

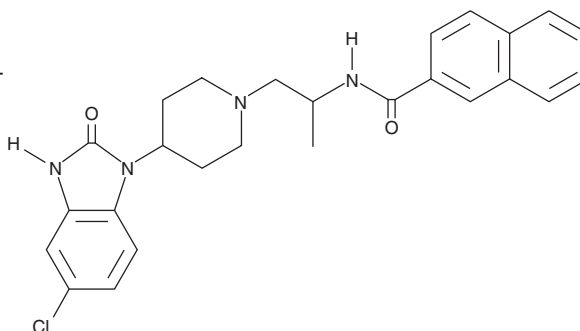
# PRODUCT INFORMATION



**VU0155069**  
Item No. 13206

**Formal Name:** N-[2-[4-(5-chloro-2,3-dihydro-2-oxo-1H-benzimidazol-1-yl)-1-piperidiny]-1-methylethyl]-2-naphthalenecarboxamide

**Synonym:** CAY10593  
**MF:** C<sub>26</sub>H<sub>27</sub>ClN<sub>4</sub>O<sub>2</sub>  
**FW:** 463.0  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 212, 232, 292 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

VU0155069 is supplied as a crystalline solid. A stock solution may be made by dissolving the VU0155069 in the solvent of choice, which should be purged with an inert gas. VU0155069 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of VU0155069 in these solvents is approximately 25 and 30 mg/ml, respectively.

VU0155069 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, VU0155069 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. VU0155069 has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMF: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, PLD<sub>1</sub> and PLD<sub>2</sub>, 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. VU0155069 is a potent and selective inhibitor of phospholipase D (PLD)1, both *in vitro* (IC<sub>50</sub> = 46 nM) and in cells (IC<sub>50</sub> = 11 nM).<sup>1</sup> It is also effective as a PLD<sub>2</sub> inhibitor at higher concentrations (IC<sub>50</sub> = 933 nM *in vitro*; 1,800 nM in cells).<sup>1</sup> VU0155069 strongly inhibits the invasive migration of several breast cancer cell lines in transwell assays, suggesting that PLD might be a useful target in blocking tumor cell invasion.<sup>1</sup>

## 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