

CERTIFICATE of ANALYSIS



Lipopolysaccharyl β -1,3-N-acetylgalactosaminyltransferase (H. influenzae F3031, recombinant)

Item No. 44171 • Batch No. 0810511 • Expires 04AUG2027

Overview

Formulation:	50 mM Tris Hydrochloric acid, pH 7.5, with 50 mM sodium chloride, and 5 mM DTT
Concentration:	4.27 mg/mL
Storage/stability:	-80°C/1yr
Biofunctional assay:	Activity confirmed by demonstrating the transfer of GalNAc (N-acetylgalactosamine) from UDP-GalNAc to globotriose (Gb3) for the synthesis of globotetraose (Gb4)
Purity:	≥90% estimated by SDS-PAGE

Reviewed and approved by: Beth Hurst

Approved date: 05AUG2026

A handwritten signature in black ink that reads "Elizabeth Hurst".

WARNING

THIS PRODUCT IS FOR RESEARCH USE - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE. IT IS THE RESPONSIBILITY OF THE PURCHASER TO DETERMINE SUITABILITY FOR OTHER APPLICATIONS.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 10/12/2018

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD

ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM