

PRODUCT INFORMATION



MLK3 Inhibitor

Item No. 42887

Formal Name: N⁴-(3-(methylsulfonyl)benzyl)-5-(trifluoromethyl)-N²-(4-(trifluoromethyl)phenyl)pyrimidine-2,4-diamine

Synonyms: Mixed-lineage Kinase 3-IN-1, Mixed-lineage Kinase 3 Inhibitor, Mixed-lineage Kinase 3 Inhibitor 1, MLK3-IN-1, MLK3 Inhibitor 1

MF: C₂₀H₁₆F₆N₄O₂S

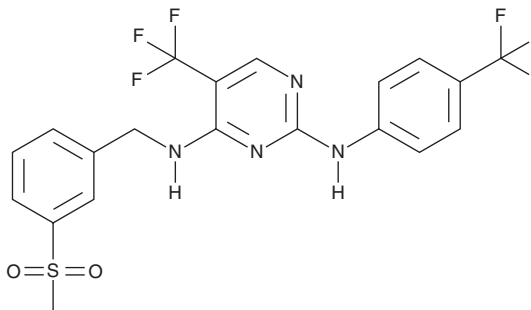
FW: 490.4

Purity: ≥98%

Supplied as: A neat 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

MLK3 inhibitor is supplied as a neat solid. A stock solution may be made by dissolving the MLK3 inhibitor in the solvent of choice, which should be purged with an inert gas. MLK3 inhibitor is soluble (≥10 mg/ml) in DMSO.

Description

MLK3 inhibitor is an inhibitor of mixed-lineage kinase 3 (MLK3; IC₅₀ = <0.001 μM).¹ It is selective for MLK3 over focal adhesion kinase (FAK; IC₅₀ = 15.5 μM).

Reference

1. Sander, P., Schwalm, M.P., Krämer, A., *et al.* Design, synthesis, and biochemical evaluation of novel MLK3 inhibitors: A target hopping example. *J. Med. Chem.* **68**(1), 674-694 (2025).

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, 05/20/2025

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