

# PRODUCT INFORMATION



**MP07-66**

Item No. 42337

**CAS Registry No.:** 1938056-90-6  
**Formal Name:** N-(2,2-diethoxyethyl)-4-(hexyloxy)-benzenemethanamine

**MF:** C<sub>19</sub>H<sub>33</sub>NO<sub>3</sub>

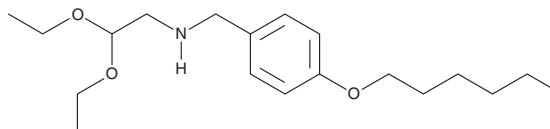
**FW:** 323.5

**Purity:** ≥98%

**Supplied as:** A solution in ethanol

**Storage:** -20°C

**Stability:** ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

## Description

MP07-66 is an inhibitor of the protein-protein interaction between SET and protein phosphatase 2A (PP2A) and a derivative of FTY720 (fingolimod; Item No. 11975).<sup>1</sup> It increases PP2A activity in isolated human B cell chronic lymphocytic leukemia (B-CLL) cells when used at concentrations of 3 and 8 μM. Unlike FTY720, MP07-66 (8 μM) does not decrease plasma membrane levels of sphingosine-1-phosphate receptor 1 (S1P<sub>1</sub>) in isolated human B-CLL cells. MP07-66 induces apoptosis in isolated human B-CLL cells when used alone or in combination with dasatinib (Item No. 11498).

## Reference

1. Zonta, F., Pagano, M.A., Trentin, L., *et al.* Lyn sustains oncogenic signaling in chronic lymphocytic leukemia by strengthening SET-mediated inhibition of PP2A. *Blood* **125**(24), 3747-3755 (2015).

### 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, 05/15/2025

## 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)