



CERTIFICATE



This is to certify that



BERTHOLD TECHNOLOGIES GmbH & Co. KG

Calmbacher Straße 22
75323 Bad Wildbad
Germany

has implemented and maintains an **Environmental 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 14001 : 2015

Certificate registration no.	000051 UM15
Date of revision	2018-12-21
Valid from	2017-12-16
Valid until	2020-12-15
Date of certification	2018-12-21



DQS GmbH

Stefan Heinloth
Managing Director