

Automated DNA clean-up.

Application Note

**AUTOMATED DNA CLEAN-UP FOR PCR AND NGS WORKFLOWS
USING OMEGA BIO-TEK'S MAG-BIND® TOTALPURE NGS KIT
ON THE FLUENT® AUTOMATION WORKSTATION**



INTRODUCTION

Next generation sequencing (NGS) technologies are increasingly being used in a variety of fields, from basic biological research to pharmacogenomics and clinical medicine. Library preparation is the first crucial step of a typical NGS workflow, and involves several DNA clean-up stages. To fully exploit the potential of NGS technologies, rapid advancements are needed in throughput and turnaround time. To meet this need, a fully automated protocol has been developed that uses Mag-Bind TotalPure NGS beads (Omega Bio-tek®) to perform the required library preparation clean-up steps on a Fluent 780 workstation. This application note demonstrates the ability of Mag-Bind TotalPure NGS beads to selectively bind DNA of different fragment lengths by altering the ratio of magnetic beads to input DNA. The performance of this high throughput solution at different bead-to-sample ratios was evaluated based on DNA recovery and quality. The results indicate that the automated workflow can efficiently purify 96 samples, with an input DNA >100 bp and up to 100 µl volume, in less than 35 minutes.

MATERIALS AND METHODS

Four 25 µl aliquots of 20x diluted 50 bp ladder were transferred to a 96-well plate in quadruplicate, and purified using Mag-Bind TotalPure NGS beads on a Fluent 780

automation workstation at bead-to-sample ratios of 0.6x, 0.8x, 1.0x and 1.2x. The Fluent workstation was configured with an Air Flexible Channel Arm™ (Air FCA), a MultiChannel Arm™ (MCA), a Robotic Gripper Arm™ (RGA), a BioShake® D30-T elm (QInstruments®) for heating and shaking, and a Magnum FLX® Enhanced Universal Magnet Plate (Alpaqua®). This worktable configuration allows for full automation of the Mag-Bind TotalPure NGS protocol, including automatic dispensing of beads as per the clean-up ratio, binding, washing and, finally, elution of the DNA in 30 µl of 10 mM Tris-HCl (pH 8.5). An aliquot of unprocessed 50 bp ladder (1:20 dilution) was included as a control for post clean-up analysis, and to evaluate the fragment sizes recovered at the four different bead-to-sample ratios.

The fragment sizes of the eluted DNA were determined using a TapeStation® 2200 (Agilent®), and compared to the unprocessed 50 bp ladder control using a High Sensitivity D1000 ScreenTape® (Agilent). TapeStation 2200 analysis software was used to estimate the percentage of DNA recovered, as well as the fragment sizes removed after clean-up at the four different ratios. The results were also compared to an unaffiliated, third-party internal report published by the Genomics and Cell Characterization Core Facility of the University of Oregon,¹ to validate the automation clean-up methodology and set-up.

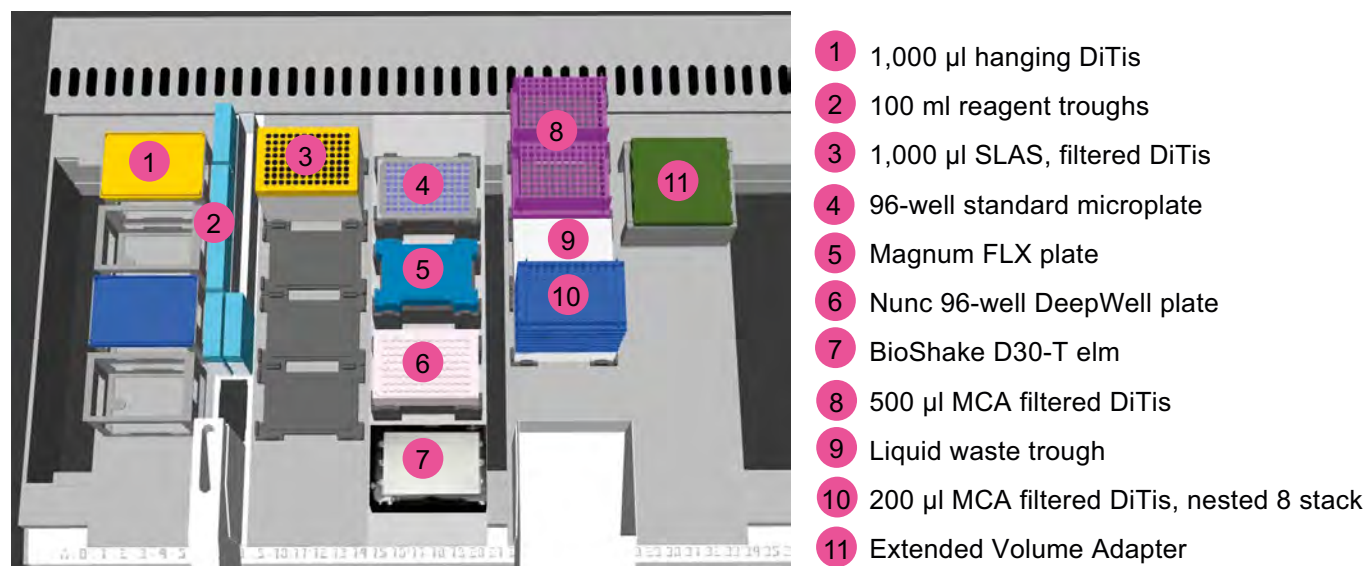


Figure 1: Fluent deck layout for extraction of 96 blood samples with 250 µl input volumes. DiTi = disposable tips.

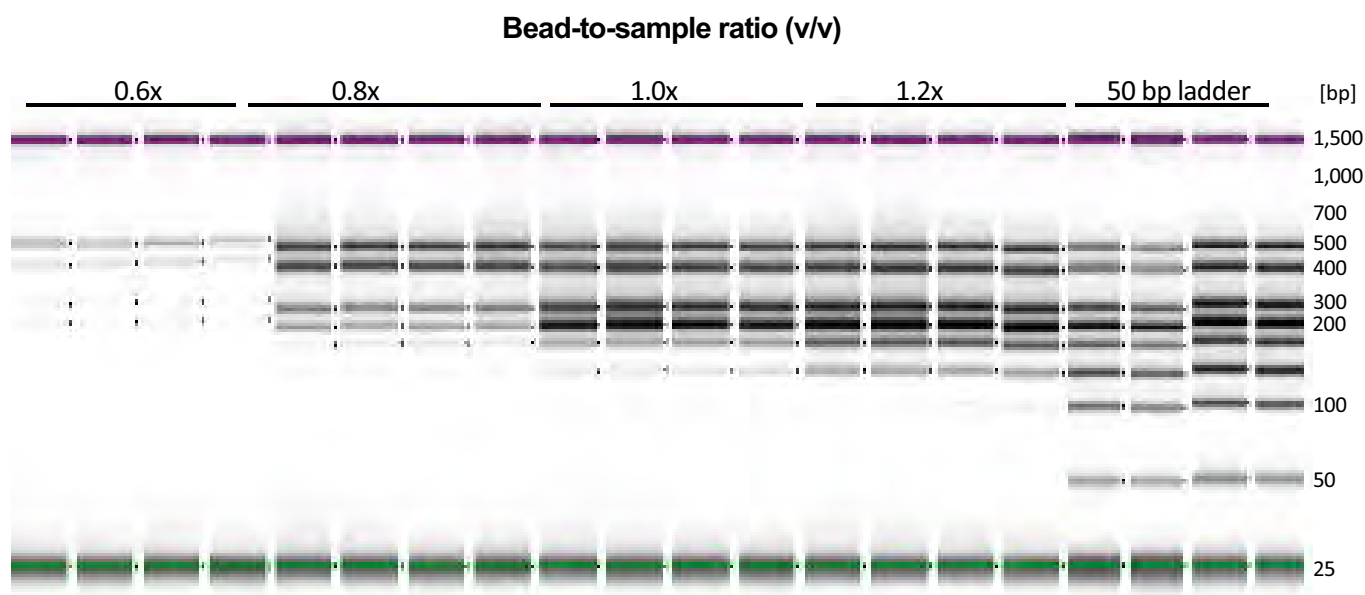


Figure 2: Fragment sizes of eluted DNA after clean-up with different bead-to-sample (v/v) ratios. TapeStation analysis was performed on 25 μ l of 20x diluted 50 bp ladder following clean-up with Mag-Bind TotalPure NGS beads on a Fluent 780 automation workstation, using unprocessed 50 bp ladder as a control.

Eight 250 μ l aliquots from the same lot of human whole blood were transferred to a 96-well deep-well plate, then placed on the Fluent workstation for automated gDNA extraction and purification using the Mag-Bind Blood & Tissue DNA HDQ 96 Kit. DNA was eluted in 100 μ l of 10 mM Tris-HCl (pH 8.5). All consumables and carriers were placed onto the Fluent workdeck (Figure 1). The workflow was fully automated, from preparation and extraction of the sample aliquots contained in the 96-well plate, to elution of the final product. Aliquots of the same lot of human whole blood were manually extracted in parallel, and the results of the two workflows compared to validate the automated methodology and instrument set-up.

RESULTS AND DATA ANALYSIS

TapeStation analysis of DNA after clean-up at various bead-to-sample ratios is shown in Figure 2. The results show that, by altering the bead-to-sample ratio, it is possible to selectively bind DNA fragment sizes. The average percentage of DNA removed at ratios of 0.6x, 0.8x, 1.0x and 1.2x (v/v) was calculated based on the unprocessed 50 bp ladder control at different fragment sizes (Figure 3).

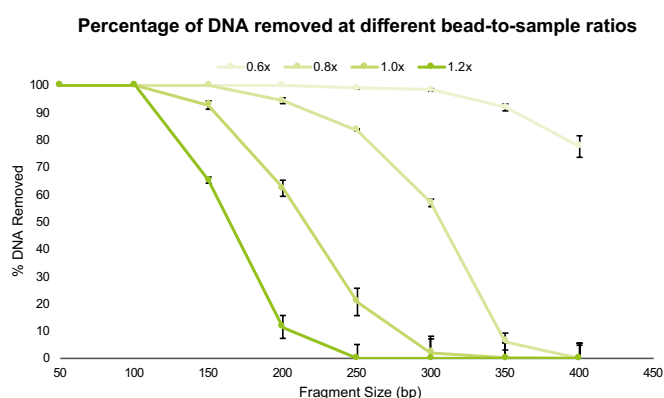


Figure 3: Average percentage of DNA removed (n = 4) at various fragment sizes using different volume ratios of Mag-Bind TotalPure NGS beads to input sample. Note: negative percentages are reported as 0 %.

The calculations were based on the concentrations of the different DNA fragments of the 50 bp ladder, as estimated by the TapeStation 2200 analysis software. The 50 bp and 100 bp fragments of the 50 bp ladder were completely removed at all the different ratios tested. The results demonstrate that higher bead volumes are capable of binding smaller fragment sizes compared to lower volumes. For example, a bead-to-sample ratio of 1.2x recovered fragments over 150 bp, whereas only fragment sizes of 300 bp and over were recovered using a 0.8x ratio. Table 1 shows that the average percentage of DNA removed using the automated protocol on a Fluent 780 workstation is in agreement with published results.¹

	100	150	200	300	400		
Bead-to-sample ratio (v/v)	0.6x	99	99	99	68	91	Published data ¹
	0.8x	98	97	96	59	17	
	1.0x	96	81	65	11	9	
	1.2x	90	51	12	0	0	
	100	150	200	300	400		
Bead-to-sample ratio (v/v)	0.6x	100	100	100	98	78	Current study
	0.8x	100	100	94	57	0	
	1.0x	100	93	62	2	0	
	1.2x	100	65	0	0	0	

Table 1: Comparison of average percentage of DNA removed (n = 4) at various bead-to-sample ratios using the automated protocol with that of published data.

¹ Note: negative percentages are reported as 0 %.

The results also show similar trends in terms of DNA binding capability at different bead-to-sample volumes. These results validate the instrument set-up and automated clean-up protocol.

SUMMARY

Integration of the Mag-Bind TotalPure NGS bead protocol with the Fluent workstation demonstrates a fully automated, high throughput DNA clean-up example for PCR and NGS applications. Using this workflow, full or partial 96-well plates containing up to 100 µl samples can be processed in less than 35 minutes. The Fluent deck configuration can easily be adapted according to the size and availability of other liquid handling arms. DNA clean-up at different bead-to-sample ratios was carried out in the same plate, demonstrating the ease and flexibility of the automated workflow. The proposed magnetic bead-based workflow is not only user-friendly, but also eliminates the need for time-consuming electrophoresis-based clean-up. The high throughput capabilities, together with the ability of the automated bead clean-up system to select the desired DNA fragment sizes, has significant potential in NGS applications.

NOTE

The third-party report published by the Genomics and Cell Characterization Core Facility of the University of Oregon evaluated and validated the performance of Mag-Bind TotalPure NGS beads at different bead-to-sample ratios (0.3x to 3x). This report serves as a guideline for the appropriate clean-up ratio(s) needed for the isolation of the DNA fragments of choice.

REFERENCES

- 1) Technical Note: Evaluation of Omega Mag-Bind® TotalPure NGS Beads for DNA Size Selection. Genomics and Cell Characterization Core Facility, University of Oregon.
https://www.omegabiotek.com/product/ngs-workflow-pcr-clean-up-mag-bind-total-pure-ngs/?mc_phishing_protection_id=55255-cbpvb4a80btra8187us0#protocols

ACKNOWLEDGEMENTS

Data provided courtesy of Kiranmai Durvasula, Evan Wolfson, Julie Baggs and Travis Butts, Omega Bio-tek, Inc, Norcross GA 30071.



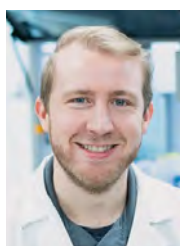
LEARN MORE

To learn more about Tecan® nucleic acid purification solutions, contact your sales representative or visit www.tecan.com/NAP.

About the authors



Kiranmai Durvasula (PhD) is a Product Manager at Omega Bio-tek, Inc., leading the efforts to develop and deliver novel nucleic acid extraction products supporting the fields of basic and translation research, diagnostics, pharmacogenomics, etc. Previously, she served as Biological Scientist at Small Business Innovation Research (SBIR/STTR) grants-funded start-up (OpenCell Technologies Inc.) developing an innovative intracellular delivery and transfection device. Kiranmai obtained her PhD in Chemical Engineering from University of Florida, Gainesville, and received postdoctoral training in cell-based therapies for treatment of insulin-dependent diabetes in Dr Athanassios Sambanis's laboratory at Georgia Institute of Technology. She has authored several peer-reviewed journal papers and is passionate about translating fundamental research platforms into practical technologies.



Evan Wolfson is a Field Applications Scientist at Omega Bio-tek, Inc., providing custom scripting solutions and automation integration for the Eastern United States and Canada. Evan has previous experience in quality systems and data management from working in

quality control for the Influenza Division of the Centers for Disease Control and Prevention, where he also acted as equipment manager and as validation engineer for the integration of a novel liquid handling solution for serological data collection and surveillance. Furthermore, he worked at Tecan as a Field Applications Engineer in the southeast United States, where he learned advanced scripting and workflow concepts that covered numerous laboratory techniques and solutions.



Adrian Cortes Sanchon (PhD) is a Product Manager at Tecan Switzerland. He studied Biochemistry at the University of Zaragoza, Spain, and did his master's in Biomedical Research at the Pompeu Fabra University, Spain, and at EMBL Heidelberg, Germany. Adrian obtained his PhD in Life Science from the University of Zurich, Switzerland, where his research focused on biochemical and -omics analysis of mistranslating systems at the Institute of Medical Microbiology, Zurich. He joined Tecan in 2022, focusing on the implementation of automated solutions for nucleic acid extraction and PCR workflows.



For Research Use Only. Not for use in diagnostic procedures.

.....
Australia +61 3 9647 4100 **Austria** +43 62 46 89 330 **Belgium** +32 15 42 13 19 **China** +86 21 220 63 206 **France** +33 4 72 76 04 80 **Germany** +49 79 51 94 170
Italy +39 02 92 44 790 **Japan** +81 44 556 73 11 **Netherlands** +31 18 34 48 17 4 **Nordic** +46 8 750 39 40 **Singapore** +65 644 41 886 **Spain** +34 93 595 25 31
Switzerland +41 44 922 89 22 **UK** +44 118 9300 300 **USA** +1 919 361 5200 **Other countries** +41 44 922 81 11

Tecan Group Ltd. makes every effort to include accurate and up-to-date information within this publication; however, it is possible that omissions or errors might have occurred. Tecan Group Ltd. cannot, therefore, make any representations or warranties, expressed or implied, as to the accuracy or completeness of the information provided in this publication. Changes in this publication can be made at any time without notice. For technical details and detailed procedures of the specifications provided in this document please contact your Tecan representative. This brochure may contain reference to applications and products which are not available in all markets. Please check with your local sales representative.

All mentioned trademarks are protected by law. In general, the trademarks and designs referenced herein are trademarks, or registered trademarks, of Tecan Group Ltd., Männedorf, Switzerland. A complete list may be found at www.tecan.com/trademarks. Product names and company names that are not contained in the list but are noted herein may be the trademarks of their respective owners.

Tecan and Fluent are registered trademarks and Air Flexible Channel Arm, MultiChannel Arm and Robotic Gripper Arm are trademarks of Tecan Group Ltd., Männedorf, Switzerland.

QInstruments and BioShake are registered trademarks of QInstruments GmbH (a BICO company), Jena, Germany. Alpaqua and Magnum FLX are registered trademarks of Alpaqua Engineering, LLC, Massachusetts, US. Agilent, TapeStation and ScreenTape are registered trademarks of Agilent Technologies, Inc., California, US. Omega Bio-tek and Mag-Bind are registered trademarks of Omega Bio-tek, Inc., Georgia, US.

© 2022, Tecan Trading AG, Switzerland, all rights reserved. For disclaimer and trademarks please visit www.tecan.com.

www.tecan.com

