

PRODUCT INFORMATION



Dasatinib N-oxide

Item No. 27845

CAS Registry No.: 910297-52-8

Formal Name: N-(2-chloro-6-methylphenyl)-
2-[[6-[4-(2-hydroxyethyl)-
4-oxido-1-piperazinyl]-2-
methyl-4-pyrimidinyl]amino]-5-
thiazolecarboxamide

MF: C₂₂H₂₆ClN₇O₃S

FW: 504.0

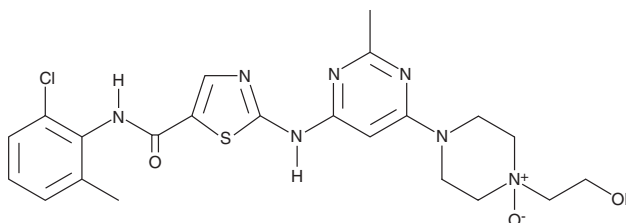
Purity: ≥95%

UV/Vis.: λ_{max}: 218, 230, 321 nm

Supplied as: A crystalline 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

Dasatinib N-oxide is supplied as a crystalline solid. A stock solution may be made by dissolving the dasatinib N-oxide in the solvent of choice, which should be purged with an inert gas. Dasatinib N-oxide is soluble in DMSO.

Description

Dasatinib N-oxide is a major metabolite of the Abl and Src kinase inhibitor dasatinib (Item No. 11498).¹ It is also a potential impurity in commercial preparations of dasatinib.²

References

1. Christopher, L.J., Cui, D., Li, W., *et al.* Biotransformation of [¹⁴C]dasatinib: In vitro studies in rat, monkey, and human and disposition after administration to rats and monkeys. *Drug Metab. Dispos.* **36**(7), 1341-1356 (2008).
2. Plachká, K., Švec, F., and Nováková, L. Ultra-high performance supercritical fluid chromatography in impurity control: Searching for generic screening approach. *Anal. Chim. Acta* **1039**, 149-161 (2018).

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, 11/18/2024

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