

Sartorius Lab Instruments GmbH & Co. KG, 37070 Göttingen

Certificate of Origin

October 25th, 2023

Shipper: Sartorius Lab Instruments GmbH & Co. KG
 Company name: Sartorius Lab Instruments GmbH & Co. KG
 Company address: Otto-Brenner-Str. 20,
 37079 Göttingen, Germany
 Phone: +49 551 308 0

Based on information and declarations from raw material manufacturers regarding the material origins and production processes, and to the best of my knowledge all components in the product are of non-animal origin and are manufactured from raw materials that do not present a risk of TSE / BSE.

Part Number	Product Name
CYK-xxxx-0012	Recombinant human LIF protein

Note: "xxxx" is a placeholder for the mass in a vial in µg, e.g. "CYK-0025-0012" for a 25 µg vial of Recombinant human LIF protein.

These products are manufactured in the UK for Sartorius Lab Instruments GmbH & Co. KG.

Best regards

Sartorius Lab Instruments GmbH & Co. KG
 Head of Quality LPS



Dr. Susanne Gerighausen

Sartorius Lab Instruments GmbH & Co. KG

Otto-Brenner-Straße 20
 37079 Göttingen, Germany
 Phone +49 551 308 0
 Fax +49 551 308 3289
 www.sartorius.com

Registered Office:
 Göttingen
 Local Court of Registration:
 Amtsgericht Göttingen
 HRA 201253

The company is a limited commercial partnership

Individually liable shareholder

Sartorius Weighing Technology GmbH
 Registered Office:
 Göttingen
 Local Court of Registration:
 Amtsgericht Göttingen
 HRB 201458
 Managing Directors:
 Mario Hesse,
 Carsten Wolter