

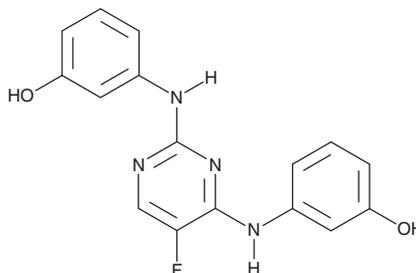
PRODUCT INFORMATION



R112

Item No. 22929

CAS Registry No.: 575474-82-7
Formal Name: 3,3'-[(5-fluoro-2,4-pyrimidinediyl)diimino]bis-phenol
MF: C₁₆H₁₃FN₄O₂
FW: 312.3
Purity: ≥98%
UV/Vis.: λ_{max}: 271, 321 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

R112 is supplied as a crystalline solid. A stock solution may be made by dissolving the R112 in the solvent of choice, which should be purged with an inert gas. R112 is soluble in the organic solvent DMSO at a concentration of approximately 30 mg/ml.

Description

R112 is an inhibitor of spleen tyrosine kinase (Syk; IC₅₀ = 226 nM in cultured human mast cells) that acts in an ATP-competitive and reversible manner (K_i = 96 nM).¹ It inhibits mast cell and basophil degranulation induced by anti-IgE cross-linking (EC₅₀s = 353 and 280 nM, respectively). It also inhibits basophil degranulation induced by an allergen with an EC₅₀ value of 490 nM. R112 inhibits the production of leukotriene C₄ (LTC₄; Item No. 20210) and decreases cytokine levels in cultured human mast cells.

References

1. Rossi, A.B., Herlaar, E., Braselmann, S., *et al.* Identification of the Syk kinase inhibitor R112 by a human mast cell screen. *J. Allergy Clin. Immunol.* **118**(3), 749-755 (2006).

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, 10/27/2022

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