

Certificate

DQS Holding has issued an IQNET recognized certificate that the organization:

BERTHOLD TECHNOLOGIES GmbH & Co. KG

**Calmbacher Straße 22
DE-75323 Bad Wildbad Germany**

has implemented and maintains an
Occupational Health and Safety Management System

for the following scope:

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

which fulfils the requirements of the following Standard:

ISO 45001:2018

Issued on: **2023-12-13**

Expires on: **2026-11-29**

Registration Number: **DE-000051 OHS18**




Alex Stoichitoiu
President of IQNet


Pedro Alves
Managing Director of IQNet

.....
This annex is only valid in connection with the original certificate number mentioned above.
.....

IQNET Members*:

AENOR Spain **AFNOR Certification** France **APCER** Portugal **CCC** Cyprus **CISQ** Italy **CQC** China **CQM** China **CQS** Czech Republic
Cro Cert Croatia **DQS Holding GmbH** Germany **EAGLE Certification Group** USA **FCAV** Brazil **FONDONORMA** Venezuela **ICONTEC**
Colombia **ICS** Bosnia and Herzegovina **INTECO** Costa Rica **IRAM** Argentina **JQA** Japan **KFQ** Korea **LSQA** Uruguay **MIRTEC** Greece
MSZT Hungary **Nemko AS** Norway **NSAI** Ireland **NYCE** México **PCBC** Poland **Quality Austria** Austria **SII** Israel **SIQ** Slovenia **SIRIM**
QAS International Malaysia **SQS** Switzerland **SRAC** Romania **TSE** Türkiye **YUQS** Serbia

* The list of IQNET Members is valid at the time of issue of this certificate. Updated information is available under www.iqnet-certification.com