



CERTIFICATE



This is to certify that



BERTHOLD TECHNOLOGIES GmbH & Co. KG

Calmbacher Straße 22
75323 Bad Wildbad
Germany

with the organizational units/sites as listed in the annex
has implemented and maintains a **Quality Management System**.

Scope:

Development and manufacturing Measuring Instruments for Process Control, Bioanalytic and
Radiation Protection

Through an audit, documented in a report, it was verified that the management system
fulfills the requirements of the following standard:

ISO 9001 : 2015

Certificate registration no.	000051 QM15
Valid from	2023-12-11
Valid until	2026-11-29
Date of certification	2023-12-11



DQS GmbH

Christian Gerling
Managing Director

DQS IS A MEMBER OF





Annex to certificate
Registration No. 000051 QM15

BERTHOLD TECHNOLOGIES GmbH & Co. KG

Calmbacher Straße 22
75323 Bad Wildbad
Germany

Location	Scope	Location	Scope
544341 BERTHOLD TECHNOLOGIES GmbH (Schweiz) Chollerstrasse 37 6300 Zug Switzerland	sales and service	533688 BERTHOLD TECHNOLOGIES (U.K.) Ltd. 6 Allied Business Centre Coldharbour Lane Harpenden AL5 4UT United Kingdom	sales and service