



Industrie Service

# CERTIFICATE

The company



**Berthold Technologies  
GmbH & Co.KG**

**Calmbacher Strasse 22  
75323 Bad Wildbad**

- Germany -

has been approved according to

**AD 2000-Codes HP0 and HP100R**

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. 600132650-HP0.

**This certificate is valid until 2024-06-30**

Filderstadt, 2021-08-03

**TÜV SÜD Industrie Service GmbH**

Anlagensicherheit - Abt. Werkstoff- und Schweißtechnik

*Ulrich Haug*  
Ulrich Haug



TÜV SÜD Industrie Service GmbH, Abteilung Anlagensicherheit  
Gottlieb-Daimler-Straße 7, 70794 Filderstadt - Germany