



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:

BS OHSAS 18001 : 2007

Certificate registration no.	000051 BSOH
Valid from	2017-12-16
Valid until	2020-12-15
Date of certification	2017-12-16



DQS GmbH

Frank Graichen
Managing Director

Accredited Body: DQS GmbH, August-Schanz-Straße 21, 60433 Frankfurt am Main, Germany