

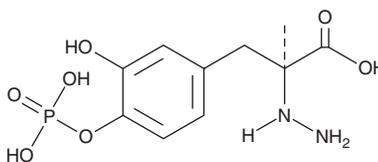
PRODUCT INFORMATION



Foscarbidopa

Item No. 43930

CAS Registry No.: 1907685-81-7
Formal Name: (αS)-α-hydrazinyl-3-hydroxy-α-methyl-4-(phosphonoxy)-benzenepropanoic acid
Synonym: ABBV-951
MF: C₁₀H₁₅N₂O₇P
FW: 306.2
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

Foscarbidopa is supplied as a solid. A stock solution may be made by dissolving the foscarbidopa in the solvent of choice, which should be purged with an inert gas. Foscarbidopa is slightly soluble (0.1-1 mg/ml) in methanol.

Description

Foscarbidopa is a monophosphate prodrug form of the dopamine decarboxylase inhibitor carbidopa (Item No. 23783).¹

Reference

1. Savaryn, J.P., Smith, R.L., Rosebraugh, M., *et al.* Metabolite profiling of foslevodopa/foscarbidopa in plasma of healthy human participants by LC-HRMS indicates no major differences compared to administration of levodopa/carbidopa intestinal gel. *Pharmacol. Res. Perspect.* **12(2)**, e1190 (2024).

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, 11/21/2025

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