



# CERTIFICATE

Nr. ISK-24-16-042

**TÜV Technische Überwachung Hessen GmbH  
Industrie Service**

we hereby certify that the company

**Sartorius Stedim Systems GmbH  
Robert – Bosch – Straße 5-7  
D-34302 Guxhagen**

as a manufacturer according to

**AD 2000-Code HP 0 / DIN EN ISO 3834-2**

was audited and accepted.

The scope of this approval and other details are described  
in our Technical Report No. 10324703 from 25. January 2017.

**Amongst others the following main preconditions are given:**

technical equipment for proper and up-to-date manufacturing and testing,  
quality procedures which assures a correct working and testing of materials  
according to the technical rules and standards,  
qualified supervisory and inspection personnel

The certificate is valid until 25. Januar 2020.

Kassel, 2.2.2017

TÜV Technische Überwachung Hessen GmbH  
Industrie Service

Department  
Systems Engineering  
Dipl.-Ing. R. Richter



Department  
Welding + Materials  
Dipl.-Ing. F. Stidronski