

# 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:

Andrew Launn, EMEA Systems Certification Director

Original Registration Date: 2012-07-02

Effective Date: 2021-01-07

Latest Revision Date: 2021-01-07

Expiry Date: 2023-07-01



003

Page: 1 of 2

...making excellence a habit.™

Certificate No: FM 665116

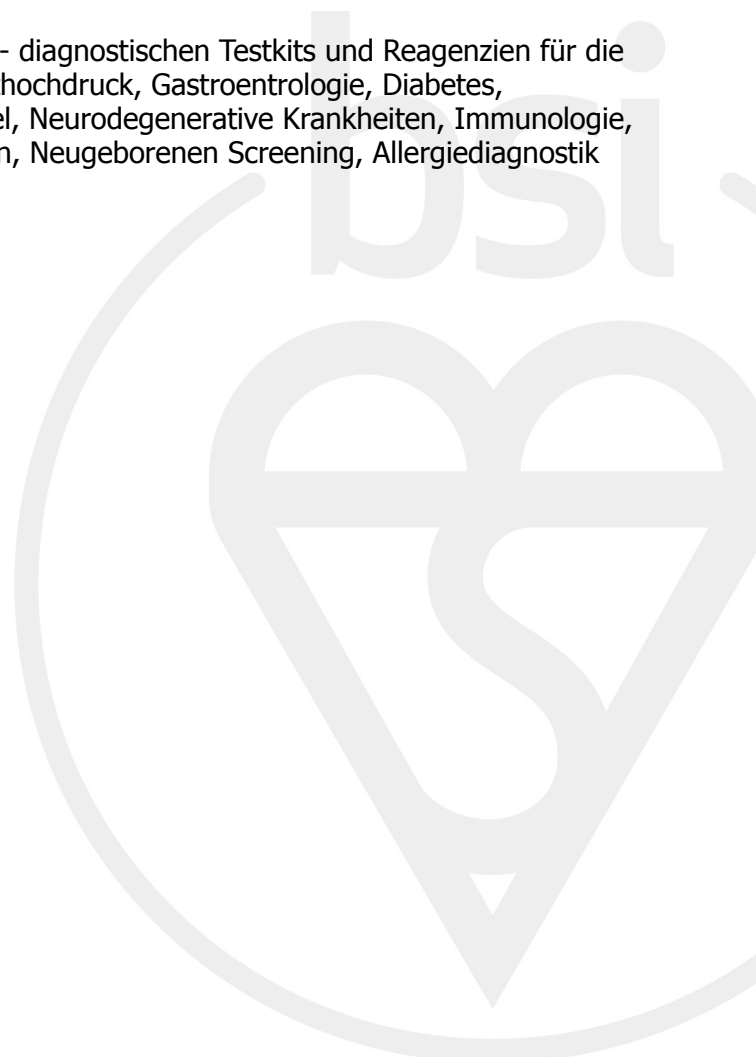
## Registered Scope:

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.

Previous certificate expires on 01.01.2021.  
Recertification audit ended 24.08.2020.

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 Nahrungsmittelnunverträglichkeit.

Der vorherigen Zertifikat ist am 01.01.2021 abgelaufen.  
Wiederzertifizierungsaudit wurde am 24.08.2020 beendet.



Original Registration Date: 2012-07-02

Latest Revision Date: 2021-01-07

Effective Date: 2021-01-07

Expiry Date: 2023-07-01

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](http://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.