

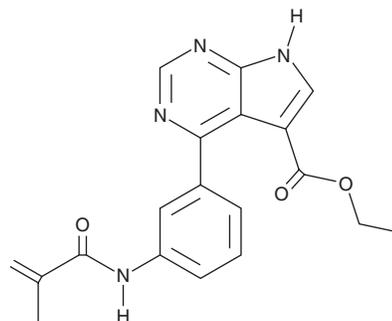
PRODUCT INFORMATION



JAK3-IN-2

Item No. 37344

CAS Registry No.: 1443235-95-7
Formal Name: 4-[3-[(2-methyl-1-oxo-2-propen-1-yl)amino]phenyl]-7H-pyrrolo[2,3-d]pyrimidine-5-carboxylic acid, ethyl ester
Synonyms: JAK3-Inhibitor 2, Janus-Associated Kinase 3 Inhibitor 2
MF: C₁₉H₁₈N₄O₃
FW: 350.4
Purity: ≥95%
UV/Vis.: λ_{max}: 250 nm
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

JAK3-IN-2 is supplied as a solid. A stock solution may be made by dissolving the JAK3-IN-2 in the solvent of choice, which should be purged with an inert gas. JAK3-IN-2 is soluble in methanol and DMSO.

Description

JAK3-IN-2 is an inhibitor of JAK3 ($K_i = 0.07$ nM).¹ It is selective for JAK3 over JAK1 and JAK2 (K_i s = 320 and 740 nM, respectively). JAK3-IN-2 inhibits IL-2-induced *STAT5* gene expression in a reporter assay ($IC_{50} = 70$ nM). *In vivo*, JAK3-IN-2 (100, 300, and 600 mg/kg) reduces paw swelling and ankle bone loss in a rat model of collagen-induced arthritis.

Reference

1. Elwood, F., Witter, D.J., Piesvaux, J., *et al.* Evaluation of JAK3 biology in autoimmune disease using a highly selective, irreversible JAK3 inhibitor. *J. Pharmacol. Exp. Ther.* **361**(2), 229-244 (2017).

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, 09/29/2023

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