

Sartorius Lab Instruments GmbH & Co. KG, 37070 Göttingen

## Certificate of Origin

October 25<sup>th</sup>, 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-0008	Recombinant human BDNF protein

Note: "xxxx" is a placeholder for the mass in a vial in µg, e.g. "CYK-0025-0008" for a 25 µg vial of Recombinant human Activin A 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