

CERTIFICATE OF REGISTRATION

This is to certify that the management system of:

Sartorius Stedim (Shanghai) Trading Co.Ltd Technical Department Validation Service

Audit Address: 2nd Floor, North Wing, Tower 1, No. 4560 Jinke Road,
Shanghai Pudong District, Shanghai

Rsgistered Address: Zone A, 3rd Floor, North Wing, Tower 1, No. 4560 Jinke
Road, China (Shanghai) Pilot Free Trade Zone

has been registered by Intertek as conforming to the requirements of:

ISO 9001:2015

The management system is applicable to:

Validation process consulting with focus on filtration and fluid management
products and filter validation testing services.

Unified Social Credit Identifier:

91310115082079882Q

Certificate Number:

111610006

Initial Certification Date:

14 July 2017

Date of Certification Decision:

26 July 2018

Issuing Date:

26 July 2018

Valid Until:

13 July 2020



014

Calin Moldovean

President, Business Assurance

Intertek Certification Limited, 10A Victory
Park, Victory Road, Derby DE24 8ZF, United
Kingdom

Intertek Certification Limited is a
UKAS accredited body under
schedule of accreditation no. 014.

