

PRODUCT INFORMATION

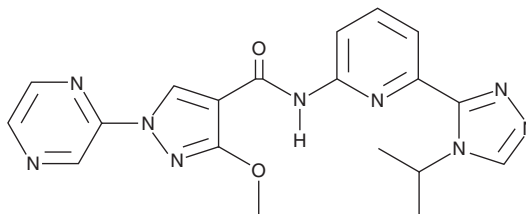


CAY10756

Item No. 30382

CAS Registry No.: 2411382-24-4
Formal Name: 3-methoxy-N-[6-[4-(1-methylethyl)-4H-1,2,4-triazol-3-yl]-2-pyridinyl]-1-(2-pyrazinyl)-1H-pyrazole-4-carboxamide

MF: C₁₉H₁₉N₉O₂
FW: 405.4
Purity: ≥98%
UV/Vis.: λ_{max}: 275, 318 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

CAY10756 is supplied as a crystalline solid. A stock solution may be made by dissolving the CAY10756 in the solvent of choice, which should be purged with an inert gas. CAY10756 is soluble in the organic solvent chloroform at a concentration of approximately 10 mg/ml.

Description

CAY10756 is a brain-penetrant inhibitor of apoptosis signal-regulating kinase 1 (ASK1; IC₅₀ = 21 nM).¹ It inhibits autophosphorylation of ASK1 in HEK293T cells (IC₅₀ = 138 nM).

Reference

1. Xin, Z., Himmelbauer, K., Jones, J.H., *et al.* Discovery of CNS-penetrant apoptosis signal-regulating kinase 1 (ASK1) inhibitors. *ACS Med. Chem. Lett.* **11**(4), 485-490 (2020).

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, 03/21/2024

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