

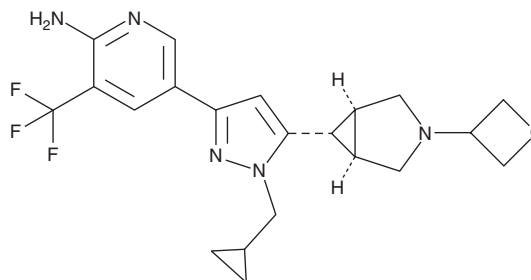
PRODUCT INFORMATION



GNE-8505

Item No. 44006

CAS Registry No.: 1620573-48-9
Formal Name: 5-[1-(cyclopropylmethyl)-5-[(1 α ,5 α ,6 α)-3-(3-oxetanyl)-3-azabicyclo[3.1.0]hex-6-yl]-1H-pyrazol-3-yl]-3-(trifluoromethyl)-2-pyridinamine
MF: C₂₁H₂₄F₃N₅O
FW: 419.4
Purity: \geq 98%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

GNE-8505 is supplied as a solid. A stock solution may be made by dissolving the GNE-8505 in the solvent of choice, which should be purged with an inert gas. GNE-8505 is slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

GNE-8505 is an inhibitor of dual leucine zipper kinase (DLK).¹ It decreases the levels of phosphorylated JNK in retinal lysates prepared from a mouse model of optic nerve crush when administered at doses of 3, 7, or 18 mg/kg.

Reference

1. Le Pichon, C.E., Meilandt, W.J., Dominguez, S., *et al.* Loss of dual leucine zipper kinase signaling is protective in animal models of neurodegenerative disease. *Sci. Transl. Med.* **9(403)**, eaag0394 (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, 09/02/2025

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