



CERTIFICATE



This is to certify that



BERTHOLD TECHNOLOGIES GmbH & Co. KG

Calmbacher Straße 22
75323 Bad Wildbad
Germany

has implemented and maintains an **Occupational Health and Safety 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 45001 : 2018

Certificate registration no.	000051 OHS18
Valid from	2023-12-13
Valid until	2026-11-29
Date of certification	2023-12-13



DQS GmbH

Christian Gerling
Managing Director

DQS IS A MEMBER OF



Accredited Body: DQS GmbH, August-Schanz-Straße 21, 60433 Frankfurt am Main, Germany
The validity of this certificate can only be verified by the QR-code.