

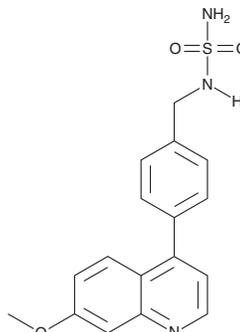
# PRODUCT INFORMATION



## ENPP1-IN-1

Item No. 31764

**CAS Registry No.:** 2289728-58-9  
**Formal Name:** N-[[4-(7-methoxy-4-quinolinyl)phenyl]methyl]-sulfamide  
**Synonyms:** N-[4-(7-Methoxyquinolin-4-yl)benzyl]sulfamide, MV 658, UUN28589  
**MF:** C<sub>17</sub>H<sub>17</sub>N<sub>3</sub>O<sub>3</sub>S  
**FW:** 343.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 234 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

ENPP1-IN-1 is supplied as a crystalline solid. A stock solution may be made by dissolving the ENPP1-IN-1 in the solvent of choice, which should be purged with an inert gas. ENPP1-IN-1 is soluble in DMSO.

### Description

ENPP1-IN-1 is an inhibitor of ectonucleotide pyrophosphatase/phosphodiesterase 1 (ENPP1).<sup>1</sup> It inhibits the hydrolysis of 2'3'-cGAMP (Item No. 19887) and the ATP analog *p*-nitrophenyl 5'-adenosine monophosphate (AMP-*p*NP; K<sub>i</sub>s = ≤100 and 100-1,000 nM, respectively, in enzyme assays). ENPP1-IN-1 enhances transcription of the gene encoding IFN-β in 2'3'-cGAMP-stimulated human foreskin fibroblast cells (HFF-1) in a concentration-dependent manner.

### Reference

1. Gallatin, W.M., Odingo, J., Dietsch, G.N., *et al.* Ectonucleotide pyrophosphatase-phosphodiesterase 1 (ENPP-1) inhibitors and uses thereof. *Mavupharma, Inc. WO2019/046778 A1* (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, 11/17/2022

#### 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