

CERTIFICATE

Quality system
according to directive 97/23/EC

Certificate No.: 07 202 1045 Z 0150/14/D0029

**Name and address of
manufacturer:**

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

We hereby certify that the manufacturer has established a quality system for the manufacturing of pressure equipment according to directive 97/23/EC. The manufacturer is entitled to mark the pressure equipment produced within the range of the quality system with the following mark:

CE 0045

Tested according to 97/23/EC:

production quality assurance (module D)

Test report No.:

1045 P 0150/14/D0029

Range of products:

**Bioprocess Equipment (Fermenter systems Biostat ®
and Filtration Systems Sartoflow ®)**

Place of manufacture:

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

valid until:

11/2017

Hannover, 2015-01-16



Zertifizierungsstelle für Druckgeräte
der TÜV NORD Systems
GmbH & Co. KG


Dipl.-Ing. Andreas Hoyer

Certification Body EC-Reg.No 0045

TÜV NORD Systems GmbH & Co. KG Tel. +49-(0) 551 3855 170
Große Bahnstr. 31 Fax +49-(0) 551 3855 121
D-22525 Hamburg e-mail goettingen@tuev-nord.de
Germany,

Member of



CONFÉDÉRATION EUROPÉENNE D'ORGANISMES DE CONTRÔLE