

PRODUCT INFORMATION

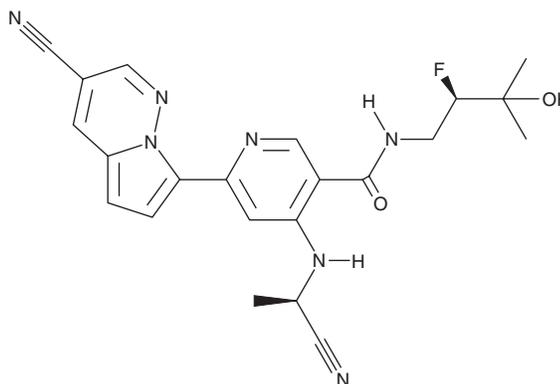


Edecesertib

Item No. 44286

CAS Registry No.: 2408839-73-4
Formal Name: 4-[[[(1R)-1-cyanoethyl]amino]-6-(3-cyanopyrrolo[1,2-b]pyridazin-7-yl)-N-[(2R)-2-fluoro-3-hydroxy-3-methylbutyl]-3-pyridinecarboxamide

Synonym: GS-5718
MF: C₂₂H₂₂FN₇O₂
FW: 435.5
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

Edecesertib is supplied as a solid. A stock solution may be made by dissolving the edecesertib in the solvent of choice, which should be purged with an inert gas. Edecesertib is soluble (≥10 mg/ml) in DMSO.

Description

Edecesertib is an inhibitor of IL-1 receptor-associated kinase 4 (IRAK4; IC₅₀ = 0.6 nM).¹ It is selective for IRAK4 over a panel of 484 additional kinases at 1,000 nM. Edecesertib inhibits toll-like receptor 7 (TLR7) agonist-induced IFN-α production in isolated human whole blood (IC₅₀ = 124 nM). Dietary administration of edecesertib (0.07% w/w in chow) reduces proteinuria and decreases plasma levels of TNF-α and IL-12 in an NZBWF1 mouse model of systemic lupus erythematosus (SLE).

Reference

1. Ammann, S.E., Cottell, J.J., Wright, N.E., *et al.* Discovery of edecesertib (GS-5718): A potent, selective inhibitor of IRAK4. *J. Med. Chem.* **68(11)**, 10619-10630 (2025).

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, 09/17/2025

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