

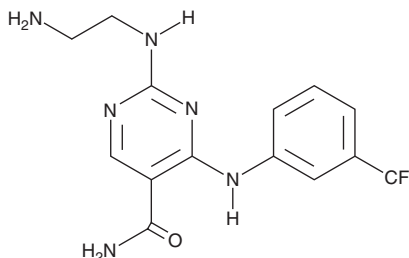
PRODUCT INFORMATION



Syk Inhibitor II

Item No. 18202

CAS Registry No.: 726695-51-8
Formal Name: 2-[(2-aminoethyl)amino]-4-[[3-(trifluoromethyl)phenyl]amino]-5-pyrimidinecarboxamide
Synonym: Spleen Tyrosine Kinase Inhibitor II
MF: C₁₄H₁₅F₃N₆O
FW: 340.3
Purity: ≥98%
UV/Vis.: λ_{max}: 260, 284 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

Spleen tyrosine kinase (Syk) inhibitor II is supplied as a crystalline solid. A stock solution may be made by dissolving the Syk inhibitor II in the solvent of choice, which should be purged with an inert gas. Syk inhibitor II is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of Syk inhibitor II in these solvents is approximately 15 and 25 mg/ml, respectively.

Description

Syk is a non-receptor tyrosine kinase that upon phosphorylation binds to immunoreceptor tyrosine-based activation motifs and mediates downstream signaling. Syk inhibitor II is a cell-permeable, pyrimidine-carboxamide compound that selectively and reversibly blocks Syk (IC₅₀ = 41 nM) in an ATP-competitive manner.¹ It is much less potent against PKCε, PKCβII, ZAP-70, Btk, and Itk (IC₅₀s = 5.1, 11, 11.2, 15.5, and 22.6 μM, respectively).¹ Syk inhibitor II has been shown to prevent FcεRI-mediated 5-HT release in RBL-2H3 cells *in vitro* (IC₅₀ = 460 nM) and to inhibit passive cutaneous anaphylaxis reactions in mice (ID₅₀ = 13.2 mg/kg, s.c.).¹

Reference

1. Hisamichi, J., Naito, R., Toyoshima, A., *et al.* Synthetic studies on novel Syk inhibitors. Part 1: Synthesis and structure-activity relationships of pyrimidine-5-carboxamide derivatives. *Bioorg. Med. Chem.* **13**(16), 4936-4951 (2005).

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, 02/05/2025

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