



Certificate of Approval

Accredited by Member of the International Accreditation Forum Multilateral Recognition Arrangement for Quality Management systems

KQA approves that the Quality Management System of the following company is conformed with the certification audit criteria.

Certificate No. : KQA-Q132740

Company : **Sartorius Korea Biotech Co., Ltd.**

Address : 8Fl., SolidSpace, 220, Pangyoyeok-ro, Bundang-gu, Seongnam-si, Gyeonggi-do, Republic of Korea

The standards for this certificate of conformity are as follows.

Management Systems Standards :
ISO 9001:2015 / KS Q ISO 9001:2015

Scope of Registration :
Design/development, Sales, Validation process consulting, Technical support and Engineering service of Pharmaceuticals Purification Filter, Pharmaceutical Fermentation, Bioprocess-related Products and the relevant Processes

Original Approval : 9 March 2012
Validity : 7 March 2021 ~ 6 March 2024
Date of issue : 18 September 2020 (Renewal of certificate)

Han Sang Sul 
Lead Auditor

Song Jong Cheol 
CEO

Korea Quality Assurance



Korea Quality Assurance(KQA) is accredited by Korea Accreditation Board(KAB) as ISO 9000 Certification Body/Register. (Accreditation No. : KAB-QC-05)

KQA Address : 2Fl., Hojeong Bldg., 49, Manan-ro, Manan-gu, Anyang-si, Gyeonggi-do, 14034, Republic of Korea