



Industrie Service

CERTIFICATE

The company



**Berthold Technologies
GmbH & Co.KG**
Calmbacher Strasse 22
75323 Bad Wildbad

- Germany -

has been approved according to

AD 2000-Merkblätter HP0 / HP100R

and DIN EN ISO 3834-2

The company has proven that the welding technical requirements to manufacture pressure equipment in accordance with

Pressure Equipment Directive 2014/68/EU

are met.

Regardless of this certificate, the process steps are to be kept in accordance with the selected module.

The company has following main preconditions:

- technical equipment for a proper and up-to-date manufacturing
- quality assurance which assures a correct working and testing of materials according to technical rules
- qualified supervising and manufacturing personnel.

The scope of this approval and other details are described in our report no. 600127752-HP0.

This certificate is valid until 2021-06-30

Filderstadt, 2018-08-01

TÜV SÜD Industrie Service GmbH
Anlagensicherheit - Abt. Werkstoff- und Schweißtechnik


Ulrich Haug

