



# CERTIFICATE



This is to certify that



## 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

Certificate registration no.	000051 QM15
Valid from	2018-11-21
Valid until	2017-12-16
Date of certification	2020-12-15
	2018-11-21



**DQS GmbH**

Stefan Heinloth  
Managing Director



**Annex to certificate**  
**Registration No. 000051 QM15**

**BERTHOLD TECHNOLOGIES GmbH & Co. KG**

Calmbacher Straße 22  
75323 Bad Wildbad  
Germany

**Location**

**Scope**

<b>533688</b> <b>BERTHOLD TECHNOLOGIES (U.K.) Ltd.</b> <b>6 Allied Business Centre</b> <b>Coldharbour Lane</b> <b>Harpenden, AL5 4UT</b> <b>United Kingdom</b>	sales and service
---	-------------------