

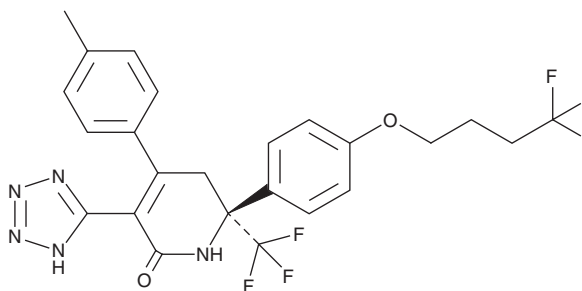
# PRODUCT INFORMATION



**BMS 963272**

Item No. 34981

**CAS Registry No.:** 1441057-15-3  
**Formal Name:** (6S)-5,6-dihydro-4-(4-methylphenyl)-3-(2H-tetrazol-5-yl)-6-[4-(4,4,4-trifluorobutoxy)phenyl]-6-(trifluoromethyl)-2(1H)-pyridinone  
**MF:** C<sub>24</sub>H<sub>21</sub>F<sub>6</sub>N<sub>5</sub>O<sub>2</sub>  
**FW:** 525.4  
**Purity:** ≥98%  
**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

BMS 963272 is supplied as a solid. A stock solution may be made by dissolving the BMS 963272 in the solvent of choice, which should be purged with an inert gas. BMS 963272 is soluble in the organic solvent DMSO.

## Description

BMS 963272 is an inhibitor of monoacylglycerol acyltransferase 2 (MGAT-2; IC<sub>50</sub> = 7.1 nM).<sup>1</sup> It is selective for MGAT-2 over MGAT-3, acyl-CoA wax alcohol acyltransferase 2 (AWAT2), and diacylglycerol acyltransferase 2 (DGAT-2; IC<sub>50</sub>s = >33 μM for all). *In vivo*, BMS 963272 (30 mg/kg) reduces body weight gain and food intake in a wild-type but not *Mogat2*<sup>-/-</sup> mouse model of high-fat diet-induced obesity. BMS 963272 (0.3 and 3 mg/kg) reduces plasma alanine transaminase (ALT), aspartate aminotransferase (AST), and bile acid levels and hepatic fibrosis in a mouse model of non-alcoholic steatohepatitis (NASH) induced by a choline-deficient and high-fat diet (CDAHFD).<sup>2</sup>

## References

1. Turdi, H., Chao, H., Hangeland, J.J., *et al.* Screening hit to clinical candidate: Discovery of BMS 963272, a potent, selective MGAT2 inhibitor for the treatment of metabolic disorders. *J. Med. Chem.* **64**(19), 14773-14792 (2021).
2. Cheng, D., Zinker, B.A., Luo, Y., *et al.* MGAT2 inhibitor decreases liver fibrosis and inflammation in murine NASH models and reduces body weight in human adults with obesity. *Cell Metab.* **34**(11), 1732-1748 (2022).

### 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, 12/06/2023

## 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