

Certificate of Registration

QUALITY MANAGEMENT SYSTEM - ISO 9001:2015

This is to certify that:

IBL International GmbH
Flughafenstrasse 52A
22335 Hamburg
Germany

Holds Certificate Number:

FM 665116

and operates a Quality Management System which complies with the requirements of ISO 9001:2015 for the following scope:

Please see scope page.

For and on behalf of BSI:



Matt Page, Managing Director Assurance - UK & Ireland

Original Registration Date: 2012-07-02

Latest Revision Date: 2023-06-23

Effective Date: 2023-07-02

Expiry Date: 2026-07-01

Page: 1 of 2



...making excellence a habit.™

Certificate No: FM 665116

Registered Scope:

The design, development, manufacture and distribution of in vitro diagnostic test kits and reagents in the field of endocrinology, cancer diagnostic, hypertension, gastric diseases, diabetes, thyroid function, bone and mineral metabolism, neurodegenerative diseases, immunology, apoptosis, autoimmune diseases, infectious diseases, new born screening, allergy and food intolerance.

Design, Entwicklung, Herstellung und Vertrieb von in-vitro- diagnostischen Testkits und Reagenzien für die Anwendungsgebiete Endokrinologie, Krebsdiagnostik, Bluthochdruck, Gastroenterologie, Diabetes, Schilddrüsenfunktionen, Knochen- und Mineralstoffwechsel, Neurodegenerative Krankheiten, Immunologie, Apoptose, Autoimmunerkrankungen, Infektionskrankheiten, Neugeborenen Screening, Allergiediagnostik und Nahrungsmittelunverträglichkeit.



Original Registration Date: 2012-07-02

Latest Revision Date: 2023-06-23

Effective Date: 2023-07-02

Expiry Date: 2026-07-01

Page: 2 of 2

This certificate was issued electronically and remains the property of BSI and is bound by the conditions of contract.

An electronic certificate can be authenticated [online](#).

Printed copies can be validated at www.bsigroup.com/ClientDirectory

Information and Contact: BSI, Kitemark Court, Davy Avenue, Knowlhill, Milton Keynes MK5 8PP. Tel: + 44 345 080 9000
BSI Assurance UK Limited, registered in England under number 7805321 at 389 Chiswick High Road, London W4 4AL, UK.
A Member of the BSI Group of Companies.