

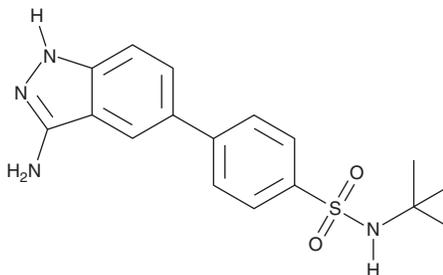
PRODUCT INFORMATION



JAK2 Inhibitor IV

Item No. 45230

CAS Registry No.: 1110502-30-1
Formal Name: 4-(3-amino-1H-indazol-5-yl)-N-(1,1-dimethylethyl)-benzenesulfonamide
Synonyms: Janus-Associated Kinase 2 Inhibitor IV
MF: C₁₇H₂₀N₄O₂S
FW: 344.4
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Special Conditions: Light sensitive



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

JAK2 inhibitor IV is supplied as a solid. A stock solution may be made by dissolving the JAK2 inhibitor IV in the solvent of choice, which should be purged with an inert gas. JAK2 inhibitor IV is slightly soluble (0.1-1 mg/ml) in acetonitrile and methanol.

Description

JAK2 inhibitor IV is an inhibitor of JAK2 (IC₅₀ = 78 nM).¹

Reference

1. Antonysamy, S., Hirst, G., Park, F., *et al.* Fragment-based discovery of JAK-2 inhibitors. *Bioorg. Med. Chem.* **19**(1), 279-282 (2009).

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, 02/12/2026

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