

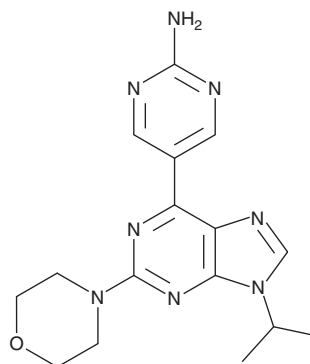
PRODUCT INFORMATION



VS-5584 analog

Item No. 30324

CAS Registry No.: 1246535-95-4
Formal Name: 5-[9-(1-methylethyl)-2-(4-morpholinyl)-9H-purin-6-yl]-2-pyrimidinamine
MF: C₁₆H₂₀N₈O
FW: 340.4
Purity: ≥98%
UV/Vis.: λ_{max}: 239, 289, 354 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

VS-5584 analog is supplied as a solid. A stock solution may be made by dissolving the VS-5584 analog in the solvent of choice, which should be purged with an inert gas. VS-5584 analog is soluble in the organic solvent chloroform at a concentration of approximately 0.5 mg/ml.

Description

VS-5584 analog is a dual inhibitor of PI3K α and mammalian target of rapamycin (mTOR; IC₅₀s = 17 and 150 nM, respectively).¹ It is a desmethyl analog of the dual PI3K/mTOR inhibitor VS-5584 (Item No. 16294).

Reference

1. Poulsen, A., Nagaraj, H., Lee, A., *et al.* Structure and ligand-based design of mTOR and PI3-kinase inhibitors leading to the clinical candidates VS-5584 (SB2343) and SB2602. *J. Chem. Inf. Model* **54**(11), 3238-3250 (2014).

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, 05/20/2020

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