

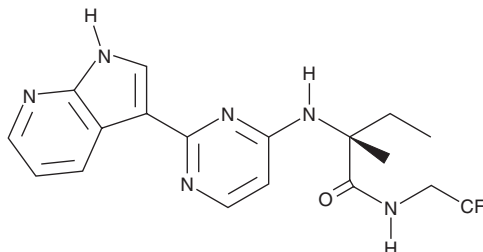
PRODUCT INFORMATION



VX-509

Item No. 19513

CAS Registry No.: 944842-54-0
Formal Name: (2R)-2-methyl-2-[[2-(1H-pyrrolo[2,3-b]pyridin-3-yl)-4-pyrimidinyl]amino]-N-(2,2,2-trifluoroethyl)-butanamide
Synonyms: Decernotinib, VRT-831509
MF: C₁₈H₁₉F₃N₆O
FW: 392.4
Purity: ≥98%
UV/Vis.: λ_{max}: 224, 262 nm
Supplied as: A crystalline 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

VX-509 is supplied as a crystalline solid. A stock solution may be made by dissolving the VX-509 in the solvent of choice. VX-509 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of VX-509 in these solvents is approximately 12, 20, and 25 mg/ml, respectively.

VX-509 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, VX-509 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. VX-509 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

VX-509 is a selective inhibitor of Janus kinase 3 (JAK3) with a K_i value of 2.5 nM in enzyme assays and an IC₅₀ value of 50-170 nM in cellular assays dependent on JAK3 activity with no measurable activity against other JAK isotypes.¹ At 10-100 mg/kg, it has been shown to dose dependently reduce proinflammatory responses in a rat collagen-induced arthritis model and a mouse model of oxazolone-induced delayed-type hypersensitivity.¹

Reference

1. Mahajan, S., Hogan, J.K., Shlyakhter, D., *et al.* VX-509 (decernotinib) is a potent and selective janus kinase 3 inhibitor that attenuates inflammation in animal models of autoimmune disease. *J. Pharmacol. Exp. Ther.* **353**(2), 405-414 (2016).

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, 06/09/2020

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