

PRODUCT INFORMATION



TC 13172

Item No. 31034

CAS Registry No.: 2093393-05-4
Formal Name: 3,7-dihydro-3-[3-(3-hydroxyphenyl)-2-propyn-1-yl]-1,7-dimethyl-8-(methylsulfonyl)-1H-purine-2,6-dione

MF: C₁₇H₁₆N₄O₅S

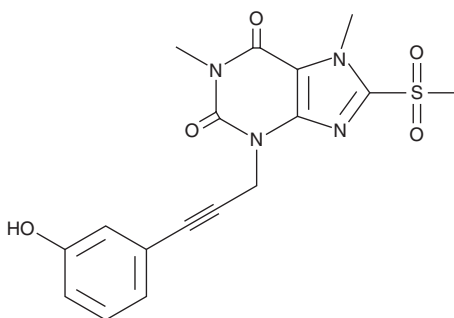
FW: 388.4

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

TC 13172 is supplied as a solid. A stock solution may be made by dissolving the TC 13172 in the solvent of choice, which should be purged with an inert gas. TC 13172 is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

TC 13172 is a covalent inhibitor of mixed lineage kinase domain-like protein (MLKL).¹ It is selective for MLKL over receptor-interacting serine/threonine kinase 1 (RIPK1) and RIPK3 at 10 μM. TC 13172 inhibits necroptosis induced by TSZ, which is a combination of TNF-α, a Smac mimetic, and Z-VAD-FMK, in HT-29 cells (EC₅₀ = 2 nM). It inhibits TSZ-induced oligomerization of MLKL and translocation of MLKL to the plasma membrane in HT-29 cells when used at a concentration of 100 nM.

Reference

1. Yan, B., Liu, L., Huang, S., *et al.* Discovery of a new class of highly potent necroptosis inhibitors targeting the mixed lineage kinase domain-like protein. *Chem. Commun. (Camb.)* **53**(26), 3637-3640 (2017).

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, 10/25/2024

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