

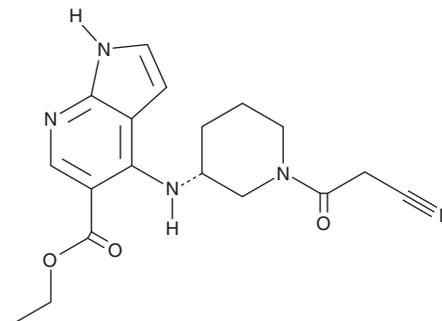
PRODUCT INFORMATION



Lepzacinib

Item No. 44278

CAS Registry No.: 2321488-47-3
Formal Name: 4-[[[(3R)-1-(2-cyanoacetyl)-3-piperidinyl] amino]-1H-pyrrolo[2,3-b]pyridine-5-carboxylic acid, ethyl ester
Synonym: ATI-1777
MF: C₁₈H₂₁N₅O₃
FW: 355.4
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

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

Description

Lepzacinib is an inhibitor of JAK1 and JAK3 (IC₅₀s = 1.5 and 3.6 nM, respectively).¹ It selectively inhibits IL-2-induced activation of JAK1/3 over IFN-γ-induced activation of JAK1/2, GM-CSF-induced activation of JAK2, and IL-12-induced activation of JAK2/tyrosine kinase 2 (TYK2) in isolated human peripheral blood mononuclear cells (PBMCs; IC₅₀s = 19, 97, 794, and 898 nM, respectively). Lepzacinib is also selective for JAK1 and JAK3 over a panel of 214 additional kinases at 1 μM. Topical application of lepzacinib (2% solution) inhibits IL-15-induced expression of mRNA encoding chemokine (C-C-motif) ligand 8 (CCL8) by 83% in a porcine model of atopic dermatitis.

Reference

1. Changelian, P., Xu, C., Mnich, S., *et al.* ATI-1777, a topical Jak1/3 inhibitor, may benefit atopic dermatitis without systemic drug exposure: Results from preclinical development and phase 2a randomized control study ATI-1777-AD-201. *JID Innov.* **4**(2), 100251 (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, 10/09/2025

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