



THE INTERNATIONAL CERTIFICATION NETWORK

# CERTIFICATE

DQS Holding GmbH has issued an IQNet recognized certificate that the organization

## BERTHOLD TECHNOLOGIES GmbH & Co. KG

Calmbacher Straße 22  
75323 Bad Wildbad  
Germany

with the organizational site 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

Issued on: 2018-11-21  
Expires on: 2020-12-15

This attestation is directly linked to the IQNet Partner's original certificate and  
shall not be used as a stand-alone document.

Registration number: DE-000051 QM15

Alex Stoichitoiu  
President of IQNet

Michael Drechsel  
Managing Director of  
DQS Holding GmbH



### IQNet Partners\*:

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 FCAV Brazil  
FONDONORMA Venezuela ICONTEC Colombia Inspecta Sertifointi Oy Finland INTECO Costa Rica  
IRAM Argentina JQA Japan KFQ Korea MIRTEC Greece MSZT Hungary Nemko AS Norway NSAI Ireland  
NYCE-SIGE México PCBC Poland Quality Austria Austria RR Russia SII Israel SIQ Slovenia  
SIRIM QAS International Malaysia SQS Switzerland SRAC Romania TEST St Petersburg Russia TSE Turkey YUQS Serbia  
IQNet is represented in the USA by: AFNOR Certification, CISQ, DQS Holding GmbH and NSAI Inc.

\* The list of IQNet partners is valid at the time of issue of this certificate. Updated information is available under [www.iqnet-certification.com](http://www.iqnet-certification.com)





## **Annex to IQNet Certificate Number: 000051 QM15**

### **BERTHOLD TECHNOLOGIES GmbH & Co. KG**

Calmbacher Straße 22  
75323 Bad Wildbad  
Germany

#### **Location**

#### **Scope**

**533688**

**BERTHOLD TECHNOLOGIES (U.K.) Ltd.**

sales and service

**6 Allied Business Centre**

**Coldharbour Lane**

**Harpenden**

**AL5 4UT**

**United Kingdom**