

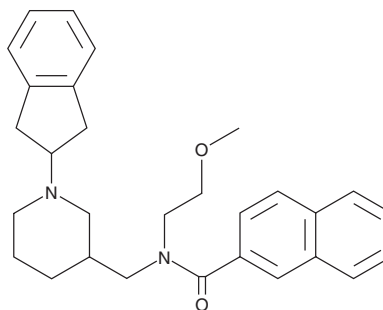
PRODUCT INFORMATION



ZINC12613047

Item No. 16975

CAS Registry No.: 1069521-64-7
Formal Name: N-[[1-(2,3-dihydro-1H-inden-2-yl)-3-piperidinyl]methyl]-N-(2-methoxyethyl)-2-naphthalenecarboxamide
MF: C₂₉H₃₄N₂O₂
FW: 442.6
Purity: ≥90%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

ZINC12613047 is supplied as a crystalline solid. A stock solution may be made by dissolving the ZINC12613047 in the solvent of choice, which should be purged with an inert gas. ZINC12613047 is soluble in organic solvents such as ethanol and DMSO. The solubility of ZINC12613047 in these solvents is approximately 5 and 10 mg/ml, respectively.

Description

ZINC12613047 is a selective butyrylcholinesterase (BChE) inhibitor with an IC₅₀ value of 21 nM for human BChE versus an IC₅₀ value of 102.1 μM for mouse AChE.¹ At 10 μM, it has been shown to inhibit Aβ₁₋₄₂ self-induced aggregation by 61.7% and to protect human neuronal SH-SY5Y cells from Aβ₁₋₄₂ peptide toxicity.¹

Reference

1. Brus, B., Kosak, U., Turk, S., *et al.* Discovery, biological evaluation, and crystal structure of a novel nanomolar selective butyrylcholinesterase inhibitor. *J. Med. Chem.* **57**(19), 8167-8179 (2014).

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, 09/09/2021

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