

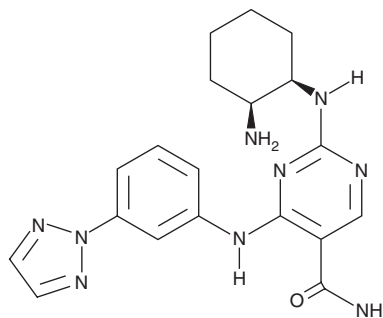
PRODUCT INFORMATION



BIIB-057

Item No. 17652

CAS Registry No.: 1370261-96-3
Formal Name: 2-[[[(1R,2S)-2-aminocyclohexyl]amino]-4-[[3-(2H-1,2,3-triazol-2-yl)phenyl]amino]-5-pyrimidinecarboxamide
Synonyms: P505-15, PRT062607, PRT2607
MF: C₁₉H₂₃N₉O
FW: 393.5
Purity: ≥98%
UV/Vis.: λ_{max}: 264 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

BIIB-057 is supplied as a crystalline solid. A stock solution may be made by dissolving the BIIB-057 in the solvent of choice, which should be purged with an inert gas. BIIB-057 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of BIIB-057 in these solvents is approximately 10, 25, and 50 mg/ml, respectively.

BIIB-057 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, BIIB-057 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. BIIB-057 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

BIIB-057 is a potent inhibitor of the non-receptor tyrosine kinase Syk (IC₅₀ = 1 nM).¹ It displays at least 80-fold selectivity for Syk over other kinases. BIIB-057 blocks B cell receptor-mediated cell signaling and activation in whole blood (IC_{50S} = 0.27 and 0.28 μM, respectively), as well as Fcε receptor 1-mediated basophil degranulation (IC₅₀ = 0.15 μM).¹ It antagonizes chemokine production, cell migration, and survival of chronic lymphocytic leukemia (CLL) cells after B cell receptor activation and synergistically enhances the action of fludarabine (Item No. 14128) in killing CLL cells.^{2,3} BIIB-057 is orally bioavailable, as it produces dose-dependent anti-inflammatory activity in two rodent models of rheumatoid arthritis.¹ It also prevents splenomegaly and inhibits non-Hodgkin lymphoma tumor growth in a xenograft model.³

References

1. Coffey, G., DeGuzman, F., Inagaki, M., *et al.* Specific inhibition of spleen tyrosine kinase suppresses leukocyte immune function and inflammation in animal models of rheumatoid arthritis. *J. Pharmacol. Exp. Ther.* **340**(2), 350-359 (2012).
2. Hoellenriegel, J., Coffey, G.P., Sinha, U., *et al.* Selective, novel spleen tyrosine kinase (Syk) inhibitors suppress chronic lymphocytic leukemia B-cell activation and migration. *Leukemia* **26**(7), 1576-1583 (2012).
3. Spurgeon, S.E., Coffey, G., Fletcher, L.B., *et al.* The selective Syk inhibitor P505-15 (PRT062607) inhibits B cell signaling and function in vitro and in vivo and augments the activity of fludarabine in chronic lymphocytic leukemia. *J. Pharmacol. Exp. Ther.* **344**(2), 378-387 (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, 11/29/2022

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