

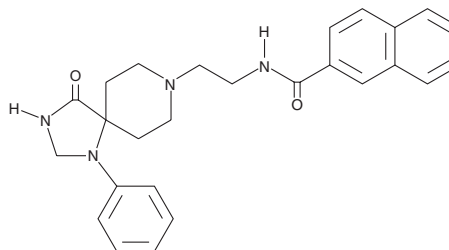
PRODUCT INFORMATION



CAY10594

Item No. 13207

CAS Registry No.: 1130067-34-3
Formal Name: N-[2-(4-oxo-1-phenyl-1,3,8-triazaspiro[4,5]dec-8-yl)ethyl]-2-naphthalenecarboxamide
MF: C₂₆H₂₈N₄O₂
FW: 428.5
Purity: ≥98%
UV/Vis.: λ_{max}: 232 nm
Supplied as: A crystalline 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

CAY10594 is supplied as a crystalline solid. A stock solution may be made by dissolving the CAY10594 in the solvent of choice, which should be purged with an inert gas. CAY10594 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of CAY10594 in these solvents is approximately 20 mg/ml.

CAY10594 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CAY10594 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. CAY10594 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Phospholipase D (PLD) is an enzyme which cleaves the head group from phospholipids, producing the second messenger phosphatidic acid. Two mammalian isoforms of PLD, PLD₁ and PLD₂, have been identified, with multiple splice variants of each. Although the two isoforms share structural and functional features, they are regulated differently and apparently subserve distinct roles. CAY10594 is a potent PLD₂ inhibitor, both *in vitro* (IC₅₀ = 140 nM) and in cells (IC₅₀ = 110 nM).¹ It is also effective as a PLD₁ inhibitor at higher concentrations (IC₅₀ = 5.1 μM *in vitro*, 1.0 μM in cells).¹ CAY10594 strongly inhibits the invasive migration of breast cancer cells in transwell assays, suggesting that PLD might be a useful target in blocking tumor cell invasion.¹

Reference

1. Scott, S.A., Selvy, P.E., Buck, J.R., *et al.* Design of isoform-selective phospholipase D inhibitors that modulate cancer cell invasiveness. *Nat. Chem. Biol.* **5**(2), 108-117 (2009).

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, 03/11/2024

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