

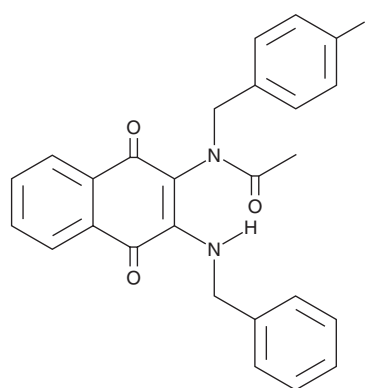
# PRODUCT INFORMATION



## RIPGBM

Item No. 27860

**CAS Registry No.:** 355406-76-7  
**Formal Name:** N-[1,4-dihydro-1,4-dioxo-3-  
[(phenylmethyl)amino]-2-naphthalenyl]-  
N-[(4-fluorophenyl)methyl]-acetamide  
**MF:** C<sub>26</sub>H<sub>21</sub>FN<sub>2</sub>O<sub>3</sub>  
**FW:** 428.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 273, 404 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

RIPGBM is supplied as a crystalline solid. A stock solution may be made by dissolving the RIPGBM in the solvent of choice, which should be purged with an inert gas. RIPGBM is soluble in organic solvents such as ethanol and DMSO.

### Description

RIPGBM is a prodrug form of the pro-apoptotic compound cRIPGBM.<sup>1</sup> It is converted to cRIPGBM in GBM-1 glioblastoma multiforme (GBM) cancer stem cells (CSCs) but not primary human lung fibroblasts (HLFs). RIPGBM induces apoptosis in GBM CSCs (EC<sub>50</sub> = ≤500 nM) and selectively reduces survival of GBM-1 cells (EC<sub>50</sub> = 0.22 μM) over human neural stem cells, primary human astrocytes, and primary HLFs (EC<sub>50</sub>s = 1.7, 2.9, and 3.5 μM, respectively). It reduces tumor growth in a GBM mouse xenograft model using patient-derived GBM-39 cells engineered to express IR fluorescent protein 720 (IRFP720) when administered at a dose of 50 mg/kg twice per day.

### Reference

1. Lucki, N.C., Villa, G.R., Vergani, N., *et al.* A cell type-selective apoptosis-inducing small molecule for the treatment of brain cancer. *Proc. Natl. Acad. Sci. U.S.A.* **116**(13), 6435-6440 (2019).

#### 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, 12/02/2022

#### 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