

# PRODUCT INFORMATION



**F1063-0967**

Item No. 36115

**CAS Registry No.:** 613225-56-2

**Formal Name:** 2-hydroxy-4-[[6-[5-[(4-methylphenyl)methylene]-4-oxo-2-thioxo-3-thiazolidinyl]-1-oxohexyl]amino]-benzoic acid

**MF:** C<sub>24</sub>H<sub>24</sub>N<sub>2</sub>O<sub>5</sub>S<sub>2</sub>

**FW:** 484.6

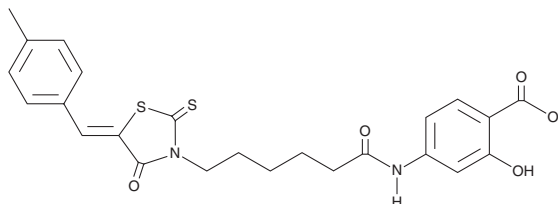
**Purity:** ≥95%

**UV/Vis.:** λ<sub>max</sub>: 275, 303, 384 nm

**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

F1063-0967 is supplied as a solid. A stock solution may be made by dissolving the F1063-0967 in the solvent of choice, which should be purged with an inert gas. F1063-0967 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of F1063-0967 in these solvents is approximately 2 mg/ml. F1063-0967 is also slightly soluble in ethanol.

## Description

F1063-0967 is an inhibitor of dual-specificity phosphatase 26 (DUSP26; IC<sub>50</sub> = 11.62 μM).<sup>1</sup> It induces IMR-32 neuroblastoma cell apoptosis (IC<sub>50</sub> = 4.13 μM).

## Reference

1. Ren, J.-X., Cheng, Z., Huang, Y.-X., *et al.* Identification of novel dual-specificity phosphatase 26 inhibitors by a hybrid virtual screening approach based on pharmacophore and molecular docking. *Biomed. Pharmacother.* **89**, 376-385 (2017).

### 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, 12/19/2022

## 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)