

# 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:** C<sub>22</sub>H<sub>26</sub>ClN<sub>7</sub>O<sub>3</sub>S

**FW:** 504.0

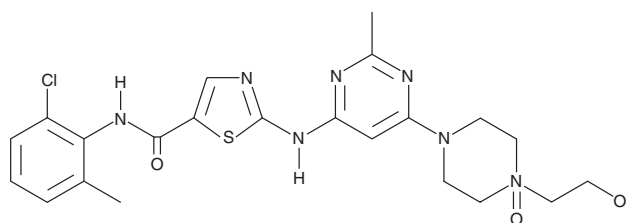
**Purity:** ≥95%

**UV/Vis.:** λ<sub>max</sub>: 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).<sup>1</sup> It is also a potential impurity in commercial preparations of dasatinib.<sup>2</sup>

### References

1. Christopher, L.J., Cui, D., Li, W., *et al.* Biotransformation of [<sup>14</sup>C]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, 10/28/2022

#### 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