌합니다.

유연한 암은 수동으로 움직일 수 있습니다. 암이 고정된 상태에서 암이 수동으로 원활하게 움직이는지 확인합니다. 암 범위 기계적 정지 위치를 포함하여 암을 단단 한 물체에 충돌하지 않도록 합니다.

7.1 오염 제거

\land 주의

오염 제거는 표준 실험실 규정에 따라 "오염 제거 고지" [▶ 36] 섹션에 기재된 상황 에서 필요합니다.

▲경고

오염!

Fluent에 물질 잔여물이 있으면 개인에게 부상을 초래할 수 있고 공정의 무결성을 저해할 수 있습니다.

• Fluent 기기와 그 부품 및 부속품을 사용하기 전에 오염을 제거하십시오.

오염 물질 유형 및 오염 정도에 따라 핵심 운영자가 오염 제거 방법을 정의해야 합 니다. 오염 제거제 및 적용 모드 선택에 대한 지침은 이 장에 나와 있습니다.



과산화수소 증기 처리에 대한 자세한 내용은 참조 설명서를 참조하십시오. 를 참조 하십시오.

\land 주의

Frida Reader의 측정 결과가 올바르지 않습니다!

삽입물이 장착되지 않은 경우 Frida Reader가 잘못된 측정 결과를 제공할 수 있습니다.

• 인서트를 제거했을 때는(예: 세척 등) 빨간색 블라인드 플러그를 사용하십시오.



7.2 세정제

7.2.1 세정제 사양

시스템 관리를 위해 특수 세정제가 필요합니다. 모든 권장 세정제는 신중하게 선택 되고 테스트되었습니다.

공지사항

효과 및 화학적 적합성이 감소합니다!

Tecan에서 권장하는 세정제 이외의 세정제를 사용할 경우 세정제의 효과와 화학 적 호환성에 대한 보장이 없습니다.

- Tecan에서 권장하는 세정제만 사용하십시오.
- 세정제는 시스템 관리대에서 사용되는 각 특정 용도에 대해 정의됩니다. 특정 작업에 사용하도록 지정되지 않은 경우 세정제를 사용하지 마십시오.

다음 표는 시스템 관리 표와 시스템 관리 활동에 명시된 대로 사용하도록 지정된 세척제를 명시합니다.

작용물	사양
탈이온수	증류수 또는 탈이온수
알코올	70% 에탄올 또는 100% 이소프로판올 (2□프로판올)
약한 세제	Liqui-Nox
표면 활성 제제	Contrad 70, Contrad 90 / Contrad 2000, Decon 90
소독제	Bacillol 플러스, SporGon
표면 소독제 (핵산 오염용)	DNAzap
약산	황산 0.3M, 10% 아세트산, 30-40% 포름산
염기	수산화나트륨 0.1M
표백제	2% 차아염소산 나트륨
시스템 액체	검사법에 지정된 경우. 시스템이 작동 하지 않는 동안(예: 야간 또는 주말) 염 분을 포함한 수용액을 씻어내야 합니 다. 시스템 지원 "일과 종료" [▶ 109]을 참조하십시오.

표 20: 시스템 관리 표에 설명된 대로 사용하는 세척제

7.2.2 상업용 세정제

세제 제조업체가 제공하거나 본 설명서에 제공된 모든 지침을 주의하여 읽고 따라 야 합니다.



다음은 시스템 관리 표 및 시스템 관리 활동에 따라 사용할 수 있도록 지정된 상용 청소제 및 소독제의 목록입니다.

표 *21:* 상업용 세정제

세정제	세정제 범주	제조업체
DNAzap	표면 소독제(핵산으로 오 염된 표면)	Ambion www.ambion.com
Decon, Contrad	표면 활성 제제	Decon Laboratories www.deconlabs.com
SporGon	소독제	Decon Laboratories www.deconlabs.com
Bacillol Plus	소독제	www.bode-chemie.com
Liqui-Nox	약한 세제	Alconox www.alconox.com

7.3 시스템 관리 모드

핵심 운영자는 "시스템 관리 표" [▶ 107]에서 제공되는 시스템 관리 표에 따라 요구 되는 시스템 관리 방법을 정의합니다. 터치스크린에서 액세스할 수 있는 System Care(시스템 관리) 모드는 시스템 관리 작업에 대한 지침을 제공합니다.

7.3.1 시스템 관리 모드로 전환

- ✔ 시스템 관리 방법을 사용할 수 있어야 합니다.
- 1. System Care(시스템 관리) (A)를 선택합니다.





2. 실행할 작업을 선택합니다.



- 3. Run(실행)을 눌러 시스템 관리 방법을 시작합니다.
- 4. 시스템 관리 작업을 수행합니다.

7.3.2 오류 재설정

메시지가 표시되면 다음과 같은 절차를 따르십시오.

- 메시지 1. 디스플레이 기능, 버튼 기능 또는 오류 메시지를 확인합니다. "작업 영역" [▶ 66] 및 "검사법 복구 버튼" [▶ 68] 섹션을 참조하십시오.
 - 2. 이 설명서 및 터치스크린의 지침에 따라 오류를 수정합니다.
 - 3. 검사법 실행을 계속합니다. "디스플레이, 옵션 및 동작 버튼" [▶ 67] 섹션을 참조 하십시오.

상태 표시등이 켜지거나 색상이 변경되면 다음과 같은 절차를 따르십시오.

- 상태 램프 1. 기기 상태를 확인하십시오. 섹션을 참조하십시오.
 - 2. Fluent에 Fluent ID 튜브 바코드 스캐너가 장착된 경우, Fluent ID 튜브 바코드 스캐너 LED 상태를 확인하십시오. "Fluent ID 상태 LED" [▶ 72] 섹션을 참조하 십시오.
 - 3. 디스플레이 기능, 버튼 기능 또는 오류 메시지를 확인합니다. "작업 영역"
 [▶ 66] 및 "검사법 복구 버튼" [▶ 68] 섹션을 참조하십시오.
 - 4. 문제 해결 표를 확인하십시오. "문제 해결 표" [▶ 144] 섹션을 참조하십시오.
 - 5. 문제를 해결할 수 없는 경우 "고객 지원" [▶ 188]을 참조하십시오.

7.4 시스템 관리 표

최적의 성능 및 신뢰성을 보장하기 위해 권장 사항에 따라 유지 관리 및 청소 작업 을 수행하십시오.



시스템 관리 테이블의 작업은 시스템 관리 모드에서만 수행할 수 있습니다. "시스 템 관리 모드" [▶ 106] 섹션을 참조하십시오.

시스템 관리 작업은 일일, 주간, 월간 등 일정한 간격으로 수행해야 합니다.



7.4.1 일일 시스템 관리

7.4.1.1 일과 시작

핵심 운영자가 수행할 수 있는 경우 DailySystemCare 방법을 실행하거나 아래 표 에 나온 시간 순서대로 Fluent 암 구성에 해당하는 개별 작업을 수행합니다.

표 22: 하루 시작 시스템 관리 표

기기/ 구성 요소	시스템 관리 작업	세정제/ 일회용 제품/ 장치	참조/ 시스템 관리 활동
피어싱 팁	관찰하여 관절 끝부분 에 침착물이 있는지 확 인합니다. 필요한 경우 세척합니다. 팁이 구부 러지지 않았는지 확인 합니다.	70% 에탄올 또는 2% 표백제 및 보풀이 없는 천	"피어싱 팁 세척" [▶ 119] 섹션 을 참조하십시오.
일회용 팁 콘 및 고정 팁	손상 및 침전물이 있는 지 검사	_	이 작업은 일일 시스템 관리 검사법에 포함되어 있습니다. 공지사항! 일회용 팁은 재사 용할 수 없습니다.
고정 팁	세정합니다. 팁이 구부 러지지 않았는지 확인 합니다. 치과 거울로 시 각적으로 확인하여 코 팅이 완전히 유지되는 지 확인하십시오.	70% 에탄올이나 100% 이소프로판올과 보풀이 없는 천을 사용하십시 오.	"고정 팁 세척" [▶ 119] 섹션을 참조하십시오.
시스템 액체 용기 (고정 팁이 있는 액체 FCA 및 MCA)	기포가 보이지 않고 깨 끗하며 가득 찼는지 확 인합니다. 튜브-용기 커넥터가 올 바르게 연결되어 있는 지 확인합니다.	_	이 작업은 일일 시스템 관리 검사법에 포함되어 있습니다.
액상 폐기물 용기 (고정 팁이 있는 액체 FCA 및 MCA)	비어 있는지 확인합니 다. 튜브-용기 커넥터가 올 바르게 재결합되었는지 확인합니다.	_	이 작업은 일일 시스템 관리 검사법에 포함되어 있습니다.
일회용 팁 폐기물 백	비어 있는지 확인합니 다.	_	"일회용 팁 폐기물 백 교체" [▶ 131] 섹션을 참조하십시오. 이 작업은 일일 시스템 관리 검사법에 포함되어 있습니다.


기기/ 구성 요소	시스템 관리 작업	세정제/ 일회용 제품/ 장치	참조/ 시스템 관리 활동
액체 시스템 (Liquid FCA)	깨끗한지 확인합니다.	시스템용 액체, 알코올, 탈이온수	이 작업은 일일 시스템 관리 방법에 포함되어 있거나 액체 FCA 정기 세척 보수 방법을 사용해 별도로 실행할 수 있 습니다. "액체 경로 세척" [▶ 133] 섹션 을 참조하십시오.
세척 시스템 (고정 팁이 있는 MCA)	플러시/프라이밍	세척액	직접 명령 Prime Wash Station(MCA384)을 실행합니 다.
액체 시스템 (Liquid FCA)	세척 후 팁 또는 DiTi 콘 에 물방울이 없는지 육 안으로 확인	-	이 작업은 일일 시스템 관리 검사법에 포함되어 있습니다.
그리퍼 핑거	핑거가 곧게 펴졌는지 확인합니다. 손상 및 오정렬 여부를 검사합니다.	_	오정렬이 발생한 경우에는 "로봇 그리퍼 암(RGA) 문제 해결" [▶ 151] 섹션을 참조하 여 해결하십시오. 변형 또는 손상. "고객 지원" [▶ 188] 섹션을 참조하십시오.
FCA 그리퍼 핑거	손상 여부를 검사합니 다.	-	손상된 경우 교체하십시오. 주문 정보는 참조 설명서를 참조하십시오. 섹션을 참조하 십시오.
Frida Reader	블라인드 플러그를 제 거하고 인서트를 Frida Reader에 장착합니다.	_	"Frida Reader" [▶ 141] 섹션을 참조하십시오.

▲ 주의

Frida Reader의 측정 결과가 올바르지 않습니다!

삽입물이 장착되지 않은 경우 Frida Reader가 잘못된 측정 결과를 제공할 수 있습니다.

• 인서트를 제거했을 때는(예: 세척 등) 빨간색 블라인드 플러그를 사용하십시오.

7.4.1.2 일과 종료

다음 표에는 일과 종료 시의 일일 시스템 관리 작업이 시간 순으로 나열되어 있습 니다.



표 23: 일과 종료 시스템 관리 테이블

기기/ 구성 요소	시스템 관리 작업	세정제/ 일회용 제품/ 장치	참조/ 시스템 관리 활동
데크 트레이	유출물이 있는지 확인 하고 필요한 경우 세척 하거나 교체합니다.	탈이온수, 알코올, 약한 세제, 소독제, 염기, 표 백제, DNAzap	"세척 데크 트레이" [▶ 124] 섹 션을 참조하십시오.
세그먼트 Fluent ID 하우징	세정	탈이온수, 알코올, 약한 세제, 소독제, 염기, 표 백제, DNAzap 공지사항! 스캐너 창에 는 세그먼트 자체와 다 른 세정제가 필요합니 다. "주간 시스템 관리" [▶ 112]를 참조하십시 오.	"러너와 세그먼트 청소" [▶ 124] 섹션을 참조하십시오. 경고! 레이저 빔을 응시하지 마십시오.
리플렉터 호일 (Fluent ID, 튜브 회전 기)	청소하고 손상 여부를 검사하십시오	알코올 공지사항! 리플렉터 호 일에는 세그먼트 자체 말고 다른 청소제가 필 요합니다.	손상. "Fluent ID 리플렉터 호 일 교체" [▶ 126] 섹션을 참조 하십시오.
러너	세정	탈이온수, 알코올, 약한 세제, 소독제, 계면 활 성제, 약한 산성, 염기 성, 표백제, DNAzap	"러너와 세그먼트 청소" [▶ 124] 섹션을 참조하십시오.
고정 팁 블럭(MCA)	파킹, 세척, 점검 및 커 버	알코올	"고정 팁 블록 세척(MCA 384)" [▶ 139] 섹션을 참조하 십시오. 주의! Tecan은 모든 작업 교 대 시간이 끝날 때 팁 블록을 파킹하고 기기에서 제거할 것 을 권장합니다. 항상 팁 블록 을 세정하고 적절히 보관하십 시오. 팁 위치를 확인합니다.
고정 팁	세정	알코올, 표백제, 보풀 없는 천	"고정 팁 세척" [▶ 119] 섹션을 참조하십시오.
일회용 팁 콘	세정	알코올, 보풀이 없는 천	"일회용 팁 콘 세척" [▶ 118] 섹 션을 참조하십시오.
세척 및 폐기물 스테이 션(액체 FCA)	세정	탈이온수, 알코올, 약한 세제, 소독제	"일회용 팁 폐기물 및 세척 스 테이션 유닛 세척" [▶ 128] 섹 션을 참조하십시오.
일회용 팁 폐기물 슬라 이드 및 폐기물 커버	세정	탈이온수, 알코올, 약한 세제, 소독제	"일회용 팁 폐기물 슬라이드 세척" [▶ 129] 섹션을 참조하 십시오.



기기/ 구성 요소	시스템 관리 작업	세정제/ 일회용 제품/ 장치	참조/ 시스템 관리 활동
액체 시스템 (Liquid FCA)	세척	시스템 액체 공지사항! 액체 시스템 에 염분 함량이 높은 경 우 탈이온수로 씻어냅 니다.	액체 FCA 정기 세척 보수 방 법을 실행합니다.
MCA 세척 스테이션 (고정 팁이 있는 MCA)	세정	알코올, 표백제, Decon 90, Contrad 공지사항! Decon 또는 Contraad와 같은 표면 활성 물질이 프로세스 에 영향을 줄 수 있습니 다. 따라서 이러한 세정 제를 사용하는 경우, 절 차를 주의 깊게 확인하 십시오.	직접 명령 Prime Wash Station(MCA384)을 실행합니 다.
일회용 팁 폐기물 백	변경 	권장 백 사양: 너비 x 길이: 300mm x 600mm; 두 께: 0.5mm 소재: 폴리프로필렌, 폴 리에틸렌 또는 공중합 체(고압 멸균 가능) 공지사항! 사용한 폐기 물 백은 현지 안전 지침 을 준수해야 합니다.	"일회용 팁 폐기물 슬라이드 세척" [▶ 129] 섹션을 참조하 십시오.
시스템 액체 용기 (Liquid FCA)	깨끗한지 확인합니다.	시스템 액체	"시스템 액체 용기 및 폐기물 용기 연결" [▶ 134] 섹션을 참 조하십시오.
폐기물 용기 (고정 팁이 있는 액체 FCA 및 MCA)	비우고 세척	탈이온수, 알코올, 약한 세제, 계면 활성제, 소 독제, 염기, 표백제	현지 검사실 규칙/규정에 따 라 매일 또는 매주 세척합니 다. "시스템 액체 용기 및 폐기물 용기 연결" [▶ 134] 섹션을 참 조하십시오.
안전 패널	세정	탈이온수, 알코올, 약한 세제	"안전 패널 세척" [▶ 128] 섹션 을 참조하십시오.
튜브 로테이터	표면, 다운홀더 및 세척 스테이션 세척	2% 표백제, 70% 에탄 올 또는 100% 이소프 로판올의 용액을 적신 보풀이 없는 천	"튜브 로테이터 세정" [▶ 119] 섹션을 참조하십시오.



기기/ 구성 요소	시스템 관리 작업	세정제/ 일회용 제품/ 장치	참조/ 시스템 관리 활동
피어싱 팁	관찰하여 관절 끝부분 에 침착물이 있는지 확 인합니다. 필요한 경우 세척합니다.	70% 에탄올 또는 2% 표백제 및 보풀이 없는 천	"피어싱 팁 세정" [▶ 119] 섹션 을 참조하십시오.
피어싱 팁 세척 스테이 션 및 폐기물 튜브	잔여 샘플 물질을 제거 한 후 세척 스테이션과 튜브를 깨끗이 청소합 니다.	탈이온수, 약한 세제, 소독제 세척 스테이션 구성품 을 물로 헹구지 않은 상 태에서 표백제를 사용 하지 마십시오.	모든 구획(세척 스테이션의 전면, 가운데, 후면)을 포함하 여 세척 스테이션의 확장된 영역 세척을 실행합니다.
Frida Reader	인서트를 분리하고 블 라인드 플러그를 Frida Reader에 장착	-	"Frida Reader" [▶ 141] 섹션을 참조하십시오.

7.4.2 주간 시스템 관리

매주 마지막 근무일에 주간 시스템 관리 작업을 수행해야 합니다.

핵심 운영자가 수행할 수 있는 경우 WeeklySystemCare 방법을 실행하거나 아래 표에 나온 시간 순서대로 일일 작업 외에도 Fluent 암 구성에 해당하는 개별 작업을 수행합니다.

표 24: 주간 시스템 관리 테이블

기기/ 구성 요소	시스템 관리 작업	세정제/ 일회용 제품/ 장치	참조/ 시스템 관리 활동
세척 시스템 (고정 팁이 있는 MCA 384)	튜브와 필터에 오물이 있는지 육안으로 확인	필터	더러운 필터. "고객 지원" [▶ 188] 섹션을 참조하십시오.
피펫팅 헤드 (MCA 384)	누출 테스트 수행	_	MCA 384 누출 및 제로 분사 검사법을 실행합니다.
플레이트 어댑터 (MCA 384)	세정	알코올 압축 공기 (건조용)	"세척 플레이트 어댑터(MCA 384)" [▶ 139] 섹션을 참조하 십시오.
액체 시스템 (Liquid FCA)	세정	Fluent에서 액체를 처 리하는 방식에 따라 Decon, Contrad, 염기 성, 약한 산성, 소독제 이후 물, 알코올 및 시 스템 액체 플러시	"액체 경로 세척" [▶ 133] 섹션 을 참조하십시오.



기기/ 구성 요소	시스템 관리 작업	세정제/ 일회용 제품/ 장치	참조/ 시스템 관리 활동
Liquid FCA	밸브 인터페이스에서 주사기의 제대로 조여 졌는지, 플런저 잠금 나 사에서 주사기 플런저 가 제대로 조여졌는지 확인합니다.		"주사기가 조여졌는지 확인" [▶ 134] 섹션을 참조하십시오.
DiTi 콘	DiTi 콘의 조임 상태를 점검합니다	-	"DiTi 콘 조이기" [▶ 140] 섹션 을 참조하십시오.
Liquid FCA	누출 테스트 수행(액체 FCA)	_	Liquid FCA 누출 테스트 를 실 행합니다.
Air FCA	공기 누출 테스트(FCA) 수행	-	AirFCA 누출 테스트 및 cLLD 자가 테스트 를 수행합니다.
Air FCA MultiSense	MultiSense Air FCA cLLD 자가 테스트 수행	_	AirFCA 누출 테스트 및 cLLD 자가 테스트 를 수행합니다.
시스템 액체 용기	세정	탈이온수, 알코올, 약한 세제, 계면 활성제, 소 독제, 염기, 표백제	"시스템 액체 용기 및 폐기물 용기 세척" [▶ 134] 섹션을 참 조하십시오.
세척 스테이션 (Liquid FCA)	세정	세제 또는 소독 용액	_
RGA 그리퍼 핑거 패드	그리퍼 핑거 패드에서 입자와 잔여물 제거	알코올을 묻힌 보풀 없 는 천	세정제로 문질러서 닦습니다.
도킹 스테이션 및 그리 퍼 핑거(연결 인터페이 스)	그리퍼 핑거 연결 인터 페이스(PCBA, 자석 및 콘체)에서 입자와 잔여 물 제거	알코올을 묻힌 보풀 없 는 천	세정제로 문질러서 닦습니다.
독립 실행형 바코드 스 캐너 창	세정	약한 세제	경고! 레이저 빔을 응시하지 마십시오. 바코드 스캐너 제조업체의 설 명서를 참조하십시오. "레이저 방사선 기기" [▶ 35] 섹션을 참조하십시오.
Fluent ID와 튜브 로테 이터 스캐너 창	먼지 및 손상 여부 점검 필요한 경우 세정	약한 세제 헹구는 데 사용할 탈이 온수	경고! 레이저 빔을 응시하지 마십시오. 부드러운 천으로 세척하고 헹 구십시오.
Fluent ID와 튜브 로테 이터 리플렉터	먼지 및 손상 여부 점검 필요한 경우 세정	약한 세제 헹구는 데 사용할 탈이 온수	경고! 레이저 빔을 응시하지 마십시오. 부드러운 천으로 세척하고 헹 구십시오.



기기/ 구성 요소	시스템 관리 작업	세정제/ 일회용 제품/ 장치	참조/ 시스템 관리 활동
FCA 그리퍼	세정	알코올	-
튜브 로테이터	잠금 및 위치 지정 핀의 존재 여부 및 조임 상태 를 점검하십시오. 필요 한 경우 핀을 조이거나 교체하십시오.	_	"잠금 핀 및 위치 지정 핀 교 체" [▶ 183] 섹션을 참조하십 시오.
MCA 96	누출 테스트 수행	-	MCA 96 누출 검사법을 구현 하여 실행합니다.
MCA 96	피펫팅 성능 테스트 수 행		다음과 같은 경우 MCA 96 피 펫팅 성능 검사법을 실행하십 시오. • 콘실 수명이 90%에 도달 함 • 10네 또는 50네 팁이 사용 됨 Fluent Control은 콘 밀봉 수 명의 90%에 도달하면 이로 인한 영향에 대한 경고를 표 시합니다.

7.4.3 월간 시스템 관리

다음 표에는 월간 시스템 관리 작업이 시간 순으로 나열되어 있습니다.

표 25: 월간 시스템 관리 테이블

기기/ 구성 요소	시스템 관리 작업	세정제/ 일회용 제품/ 장치	참조/ 시스템 관리 활동
소프트웨어	컴퓨터 재시작	-	컴퓨터를 끕니다. 10초 기다 립니다. 컴퓨터를 재시작합니 다.
암 가이드	세정	면봉이나 스크루드라이 버에 보풀이 없는 천 사 용	"암 가이드 세척" [▶ 140] 섹션 을 참조하십시오.
MCA 96	MCH 96 청소	알코올을 묻힌 보풀 없 는 천, 압축 공기	"MCH 96 청소" [▶ 142] 섹션을 참조하십시오.

7.4.4 정기 시스템 관리



이러한 작업을 수행해야 하는 간격은 핵심 운영자가 결정해야 합니다.



다음 표에는 시스템 관리 작업이 시간 순으로 나열되어 있습니다.

표 26: 정기 시스템 관리 표

기기/ 구성 요소	시스템 관리 작업	세정제/ 일회용 제품/ 장치	참조/ 시스템 관리 활동
콘 슬리브 연결부	입자 제거 표면 청소	알코올, 보풀이 없는 천	
UVC 조명	지문이 있는지 확인합 니다. 필요한 경우 세척합니 다.	알코올, 보풀이 없는 천	
MCA 96 그리퍼 핑거	특히 그리퍼 핑거의 충 돌 후에 그리퍼 핑거 육 안 검사	그리퍼 핑거에 먼지와 손상이 있는지 점검합 니다. 더러운 경우 알코 올과 보풀이 없는 천을 사용하여 닦습니다.	

7.4.5 연간 시스템 관리

연간 시스템 관리는 정확도와 정밀도를 유지하고 기기 가동 중지 시간을 최소화하는 데 도움이 됩니다. Fluent의 수명을 연장하는 데도 도움이 됩니다.

현지 Tecan 서비스 조직에 문의하여 연간 시스템 관리 일정을 예약하십시오. "고 객 지원" [▶ 188] 섹션을 참조하십시오.

7.4.6 격년제 시스템 관리

다음 예방 유지보수 작업은 2년마다 수행해야 합니다.

표 27: 격년제 시스템 관리

구성 요소	작업	참조
FCA 그리퍼	FCA 그리퍼의 핑거를 교체합 니다. FluentControl 의 카운터를 재 설정합니다.	주문 정보는 참조 설명서를 참조하십시오. "참조 문서" [▶ 10]를 참조하십시오.

7.5 시스템 관리 활동

아래 설명된 시스템 관리 작업을 수행하려면 다음과 같은 절차를 따르십시오.

- 시스템 관리 모드로 전환합니다. "시스템 관리 모드" [▶ 106] 섹션을 참조하십시 오.
- 아래 설명된 지침을 따르십시오.



7.5.1 실험실 내 캐비닛에서 기기 이동하기

▲ 주의

캐비닛이 손상됩니다!

원심 분리기 설치 또는 폐기물 카트와 같은 캐비닛 선반이 제거되었을 수 있습니 다. 장치가 설치된 캐비닛 위에 놓인 악기를 이동하면 캐비닛에 손상을 유발할 수 있고 부상을 일으킬 수도 있습니다.

- 기기를 옮기기 전에 캐비닛 선반을 설치하십시오.
- 계단이나 구멍이 없는 평평한 바닥에서만 캐비닛을 옮기십시오. 계단이나 패인 지형이 나오면 Fluent 리프팅 바를 사용하여 장애물 위로 시스템을 들어 올리 거나 서비스 담당자에게 문의하십시오.

룸 내의 캐비닛 위에서 기기를 이동하려면 다음과 같은 절차를 따르십시오.

- 캐비닛이 안전한 곳에 배치되어 있고 굴러서 이동하지 않도록 고정되어 있는지 확인합니다.
- 2. 캐비닛 선반(A)이 설치되어 있는지 확인합니다.



- 3. 오픈 엔드 스패너를 사용하여 캐비닛 피트(B)의 너트를 돌립니다.
- 잠금이 해제되고 휠이 이동 위치에 올 때까지 캐비닛 피트(C)의 빨간색 나사 (D)를 돌립니다.



5. 캐비닛의 기기를 새 위치로 이동합니다.



 캐비닛이 안전한 곳에 배치되어 있고 굴러서 이동하지 않도록 고정되어 있는지 확인합니다.

7.5.1.1 기기 수평 조정

기기 수평을 맞추려면 다음과 같은 절차를 따르십시오.

 오픈 엔드 렌치를 사용하여 캐비닛 휠을 손으로 돌릴 수 있을 때까지 조정 가능 한 모든 피트를 내립니다.



2. 해당 발의 잠금 너트(A)를 느슨하게 풀어 봅니다.



아래 나열된 그리드 위치에 따라 참조 세그먼트를 배치합니다.
 기기 크기 480: 그리드 위치 왼쪽 1 및 그리드 위치 오른쪽 21.
 기기 크기 780: 그리드 위치 왼쪽 1 및 그리드 위치 오른쪽 41.





기기 크기 1080: 그리드 위치 왼쪽 1 및 그리드 위치 오른쪽 59.

- 4. 수평 수준기를 사용하여 기기가 가로 및 세로로 정렬되었는지 확인합니다.
- 필요에 따라 캐비닛 수준을 조정합니다(올리려면 시계 방향으로, 낮추려면 시 계 반대 방향으로).



- 6. 기기를 수평으로 조정한 후 캐비넷의 발에 있는 잠금 너트를 다시 조입니다.
- 캐비닛이 안전한 곳에 배치되어 있고 굴러서 이동하지 않도록 고정되어 있는지 확인합니다.

7.5.2 일회용 팁 콘 세척

일회용 팁 콘을 세척하려면 다음과 같은 절차를 따르십시오.

- 1. 알코올과 보풀이 없는 천으로 1회용 팁 콘을 깨끗이 세척합니다.
- 시스템 관리 과정에서 일회용 팁 콘과 팁의 돌출 부분을 점검합니다.
 액체 FCA의 경우: 콘의 외부에 튀어나온 튜빙 확장 부분이 손상되지 않았는지 확인하십시오.
- 3. 튜브 확장부가 깨끗하고 침전물이 없는지 확인합니다.



7.5.3 고정 팁 세척

▲ 주의

세척 중 고정 팁으로 인한 부상 위험 있음

피펫팅에서 고정 팁을 사용할 경우 부상 위험이 있습니다.

 작업대 접근 시 적절한 보호복을 착용하여 피펫팅 팁 및 에어로졸과 접촉하지 않도록 합니다.

고정 팁을 청소하려면 다음과 같은 절차를 따르십시오.

- 1. 보풀이 없는 천을 사용하여 고정 팁을 알코올로 닦습니다.
- 2. 고정 팁이 깨끗하고 침전물이 없는지 확인합니다.

7.5.4 피어싱 팁 세정

피어싱 팁을 세정하려면 **피어싱 팁 세정 유지관리** 방법을 따르십시오. 이 검사법은 작업대 설정에 따라 조정되어야 합니다.

스크립트에는 다음 단계가 포함됩니다.

- 1. 작업대(예: 실험 기구 및 하드웨어)를 준비하십시오.
- 튜브 로테이터 또는 튜브 다운홀더 캐리어에서 마개가 닫힌 비어 있는 튜브 8 개의 Z-시작 위치를 피어싱합니다.
- 70% 에탄올 또는 2% 표백제로 희석한 용액을 묻힌 보풀이 없는 천을 사용하 여 피어싱 팁의 접근 가능한 부분을 직접 세척합니다. 피어싱 팁의 날카로운 끝 부분과 접촉을 피하여 안전하게 사용하십시오.
- 4. 수동 세척 후 세척 명령을 수행합니다.

7.5.5 튜브 로테이터 세정

일반 세척 절차

- 1. 튜브 로테이터의 부품을 세척하려면 보풀이 없는 천에 3가지 세정제, 즉 2% 표백제, 70% 에탄올, 100% 이소프로판올 중 하나로 적십니다.
- 해당 세정제를 묻힌 천으로 부품을 닦아 청소한 후 소독합니다.
 보풀이 없는 천으로 닿지 않는 부분은 면봉을 사용하여 닦습니다.
- 3. 세정액을 바른 후 5분 이내에 물에 적신 천으로 세정액을 닦아냅니다.



튜브 하향 고정판 제거 및 세척

1. 다운홀더 플레이트(A)를 해제하려면 한 손으로 다운홀더 플레이트(B)를 잡고 다른 손으로 다운홀더 잠금 핀(B)을 당깁니다.



2. 튜브 로테이터에서 하강판을 제거합니다.



 위의 일반 지침에 따라 다운홀더 플레이트를 세척하거나, 튜브 다운홀더를 최 대 2시간 동안 2% 표백제로 희석된 액체를 포함한 수조에 담가둘 수 있습니 다.

튜브 회전기 표면 세척

1. 위의 일반 지침에 따라 튜브 로테이터의 접근 가능한 표면을 세척하십시오.





2. 드럼 위치를 수동으로 변경하려면 한 손으로 드럼을 잡고 솔레노이드 분리 버 튼을 누릅니다.

- 3. 드럼을 수동으로 회전시키고 솔레노이드 분리 버튼을 해제합니다.
- 4. 드럼이 솔레노이드에 의해 잠길 때까지 돌립니다.
- 5. 위의 일반 지침에 따라 접근이 불가능한 표면을 세척하십시오.

다운 홀더 플레이트 장착

- 1. 튜브 로테이터 드럼의 위쪽면에 다운홀더 플레이트를 올려놓습니다.
- 한 손으로 기기 하단을 향해 다운홀더 플레이트를 누르고 검은색 슬라이더를 뒤로 밀어 다운홀더 플레이트를 제 위치에 고정합니다.





7.5.6 튜브 로테이터 세척 스테이션 세정

일반 세척 절차

- ✔ 세척 스테이션은 작업대에서 세척하거나 세척을 위해 해제할 수 있습니다.
- ✔ 천이 아닌 병을 닦는 브러시를 사용하면 더 편리하게 닦을 수 있습니다.
- 1. 세척 스테이션의 부품을 세척하려면 보풀이 없는 천에 3가지 세정제, 즉 2% 표백제, 70% 에탄올, 100% 이소프로판을 중 하나로 적십니다.
- 2. 적신 천으로 부품을 닦아 청소한 후 소독합니다.
- 3. 세정액을 바른 후 5분 이내에 물에 적신 천으로 세정액을 닦아냅니다.

튜브 로테이터 세척 스테이션 장착 해제

 세척 스테이션 해제 레버를 바코드 스캐너 하우징 쪽으로 누르고 다른 손으로 세척 스테이션을 들어 올립니다.







2. 폐기물 튜브를 분리하고 커넥터를 폐기물 튜브 홀더에 놓습니다.

튜브 로테이터 세척 스테이션 장착 1. 폐기물 튜브 커넥터를 연결합니다.







 세척 스테이션을 안내 축(A)에 장착하고 베이스 플레이트 쪽으로 누릅니다. 해제 레버가 제자리에 다시 맞물려 세척 스테이션을 제자리에 고정하는지 확인 합니다.

7.5.7 러너와 세그먼트 청소

러너와 세그먼트를 청소하려면 다음 단계에 따라 진행하십시오.

공지사항

액체 감지 기능 오작동(cLLD)!

러너 및 데크 세그먼트 사이에서 접촉부 오염으로 인해 액체 감지(cLLD)가 오작동 할 수 있습니다.

러너와 세그먼트가 깨끗하고 건조한지 항상 확인하십시오.

- 기기 데크에서 러너를 제거합니다.
 세그먼트 및 네스트는 제자리에서 세척합니다.
- 러너, 세그먼트 및 니들의 표면을 세정제로 닦습니다.
 러너. 세그먼트 및 네스트를 탈이온수로 헹굽니다.
- 3. 러너를 기기 데크에 다시 놓습니다.

7.5.8 세척 데크 트레이

데크 트레이를 세척하려면 다음과 같은 절차를 따르십시오.

- ✓ 데크 트레이 위의 세그먼트는 제거됩니다. "세그먼트 제거" [▶ 78]를 참조하십시 오.
- ✓ Fluent ID 또는 MCA Active Carrier와 같은 데크 세그먼트를 제거할 수 없는 경 우 데크 트레이를 열린 데크 위치로 밉니다.
- 1. 데크 트레이를 기기에서 제거합니다.



- 해당 액체에 대한 검사실 처리 프로토콜에 따라 액체를 제거하여 트레이를 비 웁니다.
- 3. 데크 트레이가 손상되었거나 분실된 경우 트레이를 교체해야 합니다.
- 4. 세정제로 데크 트레이의 표면을 닦습니다.
- 데크 트레이를 기기에 다시 장착합니다.
 아래 그림과 같이 데크 트레이의 방향을 조정합니다.
 인접한 데크 트레이는 서로 맞물려 있어야 합니다.



그림 44: 데크 트레이 배치 잘못됨



그림 45: 교차식 데크 트레이





그림 46:데크 트레이를 올바르게 배치

7.5.9 Fluent ID 리플렉터 호일 교체

- ✔ 자체 접착식 리플렉터 호일
- 1. 리플렉터 호일을 가열합니다. 히트 건을 사용하십시오.
- 2. 리플렉터 호일을 제거합니다.



3. 알코올로 모든 잔여물을 제거합니다.





4. 리플렉터의 위쪽 끝에 새 접착식 리플렉터 호일을 부착합니다.

7.5.10 DiTi 폐기물 슬라이드에 Fluent ID 리플렉터 호일 적용

- ✔ 자체 접착식 리플렉터 호일
- 1. 아래 그림에 따라 DiTi 폐기물 슬라이드에 테이프를 부착합니다.



 아래 그림에 따라 새 접착식 리플렉터 호일을 DiTi 폐기물 슬라이드에 부착합 니다.



레이저 빔은 반드시 리플렉터 호일 중앙에 있어야 합니다.



3. DiTi 폐기물 슬라이드에서 테이프를 제거합니다.

7.5.11 안전 패널 세척

안전 패널을 청소하려면 다음과 같은 절차를 따르십시오.

1. 세정제로 안전 패널의 내부 및 외부 표면을 닦습니다.

7.5.12 일회용 팁 폐기물 및 세척 스테이션 유닛 세척

일회용 팁 폐기물과 세척 스테이션 장치를 청소하려면 다음과 같은 절차를 따르십 시오.

- 1. 신속 해제 패스너 버튼(B)을 누릅니다.
- 2. 세척 스테이션을 뒤로 밉니다.



그림 47: 백 하우징 탈거용 고정 장치

3. 세척 스테이션을 일회용 팁 폐기물과 세척 스테이션 유닛에서 제거합니다.





그림 48: 세척 스테이션을 제거합니다

- 4. 세척액으로 세척 스테이션 표면을 닦아내고 쏟아진 시약을 제거합니다.
- 5. 신속 해제 패스너 버튼(B)을 누릅니다.
- 6. 세척 스테이션(A)을 제자리에 놓습니다.
- 7. 세척 스테이션을 앞으로 밉니다.



그림 49: 세척 스테이션을 재장착합니다

7.5.13 일회용 팁 폐기물 슬라이드 세척

일회용 팁 폐기물 슬라이드를 세척하려면 다음과 같은 절차를 따르십시오.

- ✔ 전면 안전 패널이 열려 있습니다.
- 1. 커버(A)를 일회용 팁 폐기물 슬라이드에서 제거합니다.





그림 50: 폐기용 팁 슬라이드의 커버를 제거합니다. 2. 홀더에서 일회용 팁 폐기물 슬라이드(B)를 제거합니다.



- *그림 51:* 일회용 팁 폐기물 슬라이드 제거
- 일회용 팁 폐기물 슬라이드 하단 개구부(C) 아래에 조직을 고정합니다.
 오염된 물질이 떨어지지 않도록 하십시오.



그림 52: 일회용 팁 폐기물 슬라이드 처리

- 4. 일회용 팁 폐기물 슬라이드 및 커버를 세정제가 들어 있는 통에 넣습니다.
- 5. 30분에서 4시간 동안 담가 둡니다.
- 일회용 팁 폐기물 슬라이드 및 커버를 통에서 꺼내 깨끗한 마른 수건 위에 놓습 니다.
- 7. 그대로 두어 건조합니다.



8. 일회용 팁 폐기물 슬라이드(B)를 홀더에 다시 설치합니다.



그림 53: 일회용 팁 폐기물 슬라이드 다시 삽입 후 다시 설치 9. 위치 지정 핀이 슬롯(D)에 올바르게 삽입되었는지 확인합니다. 10 폐기물 슬라이드 위에 커버(A)를 놓습니다.



그림 54: 위치 지정 핀 및 덮개

7.5.14 일회용 팁 폐기물 백 교체

일회용 팁 폐기물 백을 변경하려면 다음과 같은 절차를 따르십시오. 1. 패스너(A)를 들어 올리고 백 하우징을 앞쪽으로 밉니다.



그림 55: 백 하우징 탈거용 고정 장치



- 2. 일회용 백 하우징(A)을 제거합니다.
- 3. 일회용 팁 폐기물 백(B)을 제거합니다.
- 4. 연구소 지침에 따라 사용 후 폐기되는 팁 폐기물 백을 적절하게 처리하십시오.
- 5. 새 일회용 폐기물 백(B)을 빈 백 하우징(B)에 삽입합니다.



그림 56: 백 하우징 및 일회용 팁 폐기물 백



그림 57: 작업대에 폐기물 처리구를 적절하게 설치하는 방법



▲ 주의

폐기물 슈트를 잘못 배치할 경우 암이 충돌하거나 팁이 잘못 이탈할 수 있습니다. 폐기물 슈트가 아래 그림과 같이 올바른 위치에 있는지 확인하십시오. (폐기물 슈 트 배치에서 그림 삽입, 올바른 항목 및 잘못된 항목 표시)

6. 백 하우징을 제 위치로 밀어 넣고 패스너(A)로 닫습니다.





그림 58: 체결 부품

7.5.15 액체 경로 세척

- 고정 팁 안쪽에서 단백질 잔여물을 제거하려면 먼저 약한 산성 세정제를 사용 한 후 염기 세정제를 사용합니다.
- 고정 팁 내부에서 핵산 잔여물을 제거하려면 염기 세정제를 사용하십시오.
- Decon/Contraad와 같은 세정제는 절차를 영향을 줄 수 있습니다. 따라서 이러 한 세정제를 사용하는 경우, 절차를 주의 깊게 확인하십시오.
- 이소프로판올은 매우 효과적인 소독제입니다. 빠르게 증발시켜 표면을 사용 가 능하도록 준비합니다.
- 허용되는 세정제만 사용하십시오. 표백제 용액을 사용해 전체 액체 시스템을 세척하지 마십시오.

액체 경로를 세척하려면 다음과 같은 절차를 따르십시오.

- 1. 액체 용기에서 시스템 액체 튜브를 분리합니다.
- 2. 유지보수 튜브(30043739)를 시스템 튜브에 연결합니다.
- 3. 유지보수 튜브의 열린 끝을 세정제가 들어 있는 병에 놓습니다.
- 4. 세정제(RapidWash 20ml, 희석제 10ml)로 세척합니다.
- 5. 20분 동안 담가 둡니다.
- 6. 탈이온수를 가득 채운 병에 튜빙을 놓습니다.
- 7. 탈이온수로 두 번 헹굽니다(RapidWash 20ml 및 희석제 10ml).

▲ 경고

가연성 액체!

가연성 액체 또는 시스템 액체로 인한 화재 위험이 있습니다.

- 가연성 증기가 형성되거나 축적되지 않도록 하십시오.
- 데크 트레이 없이 시스템을 작동하지 마십시오.
- 8. 튜브를 알코올이 들어있는 병에 넣습니다.
- 9. 알코올로 씻습니다(RapidWash 20ml 및 희석제 10ml).
- 10 시스템 튜브에서 유지보수 튜브를 제거하고 시스템 튜브를 시스템 액체 용기에. 연결합니다.
- 11 탈이온수로 두 번 씻습니다(RapidWash 20ml 및 희석제 용량 5배).



12 튜브에 기포가 있는지 확인합니다.

13 기포가 보이면 다시 씻습니다.

7.5.16 시스템 액체 용기 및 폐기물 용기 연결

시스템 액체 용기와 폐기물 용기를 준비하려면 다음과 같은 절차를 따르십시오.



Tecan에서 시스템을 제어하는 원래 컨테이너를 사용하는 경우에만 고장 없이 작 동할 수 있습니다. 처음 사용하기 전에 시스템 액체 용기를 수동으로 완전히 헹궈서 병 내부에서 고형 이물질을 제거해야 합니다. "시스템 액체 용기 및 폐기물 용기 세척" [▶ 134] 섹션을 참조하십시오.

- ✔ 용량이 20리터 이상인 Tecan 용기
- 1. 액체 감지 시스템(A, B)이 올바르게 연결되어 있는지 확인합니다.
- 2. 튜브(C, D)가 올바르게 연결되어 있는지 확인합니다.



7.5.17 시스템 액체 용기 및 폐기물 용기 세척

액체 용기와 폐기물 용기를 청소하려면 다음과 같은 절차를 따르십시오.

- 1. 세척액 용기를 수동으로 비웁니다.
- 2. 세정제가 든 통에 액체 용기를 넣고 세척한 후 헹굽니다.
- 3. 액체 용기를 알코올로 소독합니다.
- 4. 시스템 액체 및 폐기물 용기를 연결합니다("시스템 액체 용기 및 폐기물 용기 연 결" [▶ 134] 섹션 참조).

7.5.18 주사기가 조여졌는지 확인

다음과 같이 주사기가 올바로 조여졌는지 점검합니다.





그림 59: 조임 점검

Α

플런저 잠금 나사 B 주사기 나사

- 공기를 흡입하는 유지보수 스크립트를 사용하여 플런저를 주사기 중앙으로 이 동합니다.
 참고: 유지보수 스크립트는 실험실의 FluentControl 관리자가 제공해야 합니 다.
- 2. 주사기 나사(B)를 조입니다. 즉, 오른쪽으로 돌립니다.
- 3. 플런저 잠금 나사(A)를 시계 방향으로 돌려 조입니다.

7.5.19 개스킷 점검(MCA 384)

MCA 개스킷을 점검하려면 다음 단계를 따르십시오.

1. 개스킷(A)이 손상되었는지 검사합니다.



개스킷이 손상되지 않았는지 확인합니다. 손상된 개스킷을 교체해야 합니다. "개스킷 교체(MCA 384)" [▶ 136] 섹션을 참조하십시오.



7.5.20 개스킷 교체(MCA 384)



그림 60: 부품 및 도구

- A 개스킷
- C 뭉툭한 튜브 삽입 도구
- E 블런트 튜브 제거 공구
- B 블런트 튜브
- D 개스킷 삽입 공구

개스킷(MCA)을 교체하려면 다음과 같은 절차를 따르십시오.

- 1. 어댑터 플레이트를 제거합니다.
- 2. 피펫팅 헤드를 앞으로 움직이고 가능한 한 높게 올립니다.
- 3. 블런트 튜브 제거 공구를 채널에 약 2mm 정도 밀어 넣습니다.





4. 도구를 시계 방향으로 돌려 뭉툭한 튜브를 고정합니다.

5. 뭉툭한 튜브를 채널 밖으로 당깁니다.







6. 롱 노즈 플라이어를 사용하여 채널에서 개스킷을 제거합니다.

- 7. 새 개스킷을 미네랄 오일 필름으로 윤활합니다.
- 8. 개스킷 삽입 공구를 사용하여 윤활 처리된 개스킷을 채널에 삽입합니다.



9. 새 뭉툭한 튜브의 바깥 표면을 미네랄 오일로 윤활합니다.



- 10 뭉툭한 튜브 삽입 도구를 사용하여 윤활된 뭉툭한 튜브를 채널에 천천히 삽입. 합니다.
 - 뭉툭한 튜브는 채널 내부에 완전히 삽입되어야 합니다.



11 새 개스킷을 검사하여 다른 개스킷과의 균일성을 확인합니다.

· · · · · ·

12 MCA 누출 테스트 검사법을 실행합니다.

7.5.21 고정 팁 블록 세척(MCA 384)

- 1. 보풀이 없는 천을 사용하여 알코올로 고정 팁 블록(MCA)을 세척합니다.
- 2. 보풀이 없는 티슈로 팁 블록을 닦거나 무급유 압축 공기로 팁 블록을 말립니다.
- 3. 고정 팁 블록(MCA)은 깨끗하고 침전물이 없는 상태로 유지하여야 합니다.
- 4. 팁 블록을 팁 블록 상자에 보관합니다.
- 팁 블록을 취급할 때 오염을 방지하려면 다음과 같이 수행해야 합니다.
- 팁 블록은 먼지가 없는 곳에 보관해야 합니다.
- 손가락으로 팁을 절대로 만지면 안 됩니다. 팁 블록은 항상 처리 중 PEEK 블록 에 의해 고정되어야 합니다.
- 테이블 위에 팁 블록의 팁을 놓지 말아야 합니다.

7.5.22 세척 플레이트 어댑터(MCA 384)

플레이트 어댑터를 청소하려면 다음과 같은 절차를 따르십시오.

- 1. 채널에 먼지가 있는지 확인합니다.
- 2. 알코올을 사용하여 기름을 제거합니다.
- 3. 압축 공기로 건조합니다.



7.5.23 암 가이드 세척

암 가이드를 세정하려면 다음과 같은 절차를 따르십시오.

- 암 가이드의 암 가이드 롤러(A)를 면봉이나 스크루드라이버에 보풀이 없는 천 을 덮어 닦습니다.
- 2. 보풀이 없는 천으로 암 레일(B)을 청소합니다.
- 보풀이 없는 천으로 MCA 암 가이드에 있는 가이드 레일의 상단 표면을 청소합 니다.



7.5.24 DiTi 콘 조이기

FCA DiTi 콘을 조정하려면 다음과 같은 절차를 따르십시오.

- 1. 팁 어댑터(D)와 팁 이젝터 튜브(C)를 잡습니다.
- 2. DiTi 콘 렌치(B)를 사용하여 DiTi 콘(A)을 조입니다.



3. FCA 정기 유지 관리 검사법을 실행합니다.



7.5.25 Frida Reader

삽입



그림 61: Frida Reader 삽입물 설치를 위해 인서트를 Frida Reader에 끼우고 표시를 정렬하십시오.

블라인드 플러그



그림 62: Frida Reader 블라인드 플러그

블라인드 플러그는 인서트가 제거되었을 때 Frida Reader를 보호합니다. 설치하려 면 블라인드 플러그를 Frida Reader에 장착하십시오.



7.5.26 MCH 96 청소

배출 플레이트(A)에 눈에 보이는 손상이 있는지 검사합니다. 특정 공정 용액 (DMSO 또는 아세토니트릴 등) 또는 세정제(표백제 등)를 흘릴 경우 플레이트가 손 상될 수 있습니다. 이 경우 현지 서비스 조직에 문의하여 교체하십시오.



- 1. Move Tool(이동 도구)을 사용하여 배출 축을 아래쪽으로 이동합니다. 2..3.
- 무급유식 압축 공기를 사용하여 배출 플레이트(C), 콘형 플레이트 및 콘(B)의 상단 표면을 세척합니다.
- 보푸라기가 없는 천과 알코올을 사용하여 배출판(D)의 바닥 표면을 세척합니다.





아래는 청소 전의 분리판을 보여주는 왼쪽 그림과 청소 후의 분리판을 보여주는 오 른쪽 그림입니다.





8 문제 해결

Fluent에 문제가 발생한 후 작동을 재개하는 방법에 대한 도움말은 이 장을 참조하 십시오. 자세한 내용이 필요하시거나 이 설명서에서 다루지 않은 문제나 세부 사항 이 충분하지 않은 경우 "고객 지원" [188] 섹션을 참조하십시오.

8.1 이 장의 안전 지침

▲ 주의

충돌 후 손상된 팁으로 인한 교차 오염!

팁이 구부러지거나 팁 코팅이 손상되면 피펫팅이 부정확해지고 액체 감지 오류가 발생합니다.

• 충돌 후 고정 팁을 점검하십시오. "고정 팁 확인" [▶ 164] 섹션을 참조하십시오.

8.2 문제 해결 표



문제 해결 표에는 가능한 문제, 원인 및 해결 방법이 나와 있습니다. 자세한 내용이 필요하시거나 이 설명서에서 다루지 않은 문제나 세부 사항이 충분하지 않은 경우 "고객 지원" [》 188] 섹션을 참조하십시오.

8.2.1 기기 문제 해결

표 28: 기기 문제 해결 표

문제/오류	가능한 원인	시정 조치
시스템 액체 누 출	튜브 및/또는 튜브 연결 부가 단단히 연결되어 있지 않습니다. 주사기에서 누출이 발 생하고 있습니다.	"고객 지원" [▶ 188] 섹션을 참조하십시오.
통신 오류	전원이 켜져 있지 않습 니다. 전원 또는 통신이 중단 되었습니다. 통신 없음.	기기의 전원을 끕니다. 기기 상태 램프 및 전원 공급 램프가 꺼질 때까지 기다립니다. PC 전원을 끕니다. 케이블 및 플러그를 점검합니다. 기기 및 PC의 전원을 켭니다.
	X, Y 또는 Z-드라이브 가 차단되었습니다	장애물이 있는지 확인합니다. 공지사항! 암이 자유롭게 움직일 수 있는 지 확인합니다.
초기화 오류	무기를 초기화할 수 없 습니다.	장애물이 있는지 확인합니다. 공지사항! 암이 자유롭게 움직일 수 있는 지 확인합니다.
	하드웨어 문제입니다.	"고객 지원" [▶ 188] 섹션을 참조하십시오.
•TECAN•

문제/오류	가능한 원인	시정 조치
전면 안전 패널 도어 센서 및 도 어 잠금 장치가 손상됨	도어 잠금 장치의 기계 적 고장	기기의 전원을 끕니다. "고객 지원" [▶ 188] 섹션을 참조하십시오.
안전 패널이 없 거나 손상됨	안전을 보장할 수 없습 니다.	기기의 전원을 끕니다. "고객 지원" [▶ 188] 섹션을 참조하십시오.
액체 감지(cLLD) 고장	접촉 표면이 더럽습니 다. 실험 기구 및 세그먼트 간의 접촉이 부적절합 니다.	데크를 준비합니다. 섹션을 참조하십시오. 접촉면을 세척합니다. "러너와 세그먼트 청 소" [▶ 124] 섹션을 참조하십시오.
	시스템 액체는 cLLD 호 환성을 위해 전도도가 10µS/cm 이상입니다.	핵심 운영자에게 문의하십시오.

8.2.2 유연한 채널 암(FCA) 문제 해결

표 29: 유연한 채널 암 문제 해결 표

문제/오류	가능한 원인	시정 조치
느슨한 DiTi 콘 주의! 부정확한 피 펫팅 용량!	충분히 조이지 않은 DiTi 콘.	DiTi 콘을 조입니다.
일회용 팁을 가 져올 수 없음	충분히 조이지 않은 DiTi 콘.	DiTi 콘을 조입니다.
일회용 팁을 폐 기할 수 없음	충분히 조이지 않은 DiTi 콘.	DiTi 콘을 조입니다.
	DiTi를 재사용함	팁이 새것인지 확인하십시오. DiTi는 재사 용하지 않는 것을 권장합니다. 폐기물 슈트에서 일회용 팁이 배출되지 않 음/폐기물 슈트가 잘못 배치됨
일회용 팁이 폐 기물 슈트에서 배출되지 않음	쓰레기 구멍의 위치가 잘못되어 있습니다.	폐기물 슈트의 설치 위치를 정확하게 확인 하여야 합니다. "일회용 팁 폐기물 슬라이 드 세척" [▶ 129] 섹션을 참조하십시오.
팁이 단일 캐리 어에서 실험 기 구에 맞게 정렬 되지 않음	캐리어 위치가 잘못되 었습니다. 세그먼트가 제 위치에 고정되지 않았습니다. 실험 기구가 올바르게 배치되지 않았습니다.	올바른 캐리어 위치를 확인하십시오. "표 준 러너 로드" [▶ 80] 섹션을 참조하십시오. 세그먼트를 제자리에서 잠급니다. "세그먼 트 확인" [▶ 181] 섹션을 참조하십시오.



문제/오류	가능한 원인	시정 조치
팁이 여러 캐리 어에서 실험 기 구에 맞게 정렬 되지 않음	충돌로 인해 발생한 암 정렬 결함.	"고객 지원" [▶ 188] 섹션을 참조하십시오.
팁이 실험 기구 의 바닥과 충돌 했습니다.	잘못된 실험 기구입니 다. 실험 기구가 올바르게 배치되지 않았습니다.	데크의 실험 기구가 검사법 데크 레이아웃 과 일치하는지 확인합니다.
DiTi 드립	오염된 DiTi 콘을 사용 하면 누출이 발생합니 다	DiTi 콘을 청소합니다
	DiTi를 재사용함	팁이 새것인지 확인하십시오. DiTi는 재사 용하지 않는 것을 권장합니다.
오류 메시지: 압력 범위 이탈 (에어 FCA)	잘못된 DiTi 크기로 흡 입 후 젖은 상태의 인라 인 필터.	데크의 DiTi 크기가 검사법에 정의된 크기 와 일치하는지 확인합니다. DiTi 콘이 올바 르게 조여졌는지 확인 인라인 필터를 확인합니다. "인라인 필터 (Air FCA) 확인" [▶ 157] 섹션을 참조하십시 오.
오류 메시지: DiTi를 가져오지 못함 DiTi 드롭되지 않 음	자기장이 DiTi가 있는 센서에 영향을 미침.	_
액체 감지: 팁이 액체를 감 지하지 못함	느슨한 DiTi 콘	DiTi 콘을 조입니다("DiTi 콘 조이기" [▶ 140] 참조).
마모된 O-링	MultiSense 팁 어댑터	MultiSense 팁 어댑터의 O-링과 X-링을 교체합니다.

8.2.3 Mix & Pierce

표 30: 문제 해결

증상	가능한 원인	시정 조치
소프트웨어 명령으로는 피어싱 팁을 수축할 수 없음	막힌 피어싱 팁	"막힌 피어싱 팁 수축" [▶ 175] 섹션을 참조하십시오.
피어싱의 끝 부분이 손 상됨	구부러진 피어싱 팁 팁이 손상됨	피어싱 팁을 교체합니다. "피 어싱 팁 제거" [▶ 168] 및 "피어 싱 팁 설치" [▶ 171] 섹션을 참 조하십시오.



증상	가능한 원인	시정 조치
피어싱 오류	너무 건조한 피어싱 팁	물로 윤활하기(세척장소)
	팁이 손상됨	피어싱 팁을 교체합니다. "피
	구부러진 피어싱 팁	이성 탑 세거 [108] 및 피어 싱 팁 설치"[171] 섹션을 참 조하십시오.
	잘못된 피어싱 매개 변수	핵심 운영자에게 문의하십시 오.
	잘못된 이동 유형이 사용됨	핵심 운영자에게 문의하십시 오.
	잘못된 튜브를 사용함	지원되는 튜브를 사용하십시 오. "튜브 로테이터 러너" [▶ 59] 섹션을 참조하십시오.
	암의 수명이 다함	"고객 지원" [▶ 188] 섹션을 참 조하십시오.
액체 처리 문제	막힌 피어싱 팁	피어싱 팁을 씻습니다. 일반적으로 세척 절차를 확인 하십시오.
	팁이 손상됨	피어싱 팁을 교체합니다. "피 어싱 팁 제거" [▶ 168] 및 "피어 싱 팁 설치" [▶ 171] 섹션을 참 조하십시오.
	주사기가 제대로 장착되지 않 음	주사기의 조임 상태를 점검합 니다. "주사기가 조여졌는지 확인" [▶ 134] 섹션을 참조하 십시오.
	액체 시스템 내부에 기포 발 생	세척합니다. "고객 지원" [▶ 188] 섹션을 참 조하십시오.
용혈 문제	표본 희석	과도하게 높은 볼륨 또는 분 할 볼륨
		식염수 0.9%(분할 용량)
		피펫팅 속도 감소
	팁이 손상됨	피어싱 팁을 교체합니다. "피 어싱 팁 제거" [▶ 168] 및 "피어 싱 팁 설치" [▶ 171] 섹션을 참 조하십시오.
	혼합 매개변수	튜브 로테이터에서 사용된 회 전/진동 매개 변수로 인해 용 혈을 유발하지 않는지 확인



증상	가능한 원인	시정 조치
주사기 내 검체	모두	시스템을 청소합니다. "액체 경로 세척" [▶ 133] 섹션을 참 조하십시오.
	공극이 잘못 설정되어 있습니 다.	세척 절차를 확인합니다.
	주사기는 제대로 장착되지 않 았습니다.	주사기의 조임 상태를 점검합 니다. "주사기가 조여졌는지 확인" [▶ 134] 섹션을 참조하 십시오.
		FCA 누출 검사법 을 실행합니 다.
		더 큰 선행 공극.
		흡입 속도가 느립니다.
	액체 취급에서 잘못된 샘플 준비가 이루어졌습니다. 검체 소스 튜브에는 응고, 세포 파 편 등과 같은 고형 입자가 포 함되어 있습니다.	검체 액체의 피펫팅이 가능하 도록 검체 준비가 올바른지 확인합니다.
		검체 소스 튜브에 응고, 세포 조직파편 등과 같은 고형 입 자가 포함되어 있지 않은지 확인하십시오
	액체 취급에서 잘못된 샘플 준비가 이루어졌습니다. 튜브 가 제대로 채워지지 않고 피 어싱 중 선행 공극을 줄여주 는 부분 진공이 아직 남아 있	샘플 원통은 적절한 용량으로 충전해야 합니다.
		검체 소스 튜브에 진공이 포 함되지 않았는지 확인합니다.
	습니다.	가능한 잔여 진공을 보상하기 위해 선행 공극을 늘립니다.
세척 절차 중 피어싱 팁 이 휘어집니다	피어싱 팁이 세척 스테이션 클리너 구멍 중앙에 있지 않 음	세척 스테이션의 사본을 생성 하고 피펫팅 위치를 학습합니 다.

•TECAN•

증상	가능한 원인	시정 조치
잘못된 액체 레벨 감지: 특정 채널에서만	구부러진 피어싱 팁: 피어싱 팁이 구부러져 피어싱 중 튜 브 벽에 닿습니다	피어싱 팁을 교체합니다. "피 어싱 팁 제거" [▶ 168] 및 "피어 싱 팁 설치" [▶ 171] 섹션을 참 조하십시오.
	피어싱 위치가 잘못되었으므 로 피어싱 팁이 피어싱 중에 튜브 벽에 닿습니다.	Tecan 제조업체 실험 기구를 사용하십시오. "튜브 로테이 터 러너" [▶ 59] 섹션을 참조하 십시오.
		실험 기구 피펫팅 위치 학습/ 조정
	피어싱 팁의 방향이 잘못되었 습니다.	구멍이 기기 전면을 향하도록 피어싱 팁을 장착합니다. "피 어싱 팁 설치" [▶ 171] 섹션을 참조하십시오.
잘못된 액체 레벨 감지: 예상 액체 레벨의 일정 한 편차 및 감지된 액체 레벨	피어싱 힘을 가한 상태에서 암, 튜브 로테이터 및/또는 기 기의 제조 공차로 인해 액체 레벨 감지 중에 눈에 띄는 Z 오프셋이 발생할 수 있습니 다.	튜브 실험 기구 정의에서 사 용자 지정 속성 "PercingDetectionHeightCo mpensation"을 학습/조정합 니다

8.2.4 다중 채널 암(MCA) 문제 해결

표 31: 다중 채널 암 문제 해결 표

문제/오류	가능한 원인	시정 조치
팁이 캐리어와 정렬되	기계적 고장	"고객 지원" [▶ 188] 섹션을 참조하십시오.
시 않금	암 충돌	
마이크로플레이트와 피 펫팅 헤드가 완전히 평 행하지 않음	충돌	피펫팅 헤드 및 데크의 병렬성을 점검하려 면 핵심 운영자에게 문의하십시오.
피펫팅 중에 피펫팅 헤 드 중지로 오류 발생	흡입 및 분사 가속이 속도에 비해 너 무 빠릅니다.	가속은 흡입 및 분사 속도를 기준으로 합 리적인 수준이어야 합니다.
	흡인 및 분사 감속이 속도에 비해 너 무 빠릅니다.	감속은 흡입 및 토출 속도와 합리적인 관 계에 있어야 합니다.
		문제를 해결할 수 없습니다. "고객 지원" [▶ 188] 섹션을 참조하십시오.



문제/오류	가능한 원인	시정 조치
여러 또는 모든 펌핑 채 널이 누출 중	잘못된 일회용 팁, 고정 팁 블록, 고 정 팁 어댑터, 팁 콘 밀봉 또는 개스 킷.	항상 Tecan이 공급하는 일회용 팁, 고정 팁 블록, 고정 팁 어댑터, 팁 콘 씰 또는 개 스킷을 사용하십시오.
		핵심 운영자와 연락하여 시스템을 초기화 하고 누출이 있는지 확인하십시오.
	팁 콘 씰 또는 개스킷이 오래되었거 나 결함이 있습니다.	"고객 지원" [▶ 188] 섹션을 참조하십시오. 핵심 운영자와 연락하여 시스템을 초기화 하고 누출이 있는지 확인하십시오.
	피펫팅 헤드에 결함이 있습니다.	"고객 지원" [▶ 188] 섹션을 참조하십시오.
한 개의 채널에서 유출 발생	팁 콘 씰, 개스킷 또는 피펫 헤드의 기타 씰에 결함이 있습니다.	핵심 운영자와 연락하여 시스템을 초기화 하고 누출이 있는지 확인하십시오. "고객 지원" [▶ 188] 섹션을 참조하십시오.
1개의 일회용 팁이 올 바르게 픽업되지 않음	개별 일회용 팁에 결함이 있습니다. 이 일회용 팁 위치의 팁 콘 밀봉에 결함이 있습니다.	일회용 팁을 교체합니다. 문제를 해결할 수 없습니다. "고객 지원" [▶ 188] 섹션을 참조하십시오.
1개의 일회용 팁이 드 롭되지 않음	개별 일회용 팁에 결함이 있습니다. 이 일회용 팁 위치의 팁 콘 밀봉에 결함이 있습니다.	문제를 해결할 수 없습니다. "고객 지원" [▶ 188] 섹션을 참조하십시오.
	습도 조절이 잘못됨	습도가 작동 습도 제한 범위 내에 있는지 확인합니다. "환경 조건" [▶ 41] 섹션을 참 조하십시오.
여러 개 또는 모든 일회 용 팁이 드롭되지 않음	잘못된 일회용 팁이 사용되었습니 다.	항상 Tecan에서 제공한 일회용 팁을 사용 하십시오.
		문제를 해결할 수 없습니다. "고객 지원" [▶ 188] 섹션을 참조하십시오.
일회용 팁이 들어 올릴 때 일회용 팁이 들어간	캐리어가 올바르게 조정되지 않았습 니다.	모든 캐리어(기계식)을 정확하게 조정합니 다.
일회용 팁 상자를 들어 올림	X- 및/또는 Y-오프셋이 잘못 지정 되어스니다	일회용 팁 캐리어를 교체합니다.
		문제를 해결할 수 없습니다. "고객 지원" [▶ 188] 섹션을 참조하십시오.
	일회용 팁 박스가 규격에 맞지 않습 니다.	항상 생체분자 스크리닝 표준을 준수하는 일회용 팁 상자를 사용하여 안전하게 작업 하십시오.
		문제를 해결할 수 없습니다. "고객 지원" [▶ 188] 섹션을 참조하십시오.
	일회용 팁 캐리어에 결함이 있습니 다(일회용 팁 상자 리테이너 오작 동).	항상 생체분자 스크리닝 표준을 준수하는 일회용 팁 상자를 사용하여 안전하게 작업 하십시오.
		문제를 해결할 수 없습니다. "고객 지원" [▶ 188] 섹션을 참조하십시오.



문제/오류	가능한 원인	시정 조치
부정확한 피펫팅 결과	일회용 팁이 적절하게 수거되지 않 고 있습니다.	핵심 운영자에게 문의하여 애플리케이션 스크립트 및 캐리어를 확인하십시오.
	액체 취급 매개 변수가 잘못되었습 니다.	환경 매개변수와 조제 높이를 확인하려면 핵심 운영자에게 문의하십시오.
	캐리어가 제대로 조정되지 않았습니 다.	팁 코팅을 확인하십시오. 필요한 경우 팁 블록을 교체하십시오.
	피펫팅 헤드에 결함이 있습니다.	
	표준 팁의 코팅이 손상되었습니다.	핵심 운영자에게 문의하여 애플리케이션 스크립트 및 캐리어를 확인하십시오.
		팁 코팅을 확인하십시오. 필요한 경우 팁 블록을 교체하십시오.
유출	표준 팁의 코팅이 손상되었습니다.	코팅을 확인하십시오. 고정 팁 블록 또는 고정 팁 어댑터를 교체하십시오.
	부적합한 애플리케이션 스크립트.	핵심 운영자와 연락하여 세척 또는 액체 취급 설정을 조정하고, 필요한 경우 다른 세척 완충액을 선택하여 사용할 수 있습니 다.
	세척 채널이 막혔습니다.	세척 스테이션을 세척합니다. 문제를 해결할 수 없습니다. "고객 지원" [▶ 188] 섹션을 참조하십시오.

8.2.5 로봇 그리퍼 암(RGA) 문제 해결

표 32: 로봇 그리퍼 암 문제 해결 표

문제/오류	가능한 원인	시정 조치
마이크로플레이트가 픽 업되지 않음	캐리어에 마이크로플레이트가 없습 니다. 그리퍼 핑거가 마이크로플레이트를 픽업할 수 없습니다.	마이크로플레이트를 캐리어에 놓습니다. 그리퍼 위치를 설정합니다. RGA 그리퍼 핑거를 세척합니다.
암이 움직이는 중 이상 한 소리	부품이 손상되거나 마모되었습니다.	"고객 지원" [▶ 188] 섹션을 참조하십시오.
편심 그리퍼 핑거 오정 렬	예비 핑거가 충돌합니다. 핑거 나사가 충분히 조여지지 않았 습니다.	편심 그리퍼 핑거를 정렬합니다. "그리퍼 핑거 정렬 점검" [▶ 176] 섹션을 참조하십시 오. "FES 그리퍼 핑거에 대한 기본 그리퍼 핑 거 정렬" [▶ 177] 섹션에 설명된 대로 토크 스크루드라이버를 사용하여 나사를 3Nm 로 조입니다.



8.2.5.1 긴 Z축이 있는 로봇 그리퍼 암(RGA-Z) 문제 해결

표 33: 긴 Z축이 있는 로봇 그리퍼 암 문제 해결 표

문제/오류	가능한 원인	시정 조치
마이크로플레이트가 픽 업되지 않음	캐리어에 마이크로플레이트가 없습 니다. 그리퍼 핑거가 마이크로플레이트를 픽업할 수 없습니다.	마이크로플레이트를 캐리어에 놓습니다. 그리퍼 위치를 설정합니다. RGA 그리퍼 핑거를 세척합니다.
	그리퍼 핑거가 미끄럽습니다.	RGA 그리퍼 핑거를 세척합니다.
암이 움직이는 중 이상 한 소리	부품이 손상되거나 마모되었습니다.	"고객 지원" [▶ 188] 섹션을 참조하십시오.

8.2.6 세척 시스템 문제 해결

표 34: 세척 시스템 문제 해결 테이블

문제/오류	가능한 원인	시정 조치
프로세스 실행 중 잘못 된 넘침 현상 및 빈 오 류	오버플로 센서가 깨끗하지 않습니 다. 센서가 연결되어 있지 않거나 결함 이 있습니다.	센서가 연결되어 있는지 확인하십시오. 센 서 제조업체의 설명서를 참조하십시오.
일반 세척 스테이션 물 넘침	폐기물 펌프에 결함이 있습니다. 폐기물 펌프 튜브가 꼬였거나 막혔 거나 손상되었습니다.	폐기물 튜브를 점검하십시오. 필요한 경우 폐기물 배관을 교체합니다. "고객 지원" [▶ 188] 섹션을 참조하십시오.
세척 스테이션을 통해 펌핑된 세정제가 없거 나 세척 시스템이 비어 있음	폐기물 펌프 튜브가 꼬였거나 막혔 거나 손상되었습니다.	폐기물 튜브를 점검하십시오. 필요한 경우 폐기물 펌프를 교체합니다. "고객 지원" [▶ 188] 섹션을 참조하십시오.
	세척 스테이션이 연결되어 있지 않 습니다. 세척 용기가 비어 있거나 없습니다.	세척 용기를 다시 채우거나 교체하십시오. 세척 스테이션(MCA)을 올바르게 연결합 니다. "세척 스테이션 연결(MCA 384)" [▶ 155] 섹션을 참조하십시오.
	폐기물 펌프에 결함이 있습니다.	폐기물 펌프를 확인합니다. 필요한 경우 폐기물 펌프를 교체합니다. "고객 지원" [▶ 188] 섹션을 참조하십시오.



문제/오류	가능한 원인	시정 조치
세척장의 과부하	폐기물 튜브가 폐기물 용기의 액체 표면 아래에 있습니다.	고정 세척 튜브의 입력부에 있는 세척용기 를 사용하십시오.
	일회용 팁 또는 이끼로 세척 스테이 션이 막혔습니다.	세척 스테이션을 세척합니다. "시스템 액체 용기 및 폐기물 용기 세척" [▶ 134] 섹션을 참조하십시오.
	폐기물 튜브가 꼬였습니다.	튜브에 꼬인 부분이 있는지 확인하십시오. "시스템 액체 용기 및 폐기물 용기의 튜브 점검" [▶ 85] 섹션을 참조하십시오.

8.2.7 Fluent ID 문제 해결

표 35: Fluent ID 문제 해결 표

문제/오류	가능한 원인	시정 조치
바코드가 판독되지 않 음	바코드 라벨이 스캐너를 향하지 않 습니다.	튜브 러너를 제거하고 바코드 레이블이 왼 쪽을 향하도록 튜브를 회전하십시오. Fluent에 튜브 러너를 다시 로드합니다.
	러너가 너무 빨리 적재되었습니다.	튜브 러너를 언로드하고 다시 천천히 로드 하십시오.
	라벨 품질이 불량합니다.	바코드를 수동으로 입력하거나 키 운영자 에게 문제를 보고하십시오.
	스캐너 창이 더럽습니다.	스캐너 창을 청소합니다. " <mark>주간 시스템</mark> 관 리" [▶ 112] 섹션을 참조하십시오.
	반사경이 더럽습니다.	리플렉터를 청소하십시오. " <mark>주간 시스템</mark> 관 리" [▶ 112] 섹션을 참조하십시오.
	바코드 유형 또는 바코드 길이가 검 사법에 대해 사전 정의되지 않았습 니다.	핵심 운영자에게 문제를 보고합니다.
튜브 존재 감지 안 됨	바코드 라벨이 튜브에서 너무 낮게 위치해 있습니다.	핵심 운영자에게 문제를 보고합니다.

8.2.8 소프트웨어 문제 해결

표 36: 소프트웨어 문제 해결 표

문제/오류	가능한 원인	시정 조치
비가시화 시 사용자 로 그인 화면이 표시되지 않습니다.	FluentControl에서 사용자 관리가 활성화되지 않았습니다.	핵심 운영자에게 문의하여 사용자 관리를 활성화하십시오.
사용자가 로그인할 수 없습니다.	암호가 올바르지 않거나 계정이 잠 겨 있습니다.	암호 또는 계정을 재설정하려면 핵심 운영 자에게 문의하십시오.



문제/오류	가능한 원인	시정 조치
모든 서비스 작업이 완 료되지 않았습니다. 각 FluentControl 시작 시 경고가 나타납니다.	기기 구성에서는 모든 예상되는 서 비스 조치들이 완료된 것으로 표시 되지 않을 수 있습니다.	"고객 지원" [▶ 188] 섹션을 참조하십시오.
터치스크린이 터치에 반응하지 않습니다.	소프트웨어 드라이버가 설치되지 않 았습니다.	설치 CD에 드라이버를 설치하고 터치스크 린을 구성하려면 컴퓨터 관리자에게 문의 하십시오.
	터치스크린 인터페이스가 잘못 구성 되었습니다.	터치스크린 드라이버 설정을 열고 터치스 크린이 올바르게 매핑되어 있는지 확인하 십시오.
터치 인터페이스가 터 치스크린에 표시되지 않습니다.	소프트웨어가 시작될 때 터치스크린 이 켜져 있지 않습니다.	기기를 켜고 소프트웨어를 다시 시작하거 나 FluentControl 구성 시스템에서 Touch Tool 설정을 확인하십시오.
FluentControl 시작 오 류.	FluentControl(SystemSW.exe)이 이미 백그라운드에서 실행 중입니다 (작업 관리자).	Task Manager(작업 관리자)를 열고 SystemSW.exe를 실행하고 FluentControl 을 다시 시작합니다. 또는 컴퓨터를 다시 시작합니다.
FluentControl이 연결 된 하드웨어 장치와 통 신하지 않습니다.	FluentControl이 하드웨어 장치와의 통신을 위해 올바르게 구성되지 않 았습니다.	하드웨어 장치의 I/O 상태를 활성화하도록 시스템을 구성하는 담당자에게 문의하십 시오.



8.3 문제 해결 활동

8.3.1 세척 스테이션 연결(MCA 384)

MCA 384 Wash System 세척 스테이션을 연결하려면 다음과 같은 절차를 따르십 시오.

- 1. 기기의 전원을 끕니다.
- 2. 세척 튜브(A)와 폐기물 튜브(B)를 세척 스테이션에 연결합니다.



3. 세척 튜브(A)와 폐기물 튜브(B)를 연결합니다.







4. 모든 튜브(C)를 MCA 384 세척 제어 장치에 연결합니다.

5. 핵심 운영자에게 연락하여 작동 준비 상태를 확인합니다.

8.3.2 실험 기구 배치

정확한 암 접근을 위해 네스트에 실험 기구를 올바르게 배치하려면 다음과 같은 절 차를 따르십시오.



1. 실험 기구(A)를 네스트(B)에 배치합니다.





2. 실험 기구를 정적 포지셔너(C)에 대고 천천히 밀어 넣습니다.

- 실험 기구에 정확하게 맞도록 대각선 슬라이딩 포지셔너(E)를 실험실 용기 쪽 또는 바깥쪽으로 밉니다.
- 세로 및 가로 슬라이딩 포지셔너(D, F)를 실험 기구 쪽으로 또는 먼 방향으로 밀어 실험 기구를 고정합니다.
- 실험 기구를 네스트에서 들어 올립니다.
 플레이트를 배치하거나 제거할 때 마찰이 없는지 확인합니다.

8.3.3 인라인 필터(Air FCA) 확인

제어 시스템은 액체의 과다 흡입으로부터 에어 FCA 피펫팅 채널을 보호하기 위해 각 채널에 설치됩니다.

- ✓ Air FCA 정기 유지 관리 방법에는 젖거나 손상되거나 잘못 배치된 필터 및 누 락된 필터를 감지하는 인라인 필터 점검이 포함됩니다.
- Air FCA 정기 유지 관리 방법을 실행하여 Air FCA 피펫팅 채널의 DiTi 콘 내부 에 있는 인라인 필터를 점검하여 정상 작동을 확인하십시오.
 오류가 발생하는 경우 인라인 필터를 변경해야 합니다. "인라인 필터(Air FCA) 변경" [▶ 157] 섹션을 참조하십시오.

8.3.4 인라인 필터(Air FCA) 변경

인라인 필터를 변경하려면 다음과 같은 절차를 따르십시오.

- ✔ 오염 제거된 일회용 팁 콘.
- ✓ 채널에서 DiTi 콘이 제거되었습니다. "DiTi 콘(에어 FCA) 제거" [▶ 159] 섹션을 참조하십시오.





- 필터 제거 도구를 사용하여 인라인 필터를 분리합니다. 필터는 공정 액체로 오 염되었을 수 있습니다.
- 3. 인라인 필터를 폐기합니다.



- DiTi 콘(A)을 알코올로 세척합니다.
 재장착 전에 DiTi 콘을 건조해야 합니다.
- 5. 깨끗하고 평평한 표면에 새 인라인 필터를 놓습니다.
- DiTi 콘에 인라인 필터를 밀어 넣습니다.
 인라인 필터가 DiTi 콘에서 돌출되지 않아야 합니다.
- 7. 핵심 운영자가 정의한 방법에 따라 인라인 필터를 확인합니다.



8.3.5 DiTi 콘(에어 FCA) 제거

DiTi 콘(에어 FCA)을 제거하려면 다음과 같은 절차를 따르십시오.

- ✓ DiTi 콘 렌치를 사용할 수 있습니다.
- 1. 기기의 전원을 끕니다.
- 2. 전면 안전 패널을 엽니다.
- 3. Z 로드를 모두 상위 위치까지 수동으로 올립니다.
- 4. 모든 Z 로드를 기기의 전면 쪽으로 움직입니다.
- 5. Z 로드를 최대한 넓게 펼칩니다.
- 6. 팁 어댑터(D)와 팁 이젝터 튜브(C)를 잡습니다.
- 7. DiTi 콘 렌치(B)를 사용하여 DiTi 콘(A)의 나사를 풉니다.



8. DiTi 콘을 조심스럽게 아래로 당깁니다.



경우에 따라 팁 이젝터 튜브(C) 또는 어댑터 실린더(B)가 DiTi 콘(A)에 여전히 부착되어 있을 수 있습니다. "DiTi 이젝터 튜브 조립(Air FCA)" [▶ 160] 섹션을 참 조하십시오.



8.3.6 DiTi 이젝터 튜브 조립(Air FCA)

다음과 같이 DiTi 이젝터 튜브(Air FCA)를 조립합니다.

- ✔ 지침에 따라 DiTi 이젝터 튜브가 제거되었습니다.
- ✔ DiTi 콘 렌치를 사용할 수 있습니다.
- 1. 밀봉 슬리브(B)를 어댑터 실린더(A)에 삽입합니다.





 조립된 실린더에 DiTi 콘(C)을 나사로 조입니다. 아래 그림과 같이 검은색 O-링이 보이지 않는지 확인합니다.



 다음과 같은 방법으로 조립된 실린더에 노치가 없는 팁 이젝터 튜브(D)를 삽입 하십시오.



8.3.7 DiTi 콘(Air FCA)를 설치합니다.

Air FCA DiTi 콘을 설치하려면 다음과 같은 절차를 따르십시오.

- ✓ DiTi 콘이 완전히 조립된 상태: "DiTi 이젝터 튜브 조립(Air FCA)" [▶ 160] 섹션을 참조하십시오.
- ✔ DiTi 콘 렌치를 사용할 수 있습니다.
- 1. 어댑터 실린더를 팁 이젝터 튜브(C)에 밀어 넣습니다.
- 2. 팁 어댑터(D)와 팁 이젝터 튜브(C)를 잡습니다.





3. DiTi 콘 렌치(B)를 사용하여 DiTi 콘(A)을 조입니다.

4. Air FCA 정기 보수 검사법을 실행합니다.



8.3.8 DiTi 옵션(FCA) 제거

DiTi 옵션을 제거하려면 다음과 같은 절차를 따르십시오.

- ✔ DiTi 콘 렌치
- 1. 기기의 전원을 끕니다.
- 2. 전면 안전 패널을 엽니다.
- 3. Z 로드를 모두 상위 위치까지 수동으로 올립니다.
- 4. 모든 Z 로드를 기기의 전면 쪽으로 움직입니다.
- 5. Z 로드를 최대한 넓게 펼칩니다.
- 6. 팁 어댑터(D)와 팁 이젝터 튜브(C)를 잡습니다.
- 7. DiTi 콘 렌치(B)를 사용하여 DiTi 콘(A)의 나사를 풉니다.



8. DiTi 콘을 조심스럽게 아래로 당깁니다.

8.3.9 DiTi 옵션(FCA) 설치

DiTi 옵션을 설치하려면 다음과 같은 절차를 따르십시오.

- 1. 어댑터 실린더(B)를 팁 이젝터 튜브(C)에 끼웁니다.
- 2. DiTi 콘(A)을 어댑터 실린더에 돌려 끼웁니다. DiTi 콘 렌치를 사용합니다.







3. 튜브가 DiTi 옵션에 단단히 부착될 때까지 플라스틱 바늘로 튜브를 밉니다.

- 4. 팁 어댑터(D)와 팁 이젝터 튜브(C)를 잡습니다.
- 5. DiTi 콘 렌치(B)를 사용하여 DiTi 콘(A)을 조입니다.



8.3.10 고정 팁 확인

고정 팁을 확인하려면 다음과 같은 절차를 따르십시오.

공지사항

피펫팅이 부정확한 문제 및 액체 감지 오류!

팁 코팅이 구부러지거나 손상되면 피펫팅이 부정확해지고 액체 감지 오류가 발생 합니다.

- 손상되거나 구부러진 팁을 사용해서는 안 됩니다.
- 1. 기기의 전원을 끕니다.



- 2. 전면 안전 패널을 엽니다.
- 3. 고정 팁을 검사합니다.
- 거울로 고정식 팁 코팅을 확인합니다.
 고정 팁이 구부러지지 않았는지 확인합니다. 고정 팁 코팅이 손상되었거나 고 정 팁이 구부러진 경우 교체해야 합니다. "고정 팁 제거" [▶ 165] 섹션을 참조하 십시오.

8.3.11 고정 팁 제거

고정 팁을 제거하려면 다음과 같은 절차를 따르십시오.

- ✔ 고정 팁을 청소했습니다. "시스템 관리 표" [▶ 107] 섹션을 참조하십시오.
- ✓ 고정 팁의 확인이 완료되었습니다. "고정 팁 확인" [▶ 164] 섹션을 참조하십시 오.



1. 기기의 전원을 끕니다.

- 2. 전면 안전 패널을 엽니다.
- 3. Z 로드(A)를 모두 상위 위치까지 수동으로 올려 줍니다.
- 4. Z 로드를 최대한 넓게 펼칩니다.
- 5. 조절식 고정 팁이 설치된 경우 4개의 팁 조정 나사를 풉니다.
- 6. 다른 손으로 잠금 너트 바로 아래 고정 팁을 잡고 잠금 너트(C)를 풉니다.
- 7. 잠금 너트(C)를 팁축 방향으로 슬라이딩하여 제거하십시오.
 - 잠금 너트와 팁 코팅은 접촉되지 않도록 하십시오.
- 팁(D)을 조정할 수 있는 경우 깨끗한 표면에 잠금 너트(C)를 뒤집어 놓고 O-링 과 와셔를 제거합니다.
- 채널에 저용량 옵션이 장착된 경우, 솔레노이드 밸브 상단의 플랜지를 풀어 Z 로드(A)를 통해 실행되는 피펫팅 튜브(E)를 해제합니다.



10 팁(D)을 위로 당겨 팁 어댑터(B)에서 특정 거리(a)까지 피펫팅 튜브(E)를 추출 . 합니다.

팁이 아니라 피펫팅 튜브에서 그립감을 개선하기 위해 마른 금강사 소재 천을 사용합니다.

8.3.12 고정 팁 설치 중



고정 팁을 설치하려면 다음과 같은 절차를 따르십시오.

 팁 어댑터에서 피펫팅 튜브를 약 25mm(1인치)만큼 조심스럽게 당깁니다. 그립감을 개선하고자 금강사 소재의 작은 천을 사용해 튜브 끝 근처를 잡습니 다.

이전에 팁을 설치한 경우 피펫팅 튜브에서 약 5mm(0.2인치)(b)만큼 자릅니다. 이때 날카로운 칼을 사용하여 직선으로 절단합니다.





- TE-PS 튜브 와이드너(G)를 사용하여 도구를 돌리면서 TE-PS 튜브 와이드너 를 튜브 내부 힌트까지 밀어 튜브 끝을 넓힙니다. 튜브가 넓혀진 상태에서 TE-PS 팁을 튜브 끝 안으로 약 4mm(0.16인치)까지 밀어 넣습니다.
- 집 끝에 잠금 너트를 올려놓습니다.
 집을 조정할 수 있는 경우(예: TE-PS) 잠금 너트를 와셔(H) 및 O-링(I) 위로 입니다.
 공지사항! 이때 팁의 약한 끝 및 코팅과 닿지 않도록 합니다.
- 4. 팁 및 피펫팅 튜브를 팁 어댑터에 삽입합니다.
- 장금 너트를 팁 어댑터에 끼우고 조입니다.
 팁을 조정할 수 있는 경우(예: TE-PS) 4개의 팁 조정 나사(K)가 데크의 X/Y 좌 표계에서 45각 위치에 놓이도록 잠금 너트를 조입니다.
- 6. 고정 팁을 세척합니다. "일과 종료" [▶ 109] 섹션을 참조하십시오.
- 7. 핵심 사용자가 정의한 대로 피펫팅 정밀도 검사를 실행합니다.





8.3.13 피어싱 팁 제거



피어싱 팁을 제거하려면 다음과 같은 절차를 따르십시오.

- ✔ 기기가 꺼져 있습니다.
- 1. 전면 안전 패널을 엽니다.
- 2. Z 로드를 모두 상위 위치까지 수동으로 올립니다.
- 3. 모든 Z 로드를 기기의 전면 쪽으로 움직입니다.
- 4. Z 로드를 최대한 넓게 펼칩니다.





되어싱 팁을 피어싱 팁 보호 장치로 덮습니다. 맨 뒤쪽 피어싱 팁부터 시작합니다.

6. 다른 손으로 잠금 너트 바로 아래 피어싱 팁을 잡고 잠금 너트를 풉니다.





 1. 팁을 위로 당겨 팁 어댑터에서 약 25mm 떨어진 위치까지 피펫팅 튜브를 당깁 니다. 당길 때 피어싱 팁의 위쪽 끝을 잡습니다.

- 8. 다른 손으로 튜브를 잡은 상태로 피어싱 팁을 튜브에서 당깁니다.
- 피어싱 팁 보호 장치를 제거하지 마십시오. 피어싱 팁이 있는 상태에서 생물학 적 폐기물 용기에 폐기합니다.





8.3.14 피어싱 팁 설치



✔ 기기가 꺼져 있습니다.

- ✔ 핵심 운영자가 수행할 수 있습니다.
- 1. 전면 안전 패널을 엽니다.
- 2. Z 로드를 모두 상위 위치까지 수동으로 올립니다.
- 3. 모든 Z 로드를 기기의 전면 쪽으로 움직입니다.
- 4. Z 로드를 최대한 넓게 펼칩니다.





되어싱 팁 포장을 엽니다.
 팁 보호 장치(F)를 제거하지 마십시오.

피어싱 팁의 장착 순서: 뒤에서 앞으로

- 6. 팁 어댑터에서 약 25 mm 정도 조심스럽게 피펫 튜브를 빼냅니다.
- 7. 피어싱 팁의 비어 있는 콘형 끝을 튜브 끝에 밀어 넣습니다



8. 피어싱 팁 및 피펫팅 튜브를 팁 어댑터에 삽입합니다.





9. 잠금 너트를 팁 어댑터에 끼우고 손으로 조입니다.

- 10 잠금 너트를 약간 여십시오. 구멍 끝 보호막을 약간 내려놓으면 구멍 끝 막대에. 접근할 수 있습니다. 아직 팁 보호 장치를 완전히 제거하지 마십시오.
- 11 피어싱 팁 개구부가 기기 전면부에 닿을 때까지 피어싱 팁을 돌립니다. 한 손으. 로 피어싱 팁을 이 방향으로 유지하고 다른 손으로 잠금 너트를 조입니다.
- 12 모든 팁 개방구가 기기의 앞쪽을 향하는지 확인합니다.







13 모든 피어싱 팁을 설치한 후 모든 팁 보호 장치를 제거합니다. 맨 뒤쪽 피어싱. 팁부터 시작합니다.

- 14 카운터를 다시 설정하려면 핵심 운영자와 연락하여 도움을 FluentControl요청 . 합니다.
- 15 QC 키트 검사 수행에 대해서는 핵심 운영자에게 문의하십시오. 를 참조하십시 . 오.
- 16 **피어싱 FCA 누출** 검사법을 실행합니다.
- 17 핵심 사용자가 정의한 대로 피펫팅 정밀도 검사(권장: QC 키트 사용)를 실행합 . 니다.



8.3.15 막힌 피어싱 팁 수축



막힌 피어싱 팁을 소프트웨어 명령으로 수축할 수 없는 경우 수동으로 제거해야 합 니다.

막힌 피어싱 팁을 수축하려면 다음과 같은 절차를 따르십시오.

- ✔ 기기가 꺼져 있습니다.
- 1. 전면 안전 패널을 엽니다.
- 2. Z 로드의 모든 수축 부분을 수동으로 상위 위치까지 올려주십시오.
- 수축된 모든 피어싱 팁을 피어싱 팁 보호 장치(A)로 덮습니다. 맨 뒤쪽 피어싱 팁부터 시작합니다.





 안정적인 표면에 놓인 막힌 팁 옆에 피어싱 팁 제거 도구를 놓고 잠금 너트 아 래에 맞춥니다.







- 5. 팁이 완전히 수축될 때까지 피어싱 팁 제거 도구의 손잡이를 돌립니다.
- 6. 손잡이를 반대 방향으로 돌리고 견인기를 약 1cm 아래로 내립니다.
- 피어싱 팁 제거 공구를 탈거합니다.
 이제 막힌 피어싱 팁이 수축됩니다.
- 8. 알코올로 피어싱 팁 제거 도구를 깨끗이 세정합니다.
- 9. 피어싱 팁을 점검하여 손상(예: 구부러진 피어싱 팁, 손상된 팁)이 있는지 확인 합니다.
- 10 피어싱 팁이 손상된 경우 교체합니다. "피어싱 팁 제거" [▶ 168] 섹션 및 "피어싱 . 팁 설치" [▶ 171] 섹션을 참조하십시오.
- 11 한 손으로 잠금 너트를 잡고 다른 손으로 팁 보호 장치를 제거하여 모든 피어싱. 팁 보호 장치를 제거합니다. 맨 뒤쪽 피어싱 팁부터 시작합니다.
- 12 알코올로 피어싱 팁 제거 도구를 깨끗이 세정합니다.

Sehen Sie dazu auch

2 "피어싱 팁 제거" } 168 2 "피어싱 팁 설치" } 171

8.3.16 그리퍼 핑거 정렬 점검

충돌 후 또는 예비 그리퍼 핑거를 구현할 때 그리퍼 핑거를 재정렬해야 할 수 있습 니다. 그리퍼 핑거가 있는 그리퍼를 사용하는 모든 암에 적용됩니다.



충돌 후 오정렬:

- 상황을 분석합니다.
- 충돌의 잠재적 원인 또는 핑거 오정렬의 원인(예: 판독기, 와셔에서 드로어 오 정렬, 잘못된 인식/배치된 호텔이나 다른 세그먼트)을 평가합니다.
- 정밀도 요건을 토대로 아래 절차를 선택합니다.
- 그리퍼 핑거가 평균 이상의 정밀도를 충족할 필요가 없는 경우 기본 정렬을 수 행합니다. "FES 그리퍼 핑거에 대한 기본 그리퍼 핑거 정렬" [▶ 177] 섹션 또는 "고정 그리퍼 핑거에 대한 기본 그리퍼 핑거 정렬" [▶ 178] 섹션을 참조하십시오.



- 2. 그리퍼 핑거가 고급 요구 사항(Z-편차 < ± 0.2 mm)을 충족해야 하는 경우, 고급 그리퍼 핑거 정렬 절차를 수행하십시오. "FES 그리퍼 핑거에 대한 고급 그 리퍼 핑거 정렬" [▶ 179] 또는 "고정 그리퍼 핑거에 대한 고급 그리퍼 핑거 정렬"
 [▶ 179] 섹션을 참조하십시오.
- 3. 그리퍼 핑거는 두 개의 나사로 장착할 수 있습니다.a) Torx 나사 M4x12의 경우 3Nm의 토크로 조입니다.



b) 육각 나사 M4x12의 경우 장력 와셔(아래 그림에 따라 위치 확인)를 함께 사 용하여 3.5Nm의 토크로 조입니다.





토크 드라이버를 사용할 수 없는 경우, 와셔를 평평하게 누르고 저항이 증가할 때까지 나사를 조입니다. 그런 다음 ¹/₁₂만큼 돌려 더 조입니다. 이는 약 3.5Nm에 해당합니다.

8.3.17 FES 그리퍼 핑거에 대한 기본 그리퍼 핑거 정렬

그리퍼 핑거가 있는 그리퍼를 사용하는 모든 암에 적용됩니다.

기본 정렬의 경우 다음과 같은 절차를 따르십시오.

- ✔ 정렬 불량이 명확하게 보입니다.
- ✔ 평균 초과 정밀도는 필요 없습니다.
- ✓ 토크 드라이버 사용 가능(3Nm 또는 3.5Nm 옵션).
 토크 드라이버를 사용할 수 없는 경우:
 토크 나사: 나사를 단단히 조이되, 과도한 힘을 가하지 마십시오.
 육각 나사: "그리퍼 핑거 정렬 점검" [▶ 176]을 참조하십시오.
- 1. 그리퍼 헤드에서 그리퍼 핑거를 제거합니다.
- 2. 그리퍼 핑거 및 FES 핑거 어댑터 사이의 나사를 풉니다.





3. 아래 그림과 같이 그리퍼 핑거를 어댑터의 상부 및 후방 정지 위치에 대고 누르 고 스크루드라이버를 사용해 3Nm 또는 3.5Nm의 토크로 나사를 조입니다.

8.3.18 고정 그리퍼 핑거에 대한 기본 그리퍼 핑거 정렬

기본 정렬의 경우 다음과 같은 절차를 따르십시오.

- ✔ 정렬 불량이 명확하게 보입니다.
- ✔ 평균 초과 정밀도는 필요 없습니다.
- ✔ 토크 스크류 드라이버가 사용 가능합니다(3 또는 3.5Nm 옵션으로).
- 1. 나사 A를 느슨하게 풀어보십시오.



 그리퍼 핑거를 그리퍼 헤드 장착부의 상부 및 후방 정지 위치에 대고 누르고 스 크루드라이버를 사용해 3Nm 또는 3.5Nm의 토크로 나사를 조입니다.



8.3.19 FES 그리퍼 핑거에 대한 고급 그리퍼 핑거 정렬

그리퍼 핑거가 있는 그리퍼를 사용하는 모든 암에 적용됩니다.

- ✓ 토크 드라이버 사용 가능(3Nm 또는 3.5Nm 옵션).
 토크 드라이버를 사용할 수 없는 경우:
 토크 나사: 나사를 단단히 조이되, 과도한 힘을 가하지 마십시오.
 육각 나사: "그리퍼 핑거 정렬 점검" [▶ 176]을 참조하십시오.
- 1. 이동 도구를 사용하여 Z 높이를 작업대 위 약 3mm 높이로 이동합니다.
- 2. 그리퍼 핑거 및 FES 핑거 어댑터 사이의 나사를 풉니다.



- FES 핑거 어댑터가 그리퍼 헤드에 단단히 연결되어 있는지 확인합니다. 핑거 어댑터는 자석에 의해 한 방향으로 고정됩니다.
- 4. 작업대 위 0mm 높이로 Z-고도를 이동하려면 Move Tool을 사용하십시오.
 마지막 0.1mm까지 천천히 움직여보십시오.
 참고: 이동 도구에 액세스하거나 작동하는 방법을 모르는 경우 핵심 운영자에 게 문의하십시오.
- 5. 그림과 같이 그리퍼 핑거를 그리퍼 헤드 및 기준 표면에 대고 누르고 3Nm 또 는 3.5Nm의 토크로 나사를 조입니다.
- 손으로 헤드를 90°, 180°, 270°로 돌려 조정을 확인합니다. 서로 다른 위치에 서 오정렬은 헤드 또는 암의 오정렬을 나타냅니다. 이 경우 FSE는 정렬을 확인 해야 합니다.

8.3.20 고정 그리퍼 핑거에 대한 고급 그리퍼 핑거 정렬

- ✓ 토크 스크류 드라이버가 사용 가능합니다(3 또는 3.5Nm 옵션으로).
- 1. 이동 도구를 사용하여 Z 높이를 약 3mm 높이로 이동합니다.
- 2. 그리퍼 핑거 및 그리퍼 헤드 사이의 나사를 풉니다.
- 이동 도구를 사용하여 Z 높이를 0mm의 높이로 이동합니다.
 마지막 0.1mm까지 천천히 움직여보십시오.
 참고: 이동 도구에 액세스하거나 작동하는 방법을 모르는 경우 핵심 운영자에 게 문의하십시오.
- 4. 그리퍼 핑거를 그리퍼 헤드 및 기준 표면에 대고 누르고 3Nm 또는 3.5Nm의 토크로 나사를 조입니다.



5. 손으로 헤드를 90°, 180°, 270°로 돌려 조정을 확인합니다. 서로 다른 위치에 서 오정렬은 헤드 또는 암의 오정렬을 나타냅니다. 이 경우 FSE는 정렬을 확인 해야 합니다.


8.3.21 세그먼트 확인

세그먼트가 닫혀 있는지 확인하십시오.



그림 69: 세그먼트 닫힘



그림 70: 세그먼트 열림

8 - 문제 해결 문제 해결 활동



8.3.22 위치 지정 핀 제거

위치 지정 핀을 제거하려면 다음과 같은 절차를 따르십시오.

공지사항

충돌 또는 프로세스 오류!

배치 핀이 느슨하여 데크 세그먼트에 요소를 잘못 배치하면 충돌 및 기타 프로세스 오류가 발생할 수 있습니다.

- 위치 지정 핀이 없는 경우에는 Fluent를 사용하지 말아야 합니다.
- ✔ 위치 지정 핀이 파손되었습니다.
- 1. 핀 제거기를 위치 지정 핀 위로 밉니다.







2. 핀 제거기 핸들을 들어 올리고 데크 세그먼트에서 핀을 당겨 꺼냅니다.

8.3.23 잠금 핀 및 위치 지정 핀 교체

공지사항

충돌 또는 프로세스 오류!

배치 핀이 느슨하여 데크 세그먼트에 요소를 잘못 배치하면 충돌 및 기타 프로세스 오류가 발생할 수 있습니다.

- 위치 지정 핀이 없는 경우에는 Fluent를 사용하지 말아야 합니다.
- ✔ 시스템 관리 테이블에 표시된 위치 지정 핀을 사용할 수 있습니다.
- 데크에서 세그먼트 제거:
 "세그먼트 제거" [▶ 78] 섹션을 참조하십시오.
- 2. 구멍(A)에 있는 새 잠금 핀(B)을 누릅니다.





3. 구멍(A)에 있는 새 위치 지정 핀(C)을 누릅니다.

데크 위에 세그먼트 배치:
 "세그먼트 배치" [▶ 77] 섹션을 참조하십시오.



9 포장, 포장 풀기, 운반, 보관 및 폐기

이 장에는 반드시 준수해야 하는 재활용 및 포장 라벨에 대한 규제 정보가 포함되 어 있습니다.

공지사항

자격이 없는 직원 및 허가받지 않은 직원에 의한 손상을 방지하십시오!

포장, 포장 풀기, 운송 및 보관은 Tecan 직원 또는 Tecan이 승인한 직원만이 수행 할 수 있습니다!

• "고객 지원" [▶ 188]을 참조하십시오.

기기 이동에 대한 자세한 내용은 "실험실 내 캐비닛에서 기기 이동하기" [▶ 116] 섹 션을 참조하십시오.

9.1 포장 라벨

정확하고 완전한 포장 표시는 잘못된 취급, 사고, 잘못된 배송, 중량 감소 및 보관 중 손상을 방지하는 데 도움이 됩니다.

표 37: 포장 기호

기호	의미	설명
	재활용	포장재는 재활용할 수 있습니다. 가정용 폐기물 로 폐기하지 마십시오. 이 포장에 사용된 재료에 대한 정보는 기호 아
		래에 나와 있습니다.
	이쪽이 윗면	포장이 상단에 화살표가 표시된 상태로 운반되 고 보관되었는지 확인합니다. 뒤집지 마십시오.
	건조한 상태 유지	운송 및 보관 중에 포장이 젖지 않도록 하십시 오.
	깨지기 쉬움	포장을 조심스럽게 다루십시오. 내부에 깨지기 쉬운 물품이 있습니다.
*	직사광선이 닿지 않는 곳에 보관	운반 및 보관 중에 포장이 열에 노출되지 않도 록 하십시오. 강한 직사광선으로부터 보호하십 시오.
	쌓지 마십시오	포장을 쌓지 마십시오. 이 포장은 추가 중량을 견딜 수 있도록 설계되지 않았습니다.





9.2 폐기

이 섹션에는 반드시 준수해야 하는 재활용에 대한 규제 정보가 포함되어 있습니다.

공지사항

해당 법률 규정에 따라 재활용하십시오!

해당 국가의 재활용 관련 법률을 준수하십시오.

9.2.1 유럽 연합 현지 요구 사항

유럽 연합 집행위원회는 전기전자폐기물처리에 관한 지침(WEEE; 2012/19/EU)을 발표했습니다.

2005년 8월 이후, 생산자는 전기 및 전자 장비의 재사용과 재활용에 대한 책임이 있습니다.



9.2.2 중국 현지 요구 사항

전자 및 전기 제품의 유해 물질 사용 제한 표시

중화인민공화국 전자산업 표준 SJ/T11364-2014 **전자 및 전기 제품에 함유된 유 해물질 사용 제한 표시**에서 요구하는 바에 따라 전자 및 전기 제품에 함유된 유해 물질 사용 제한 표시가 필요합니다.

SJ/T11364-2014에 명시된 요구 사항에 따라 중국 내에서 판매되는 모든 전자 및 전기 Tecan 제품에는 위험 물질 사용 제한 표시가 있습니다.

표시	설명
25	이 표시는 이 전자 제품에 특정 위험 물질이 포함되어 있고 환경 친화적 사용 기간 동안 안전하게 사용할 수 있음을 나 타내지만, 환경 친화적 사용 기간이 지난 후에는 재활용 시 스템에 들어가야 합니다.



9.2.3 기타 요구 사항

표시	설명
Hg	이 램프에는 수은이 함유되어 있습니다. • 해당 현지 법률에서 요구하는 대로 재활용하거나 폐기합 니다.



10 고객 지원

이 장에서는 Tecan이 문제에 대한 첫 번째 평가를 수행하는 데 필요한 파일 및 정 보를 설명합니다.

10.1 연락처

해당 지역 대리점이나 수입업체 또는 아래 주소 중 하나에 문의하십시오. 웹의 홈 페이지도 참조하십시오(www.tecan.com).

표 38: 고객 지원 담당자

국가/지역	주소	전화/팩스/이메일	
아시아	Tecan Asia Pte Ltd. 18 Boon Lay Way, #10-106 TradeHub 21 Singapore 609966 Singapore	전화 팩스 이메일	+65 6444 1886 +65 6444 1836 tecan@tecan.com.sg
Australia 뉴질랜드 태평양 제도	Tecan Australia Pty Ltd Unit 2, 475 Blackburn Road Mount Waverly VIC 3149 Australia	전화 전화 팩스 이메일	무료 전화: 1300 808 403 +61 3 9647 4100 +61 3 9647 4199 helpdesk-aus@tecan.com
Austria	Tecan Sales Austria GmbH Untersberstrasse 1a 5082 Grödig Austria	전화 팩스 이메일	+43 6246 8933 256 +43 6246 72770 helpdesk-at@tecan.com
Belgium	Tecan Benelux B.V.B.A. Mechelen Campus Schaliënhoevedreef 20A 2800 Mechelen Belgium	전화 팩스 이메일	+32 15 42 13 19 +32 15 42 16 12 tecan-be@tecan.com
중국	Tecan (Shanghai) Laboratory Equipment Co., Ltd. 1F, T 15-4, #999, Ningqiao Road, Pilot Free Trade Zone, Shanghai, PRC, 201206	전화 이메일	+86 40 0821 38 88 helpdesk-cn@tecan.com



국가/지역	주소	전화/팩스/이메일	
Denmark	Tecan Denmark, Filial af Tecan Nordic AB, Sverige Lejrvej 29 3500 Værløse Denmark	전화 이메일	+46 8 7503940 info-dk@tecan.com
France	Tecan France S.A.S.U Tour Swiss Life 1 bd Marius Vivier Merle F- 69 003 Lyon France	전화 팩스 이메일	+33 4 72 76 04 80 +33 4 72 76 04 99 helpdesk-fr@tecan.com
Germany	Tecan Deutschland GmbH Werner-von-Siemens-Straße 23 74564 Crailsheim Germany	전화 팩스 이메일	+49 1805 8322 633 또는 +49 1805 TECAN DE +49 7951 9417 92 helpdesk-de@tecan.com
Italy	Tecan Italia, S.r.I. Via Brescia, 39 20063 Cernusco Sul Naviglio (MI) Italy	전화 팩스 이메일	+39 800 11 22 91 +39 (02) 92 72 90 47 helpdesk-it@tecan.com
Japan	Tecan Japan Co., Ltd. Kawasaki Tech Center 580-16, Horikawa-cho, Saiwai-ku Kawasaki, Kanagawa 212-0013 Japan	전화 팩스 전화 이메일	+81 44 556 7311(가와사키) +81 44 556 7312(가와사키) +81(0)6305 8511(오사카) helpdesk-jp@tecan.com
Netherlands	Tecan Benelux B.V.B.A. Industrieweg 30 NL-4283 GZ Giessen Netherlands	전화 팩스 이메일	+31 20 708 4773 +31 183 44 80 67 helpdesk.benelux @tecan.com
스칸디나비아	Tecan Nordic AB Sveavägen 159, 1tr SE-113 46 Stockholm Sweden	전화 팩스 이메일	+46 8 750 39 40 +46 8 750 39 56 info@tecan.se
가산디지털 1로 149	Tecan Korea Ltd. 대한민국 서울 금천구 가산디지털 1로 149	전화 이메일	+82-2-818-3301 helpdesk-kr@tecan.com



국가/지역	주소	전화/팩스/이메일	
Spain 포르투갈	Tecan Ibérica Instrumentación S.L. C/ Lepanto 151 Bajos E-08013 Barcelona Spain	전화 이메일	+34 93 595 25 31 helpdesk-sp@tecan.com
Switzerland	Tecan Schweiz AG Seestrasse 103 8708 Männedorf Switzerland	전화 팩스 이메일	+41 44 922 82 82 +41 44 922 89 23 helpdesk-ch@tecan.com
United Kingdom	Tecan UK Ltd. Theale Court 11-13 High Street Theale, Reading, RG7 5AH United Kingdom	전화 팩스 이메일	+44 118 930 0300 +44 118 930 5671 helpdesk-uk@tecan.com
USA	Tecan US, Inc. 9401 Globe Center Drive, Suite 140, Morrisville, NC 27560 USA	전화 팩스 전화 이메일	+1 919 361 5200 +1 919 361 5201 미국 내 무료 전화: +1 800 TECAN US 또는 +1 800 832 2687 helpdesk-us@tecan.com
USA (Tecan Systems)	Tecan Systems, Inc. 2450 Zanker Road San Jose, CA 95131 USA	전화 팩스 이메일	+1 408 953 3100 무료 전화: +1 800 231 0711 +1 408 953 3101 helpdesk-sy@tecan.com



약어

ADT

Air Displacement Technology(공기 변위 기술)

Air FCA

유연한 채널 암(에어 시스템 포함)

ASM

Application Software Manual(애플리 케이션 소프트웨어 설명서)

CE

Conformité Européenne(유럽 적합 성 준수)

cLLD

Capacitive Liquid Level Detection(정전식 액체 레벨 감지)

CNS

공통 알림 시스템(Common Notification System)

DiTi

Disposable Tip(일회용 팁)

EMC

Electromagnetic Compatibility(전자 기 호환성)

ΕN

European Norm(유럽 규격)

FCA

유연한 채널 암

FES

핑거 교체 시스템

FSE

현장 서비스 엔지니어

GLP

비임상시험관리기준

HEPA

고효율 미립자 방지

IEC

International Electrotechnical Commission(국제 전기기술 위원회)

IQ

Installation Qualification(설치 자격 평가)

ISO

International Organization for Standardization(국제 표준화 기구)

LED

Light Emitting Diode(발광 다이오드)

Liquid FCA

유연한 채널 암(액체 시스템 포함)

MCA

다중 채널 암

MCA

다중 채널 암

MCH

Multiple Channel Head(다중 채널 헤 드)

MET

국가 인정 시험 연구소로서 Eurofins EE의 등록 마크

MIO

모니터링 배양기 옵션

MP

Microplate(마이크로플레이트)



NRTL

Nationally Recognized Test Laboratory(국가적으로 인정되는 검 사 실험실)

OM

Operating Manual(작동 설명서)

OQ

Operating Qualification(작동 적합성 평가)

PC

Personal Computer(개인용 컴퓨터)

PP

Polypropylene(폴리프로필렌)

rcf

상대적 원심력

RF

Radio Frequency(무선 주파수)

RGA

로봇 그리퍼 암

RGA 롱 Z

로봇 그리퍼 암 긴 높이

RGA 표준 Z

로봇 그리퍼 암 표준 높이

RUO

연구 전용

RWP

RapidWash Pump(RapidWash 펌프)

SN

일련번호

Te-Shake

Tecan 쉐이커

Te-VacS

Tecan 진공 분리기

USB

Universal Serial Bus(범용 직렬 버스)

WEEE

Waste Electrical and Electronic Equipment(전기 및 전자 장비 폐기 물)

WRC

세척 및 리필 센터