

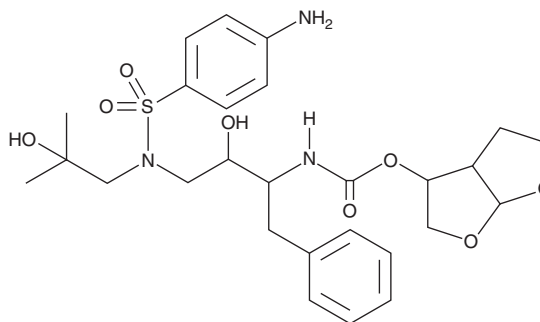
PRODUCT INFORMATION



Hydroxy Darunavir

Item No. 40225

CAS Registry No.: 1130635-75-4
Formal Name: N-[3-[[[(4-aminophenyl)sulfonyl](2-hydroxy-2-methylpropyl)amino]-2-hydroxy-1-(phenylmethyl)propyl]-carbamic acid, hexahydrofuro[2,3-b]furan-3-yl ester
Synonym: R-426857
MF: C₂₇H₃₇N₃O₈S
FW: 563.7
Purity: ≥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

Hydroxy darunavir is supplied as a solid. A stock solution may be made by dissolving the hydroxy darunavir in the solvent of choice, which should be purged with an inert gas. Hydroxy darunavir is soluble in organic solvents such as methanol and DMSO.

Description

Hydroxy darunavir is a metabolite of the HIV-1 protease inhibitor darunavir (Item Nos. 15866 | 31083).¹ It is formed from darunavir by isobutyl aliphatic hydroxylation.

Reference

1. Vermeir, M., Lachau-Durand, S., Mannens, G., *et al.* Absorption, metabolism, and excretion of darunavir, a new protease inhibitor, administered alone and with low-dose ritonavir in healthy subjects. *Drug Metab. Dispos.* **37**(4), 809-820 (2009).

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, 12/13/2023

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