

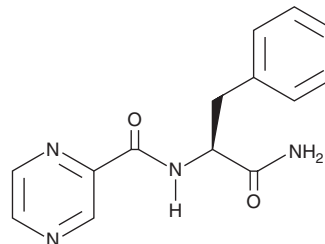
PRODUCT INFORMATION



Bortezomib Impurity A

Item No. 36499

CAS Registry No.: 289472-80-6
Formal Name: N-[(1S)-2-amino-2-oxo-1-(phenylmethyl)ethyl]-2-pyrazinecarboxamide
Synonym: Bortezomib Amide Impurity
MF: C₁₄H₁₄N₄O₂
FW: 270.3
Purity: ≥98%
UV/Vis.: λ_{max}: 269 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

Bortezomib impurity A is supplied as a solid. A stock solution may be made by dissolving the bortezomib impurity A in the solvent of choice, which should be purged with an inert gas. Bortezomib impurity A is slightly soluble in ethanol, methanol, chloroform, and DMSO.

Description

Bortezomib impurity A is a metabolite of the 20S proteasome inhibitor bortezomib (Item No. 10008822).¹ It is also a degradation product of bortezomib formed under oxidative conditions.²

References

1. Pekol, T., Daniels, J.S., Labutti, J., *et al.* Human metabolism of the proteasome inhibitor bortezomib: Identification of circulating metabolites. *Drug Metab. Dispos.* **33**(6), 771-777 (2005).
2. Martignac, M., Balayssac, S., Gilard, V., *et al.* Photochemical degradation of the anticancer drug bortezomib by V-UV/UV (185/254 nm) investigated by ¹H NMR fingerprinting: A way to follow aromaticity evolution. *J. Phys. Chem. A* **119**(24), 6215-6222 (2015).

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, 06/08/2023

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