

PRODUCT INFORMATION



GLPG2534

Item No. 42048

CAS Registry No.: 2095615-97-5

Formal Name: 6-[6-[2-(2-hydroxyethoxy)ethoxy]-5-[(tetrahydro-2H-pyran-4-yl)amino]-3H-imidazo[4,5-b]pyridin-3-yl]-3-pyridinecarbonitrile

MF: $C_{21}H_{24}N_6O_4$

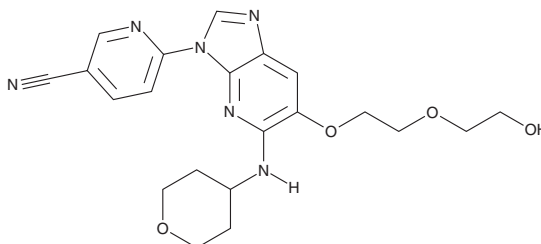
FW: 424.5

Purity: $\geq 98\%$

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

GLPG2534 is supplied as a solid. A stock solution may be made by dissolving the GLPG2534 in the solvent of choice, which should be purged with an inert gas. GLPG2534 is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in ethanol.

Description

GLPG2534 is an inhibitor of IL-1 receptor-associated kinase 4 (IRAK4; IC_{50} s = 6.4 and 3.5 nM for the human and mouse enzymes, respectively).¹ It is selective for IRAK4 over IRAK1 (IC_{50} = 179 nM) and a panel of 151 additional kinases at 1 μM but does inhibit FMS-related tyrosine kinase 3 (FLT3), tropomyosin-related kinase B (TrkB), and VEGFR3 (IC_{50} s = 2.2, 5.8, and 57 nM, respectively). GLPG2534 inhibits TNF- α release induced by CL097 (Item No. 39513) in isolated human peripheral blood mononuclear cells (PBMCs) and human whole blood (IC_{50} s = 122.8 and 568 nM, respectively). *In vivo*, GLPG2534 (30 mg/kg, p.o.) reduces ear thickness in mouse models of psoriasis induced by recombinant mouse IL-23 or imiquimod (Item No. 14956) and mouse models of atopic dermatitis induced by recombinant mouse IL-33 or MC903.

Reference

1. Lavazais, S., Jargosch, M., Dupont, S., *et al.* IRAK4 inhibition dampens pathogenic processes driving inflammatory skin diseases. *Sci. Transl. Med.* **15**(683), 683 (2023).

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, 01/21/2025

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