

Certificate of Approval

This is to certify that the Management System of:

Sartorius Stedim Chromatography Resins

48 avenue des Genottes, 95000 CERGY PONTOISE, France

has been approved by Lloyd's Register to the following standards:

ISO 9001:2015

Approval number(s): ISO 9001 – 0033286

The scope of this approval is applicable to:

Manufacture of chromatography sorbents, reagents and products.



Daniel Oliva Marcilio de Souza

Area Operations Manager - South Europe

Issued by: LRQA France SAS



Certificat d'Approbation

Nous certifions que le Système de Management de la société :

Sartorius Stedim Chromatography Resins

48 avenue des Genottes, 95000 CERGY PONTOISE, France

a été approuvé par la société Lloyd's Register selon les normes suivantes :

ISO 9001:2015

Numéro(s) d'approbation : ISO 9001 – 0033286

Le Système de Management concerne :

Réalisation de supports de chromatographie, réactifs et produits.



Daniel Oliva Marcilio de Souza

Area Operations Manager - South Europe

Emis par : LRQA France SAS

