

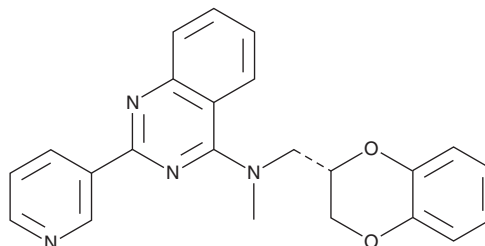
PRODUCT INFORMATION



GCase Activator 3

Item No. 45307

CAS Registry No.: 2270984-21-7
Formal Name: N-[[[(2S)-2,3-dihydro-1,4-benzodioxin-2-yl]methyl]-N-methyl-2-(3-pyridinyl)-4-quinazolinamine
Synonym: β -Glucocerebrosidase Activator 3
MF: C₂₃H₂₀N₄O₂
FW: 384.4
Purity: \geq 98%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

GCase activator 3 is supplied as a solid. A stock solution may be made by dissolving the GCase activator 3 in the solvent of choice, which should be purged with an inert gas. GCase activator 3 is slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

GCase activator 3 is an activator of β -glucocerebrosidase (GCase) with a half-maximal activation concentration (AC_{50}) value of 1.49 μ M.¹ It increases GCase protein levels and activity and reduces mutant GCase protein misfolding in Gaucher disease patient-derived fibroblasts expressing the reduction-in-function GBA1^{N370S} mutation and the loss-of-function GBA1^{84GG} mutation. GCase activator 3 (5 and 15 μ M) also increases GCase protein levels and activity in midbrain dopaminergic neurons derived from Parkinson's disease patient-derived induced pluripotent stem cells (iPSCs) and GBA1^{N370S}-expressing iPSCs, both generated from skin fibroblasts.

Reference

1. Zheng, J., Jeon, S., Jiang, W., et al. Correction to conversion of quinazoline modulators from inhibitors to activators of β -glucocerebrosidase. *J. Med. Chem.* **62**(5), 2850 (2019).

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

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, 03/11/2026

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