

PRODUCT INFORMATION

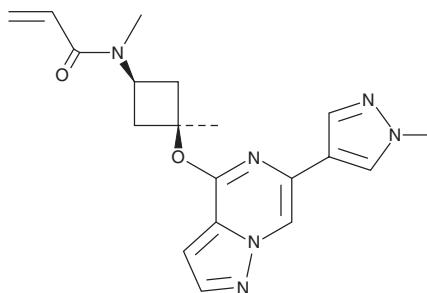


BIIB129

Item No. 43712

CAS Registry No.: 2770960-52-4
Formal Name: N-methyl-N-[cis-3-methyl-3-[[6-(1-methyl-1Hpyrazol-4-yl)pyrazolo[1,5-a]pyrazin-4-yl]oxy]cyclobutyl]-2-propenamide

MF: C₁₉H₂₂N₆O₂
FW: 366.4
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

BIIB129 is supplied as a solid. A stock solution may be made by dissolving the BIIB129 in the solvent of choice, which should be purged with an inert gas. BIIB129 is soluble (≥10 mg/ml) in DMSO. BIIB129 is sparingly soluble (1-10 mg/ml) in ethanol.

Description

BIIB129 is a covalent inhibitor of Bruton's tyrosine kinase (BTK).¹ It is selective for BTK over a panel of 403 additional kinases at 1 μM. BIIB129 inhibits CD69 activation in isolated human whole blood (IC₅₀ = 7.9 nM). *In vivo*, BIIB129 inhibits anti-myelin oligodendrocyte glycoprotein (MOG) antibody-induced microglia activation in mice (ED₅₀ = 1.5 mg/kg).

Reference

- Himmelbauer, M.K., Bajrami, B., Basile, R., et al. Discovery and preclinical characterization of BIIB129, a covalent, selective, and brain-penetrant BTK inhibitor for the treatment of multiple sclerosis. *J. Med. Chem.* **67**(10), 8122-8140 (2024).

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, 07/25/2025

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