

PRODUCT INFORMATION



GSK2250665A

Item No. 45942

CAS Registry No.: 1246030-96-5
Formal Name: *trans*-4-[[4-[(6-ethyl-2-benzothiazolyl)amino]-6-(phenylmethyl)-2-pyrimidinyl]amino]-cyclohexanol

MF: C₂₆H₂₉N₅OS

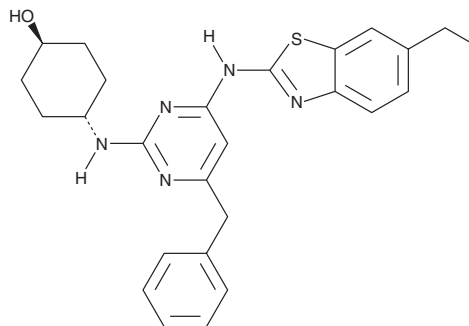
FW: 459.6

Purity: ≥98%

Supplied as: A 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

GSK2250665A is supplied as a solid. A stock solution may be made by dissolving the GSK2250665A in the solvent of choice, which should be purged with an inert gas. GSK2250665A is slightly soluble (0.1-1 mg/ml) in DMSO and in a 1:1 solution of acetonitrile:water.

Description

GSK2250665A is an IL-2-inducible T cell kinase (Itk) inhibitor (IC₅₀ = 0.63 nM).¹ It is selective for Itk over Aurora B kinase (IC₅₀ = 398 nM) and a panel of 11 kinases (IC₅₀s = 4-631 nM). GSK2250665A reduces the levels of IFN-γ in isolated human peripheral blood mononuclear cells (PBMCs; IC₅₀ = 25 nM).

Reference

1. Alder, C.M., Ambler, M., Campbell, A.J., *et al.* Identification of a novel and selective series of Itk inhibitors via a template-hopping strategy. *ACS Med. Chem. Lett.* **4**(10), 948-942 (2013).

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, 09/01/2026

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