

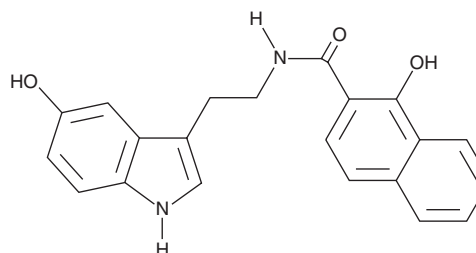
PRODUCT INFORMATION



MS-1020

Item No. 14273

CAS Registry No.: 1255516-86-9
Formal Name: 1-hydroxy-N-[2-(5-hydroxy-1H-indol-3-yl)ethyl]-2-naphthalenecarboxamide
MF: $C_{21}H_{18}N_2O_3$
FW: 346.4
Purity: $\geq 95\%$
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

MS-1020 is supplied as a solid. A stock solution may be made by dissolving the MS-1020 in the solvent of choice, which should be purged with an inert gas. MS-1020 is soluble in the organic solvent DMSO.

Description

MS-1020 is a cell-permeable inhibitor of janus kinase 3 (JAK3), strongly inhibiting constitutive autophosphorylation of JAK3 in L540 cells when used at 30-50 μM .¹ It is without effect on other JAK isoforms and several other kinases, including Src, Lyn, Akt, EGFR, and ERK1/2. Through its effects on JAK3, MS-1020 blocks phosphorylation of downstream signal transduction and activators of transcription isoforms, reducing the expression of anti-apoptotic genes, leading to cell death.¹

References

1. Kim, B.H., Oh, S.-R., Yin, C.-H., *et al.* MS-1020 is a novel small molecule that selectively inhibits JAK3 activity. *Br. J. Haematol.* **148**(1), 132-143 (2010).

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, 12/13/2022

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