

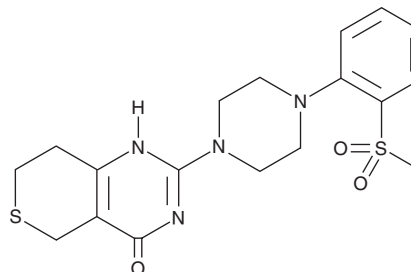
# PRODUCT INFORMATION



## G244-LM

Item No. 21876

**CAS Registry No.:** 1563007-08-8  
**Formal Name:** 3,5,7,8-tetrahydro-2-[4-[2-(methylsulfonyl)phenyl]-1-piperazinyl]-4H-thiopyrano[4,3-d]pyrimidin-4-one  
**MF:** C<sub>18</sub>H<sub>22</sub>N<sub>4</sub>O<sub>3</sub>S<sub>2</sub>  
**FW:** 406.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 213, 235, 300 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

G244-LM is supplied as a crystalline solid. A stock solution may be made by dissolving the G244-LM in the solvent of choice, which should be purged with an inert gas. G244-LM is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of G244-LM in these solvents is approximately 0.25, 5, and 2 mg/ml, respectively.

G244-LM is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, G244-LM should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. G244-LM has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

G244-LM is an inhibitor of Wnt signaling.<sup>1</sup> It inhibits Wnt3a-induced signaling in HEK293 human and 10T1/2 mouse cells in a reporter assay. It also inhibits Wnt signaling and colony formation in COLO-320DM and SW403, but not HCT-15 or DLD-1, colorectal cancer cells when used at a concentration of 0.2 μM for seven to 13 days. G244-LM inhibits growth of mouse intestinal organoids (IC<sub>50</sub> = 0.11 μM) and organoid spheroids cultured from small intestine adenomas isolated from Apc<sup>Min</sup> mice, which exhibit constitutively activated Wnt/β-catenin signaling.

### Reference

1. Lau, T., Chan, E., Callow, M., *et al.* A novel tankyrase small-molecule inhibitor suppresses APC mutation-driven colorectal tumor growth. *Cancer Res.* **73**(10), 3132-3144 (2013).

#### 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, 10/03/2022

#### 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)