

PRODUCT INFORMATION



BIIB068

Item No. 31619

CAS Registry No.: 1798787-27-5
Formal Name: 3-(1-methylethoxy)-N-[[2-methyl-4-[[2-[(1-methyl-1H-pyrazol-4-yl)amino]-4-pyrimidinyl]phenyl)methyl]-1-azetidinecarboxamide

MF: C₂₃H₂₉N₇O₂

FW: 435.5

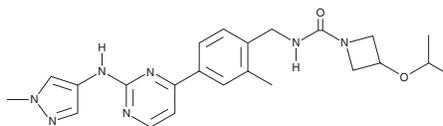
Purity: ≥98%

UV/Vis.: λ_{max}: 274 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

BIIB068 is supplied as a crystalline solid. A stock solution may be made by dissolving the BIIB068 in the solvent of choice, which should be purged with an inert gas. BIIB068 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of BIIB068 is approximately 10 mg/ml in ethanol and 30 mg/ml in DMSO and DMF.

BIIB068 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, BIIB068 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. BIIB068 has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

BIIB068 is an inhibitor of Bruton's tyrosine kinase (BTK; IC₅₀ = <10 nM).^{1,2}

References

1. Li, C., Franklin, L., Chen, R., *et al.* Process development and large-scale synthesis of BTK inhibitor BIIB068. *Org. Process Res. Dev.* **24(6)**, 1199-1206 (2020).
2. Hopkins, B.T., Ma, B., Chan, T.R., *et al.* Biaryl compounds useful for the treatment of human diseases in oncology, neurology, and immunology. *Biogen Idec and Sunesis Pharmaceuticals*. **WO2015089337A1** (2015).

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