

PRODUCT INFORMATION



NNC-0640

Item No. 41234

CAS Registry No.: 307986-98-7

Formal Name: 4-[[[3-(4-cyclohexylphenyl)-N-2H-tetrazol-5-yl-benzamido]methyl]-N-2H-tetrazol-5-yl-benzamide

MF: C₂₉H₃₁N₇O₄S

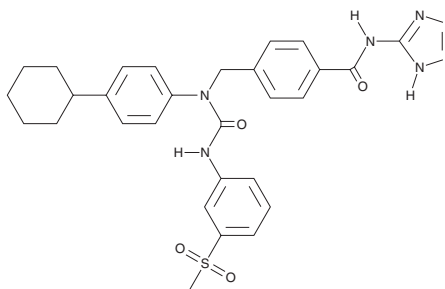
FW: 573.7

Purity: ≥98%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years



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

Laboratory Procedures

NNC-0640 is supplied as a solid. A stock solution may be made by dissolving the NNC-0640 in the solvent of choice, which should be purged with an inert gas. NNC-0640 is soluble (≥10 mg/ml) in DMSO.

Description

NNC-0640 is a negative allosteric modulator of the glucagon receptor (GCGR; K_i = 39.8 nM).¹

Reference

1. Jazayeri, A., Doré, A.S., Lamb, D., *et al.* Extra-helical binding site of a glucagon receptor antagonist. *Nature* **533(7602)**, 274-277 (2016).

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, 07/25/2024

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