

# PRODUCT INFORMATION



## Des(benzylpyridyl) Atazanavir

Item No. 27372

**CAS Registry No.:** 1192224-24-0  
**Formal Name:** (3S,8S,9S,12S)-3,12-bis(1,1-dimethylethyl)-8-hydroxy-4,11-dioxo-9-(phenylmethyl)-2,5,6,10,13-pentaazatetradecanedioic acid 1,14-dimethyl ester

**Synonym:** Atazanavir metabolite M1

**MF:** C<sub>26</sub>H<sub>43</sub>N<sub>5</sub>O<sub>7</sub>

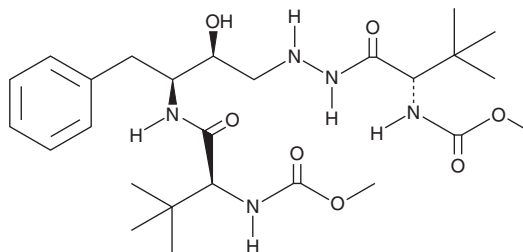
**FW:** 537.7

**Purity:** ≥98%

**Supplied as:** A crystalline solid

**Storage:** -20°C

**Stability:** ≥2 years



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

### Laboratory Procedures

Des(benzylpyridyl) atazanavir is supplied as a crystalline solid. A stock solution may be made by dissolving the des(benzylpyridyl) atazanavir in the solvent of choice, which should be purged with an inert gas. Des(benzylpyridyl) atazanavir is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of des(benzylpyridyl) atazanavir in these solvents is approximately 30 mg/ml.

### Description

Des(benzylpyridyl) atazanavir is a major metabolite of the HIV-1 protease inhibitor atazanavir (Item No. 11733).<sup>1,2</sup> It is formed from atazanavir by the cytochrome P450 (CYP) isoforms CYP3A4 and CYP2D6.<sup>2</sup>

### References

1. ter Heine, R., Hillebrand, M.J.X., Rosing, H., *et al.* Identification and profiling of circulating metabolites of atazanavir, a HIV protease inhibitor. *Drug Metab. Dispos.* **37(9)**, 1826-1840 (2009).
2. Li, F., Lu, J., Wang, L., *et al.* CYP3A-mediated generation of aldehyde and hydrazine in atazanavir metabolism. *Drug Metab. Dispos.* **39(3)**, 394-401 (2011).

#### 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, 11/13/2019

#### 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