

CERTIFICATE of ANALYSIS



β -(1 $\rightarrow$ 4)-N-Acetylgalactosaminyltransferase (CgtA-II) (C. jejuni, recombinant)

Item No. 40540 • Batch No. 0820637 • Expires 16JAN2027

Overview

Formulation:	50 mM HEPES, with 20 mM magnesium dichloride and 5% glycerol
Concentration:	24.7 mg/mL
Storage/stability:	-80°C/9mo
Biofunctional assay:	Activity confirmed by demonstrating the transfer of GalNAc (N-acetylgalactosamine) from UDP-GalNAc to ganglioside GD3 for the synthesis of ganglioside GD2
Purity:	Clarified lysate

Reviewed and approved by: Beth Hurst

Approved date: 16APR2026

Beth 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