

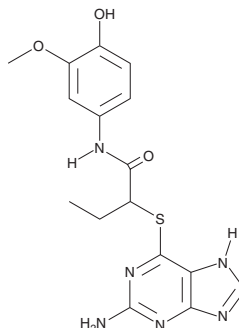
PRODUCT INFORMATION



ENPP1 Inhibitor 43

Item No. 36836

CAS Registry No.: 2631703-41-6
Formal Name: 2-[(2-amino-9H-purin-6-yl)thio]-N-(4-hydroxy-3-methoxyphenyl)-butanamide
MF: C₁₆H₁₈N₆O₃S
FW: 374.4
Purity: ≥98%
UV/Vis.: λ_{max}: 216, 242, 310 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

ENPP1 inhibitor 43 is supplied as a crystalline solid. A stock solution may be made by dissolving the ENPP1 inhibitor 43 in the solvent of choice, which should be purged with an inert gas. ENPP1 inhibitor 43 is soluble in the organic solvent DMSO at a concentration of approximately 1 mg/ml.

Description

ENPP1 inhibitor 43 is an inhibitor of ectonucleotide pyrophosphatase/phosphodiesterase family member 1 (ENPP1; K_i = 41 nM).¹ It inhibits tumor growth in an LLC1 murine lung cancer model when administered orally or intravenously at doses of 10 and 100 mg/kg, respectively.

Reference

1. Gangar, M., Goyal, S., Raykar, D., *et al.* Design, synthesis and biological evaluation studies of novel small molecule ENPP1 inhibitors for cancer immunotherapy. *Bioorg. Chem.* **119**, 105549 (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, 09/09/2022

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