

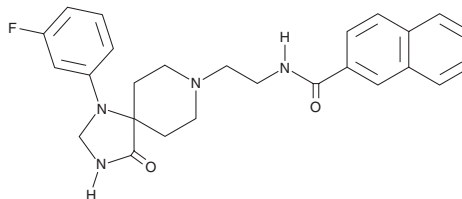
PRODUCT INFORMATION



VU0364739

Item No. 42914

CAS Registry No.: 1244639-78-8
Formal Name: N-[2-[1-(3-fluorophenyl)-4-oxo-1,3,8-triazaspiro[4.5]dec-8-yl]ethyl]-2-naphthalenecarboxamide
MF: $C_{26}H_{27}FN_4O_2$
FW: 446.5
Purity: $\geq 98\%$
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

VU0364739 is supplied as a solid. A stock solution may be made by dissolving the VU0364739 in the solvent of choice, which should be purged with an inert gas. VU0364739 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

VU0364739 is an inhibitor of phospholipase D_2 (PLD $_2$; IC $_{50}$ = 20 nM).¹ It is selective for PLD $_2$ over PLD $_1$ (IC $_{50}$ = 1,500 nM). VU0364739 (10 μM) induces apoptosis in MDA-MB-231 breast cancer cells. It decreases levels of deoxyribonucleoside triphosphates in serum-starved U87MG glioma cells when used at a concentration of 5 μM .² VU0364739 (10 μM) inhibits arachidonic acid-induced increases in the production of prostaglandin E_2 (PGE $_2$) in MC3T3-E1 osteoblasts.³

References

1. Lavieri, R.R., Scott, S.A., Selvy, P.E., et al. Design, synthesis, and biological evaluation of halogenated N-(2-(4-oxo-1-phenyl-1,3,8-triazaspiro[4.5]decan-8-yl)ethyl)benzamides: Discovery of an isoform-selective small molecule phospholipase D2 inhibitor. *J. Med. Chem.* **53**(18), 6706-6719 (2010).
2. Mathews, T.P., Hill, S., Rose, K.L., et al. Human phospholipase D activity transiently regulates pyrimidine biosynthesis in malignant gliomas. *ACS Chem. Biol.* **10**(5), 1258-1268 (2015).
3. Leis, H.J., and Windischhofer, W. Phospholipase D1 activity is crucial for cytosolic phospholipase A $_2$ -dependent prostaglandin E_2 formation in murine osteoblastic MC3T3-E1 cells. *Prostaglandins Leukot. Essent. Fatty Acids* **198-199**, 102592 (2023).

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, 03/03/2025

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