

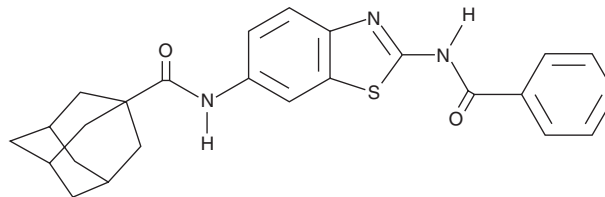
# PRODUCT INFORMATION



## NVP-231

Item No. 13858

**CAS Registry No.:** 362003-83-6  
**Formal Name:** N-[2-(benzoylamino)-6-benzothiazolyl]-tricyclo[3.3.1.1<sup>3,7</sup>]decane-1-carboxamide  
**MF:** C<sub>25</sub>H<sub>25</sub>N<sub>3</sub>O<sub>2</sub>S  
**FW:** 431.6  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 224, 316 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

NVP-231 is supplied as a crystalline solid. A stock solution may be made by dissolving the NVP-231 in the solvent of choice, which should be purged with an inert gas. NVP-231 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of NVP-231 in these solvents is approximately 1, 12, and 20 mg/ml, respectively.

### Description

NVP-231 is a potent and reversible inhibitor of ceramide kinase (IC<sub>50</sub> = 12 nM).<sup>1</sup> It is selective for ceramide kinase over SPHK1, SPHK2, DAGKa, PI3Ka, PI4Kb, VPS34, GCS, SMS-1, and CERT (IC<sub>50</sub>s = >5 mM). NVP-231 inhibits formation of ceramide-1-phosphate (C1P; IC<sub>50</sub> = 10 nM) and cell death in ceramide kinase overexpressing COS cells (COS-CerK). It also inhibits C1P formation and tube formation in murine dermal microvascular endothelial cells when used at a concentration of 100 nM.<sup>2</sup>

### References

1. Graf, C., Klumpp, M., Habig, M., *et al.* Targeting ceramide metabolism with a potent and specific ceramide kinase inhibitor. *Mol. Pharm.* **74**(4), 925-932 (2008).
2. Niwa, S., Graf, C. and Bornancin, F. Ceramide kinase deficiency impairs microendothelial cell angiogenesis in vitro. *Microvasc. Res.* **77**(3), 389-393 (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, 11/29/2022

#### 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](http://WWW.CAYMANCHEM.COM)