

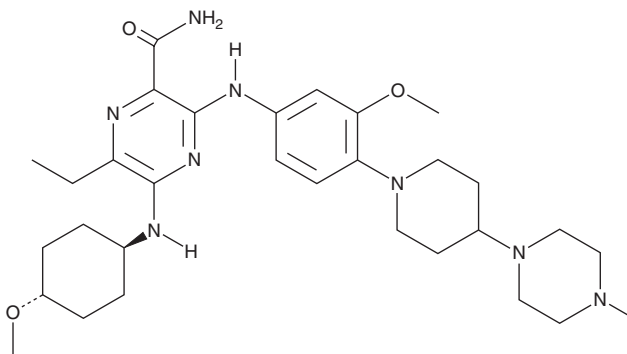
PRODUCT INFORMATION



EML4-ALK Kinase Inhibitor 1

Item No. 45387

CAS Registry No.: 1373409-08-5
Formal Name: 6-ethyl-5-[(*trans*-4-methoxycyclohexyl)amino]-3-[[[3-methoxy-4-[4-(4-methyl-1-piperazinyl)-1-piperidinyl]phenyl]amino]-2-pyrazinecarboxamide
MF: C₃₁H₄₈N₈O₃
FW: 580.8
Purity: ≥95%
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

EML4-ALK kinase inhibitor 1 is supplied as a solid. A stock solution may be made by dissolving the EML4-ALK kinase inhibitor 1 in the solvent of choice, which should be purged with an inert gas. EML4-ALK kinase inhibitor 1 is slightly soluble (0.1-1 mg/ml) in ethanol.

Description

EML4-ALK kinase inhibitor 1 is an inhibitor of the anaplastic lymphoma kinase (ALK) fusion protein EML4-ALK (IC₅₀ = 1 nM).¹ It inhibits the growth of Ba/F3 cells expressing EML4-ALK (IC₅₀ = 9.7 nM). *In vivo*, EML4-ALK kinase inhibitor 1 (10 mg/kg) induces tumor regression in mice xenografted with NIH3T3 cells expressing EML4-ALK.

Reference

1. Iikubo, K., Kurosawa, K., Matsuya, T., *et al.* Synthesis and structure-activity relationships of pyrazine-2-carboxamide derivatives as novel echinoderm microtubule-associated protein-like 4 (EML4)-anaplastic lymphoma kinase (ALK) inhibitors. *Bioorg. Med. Chem.* **27(8)**, 1683-1692 (2019).

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, 04/14/2026

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