

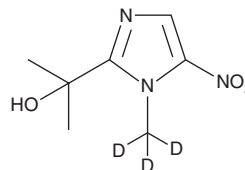
PRODUCT INFORMATION



Hydroxy Ipronidazole-d₃

Item No. 39166

CAS Registry No.: 1156508-86-9
Formal Name: α,α -dimethyl-1-(methyl-d₃)-5-nitro-1H-imidazole-2-methanol
MF: C₇H₈D₃N₃O₃
FW: 188.2
Chemical Purity: ≥98% (Hydroxy Ipronidazole)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₃); ≤1% d₀
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 ipronidazole-d₃ is intended for use as an internal standard for the quantification of hydroxy ipronidazole by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Hydroxy ipronidazole-d₃ is supplied as a solid. A stock solution may be made by dissolving the hydroxy ipronidazole-d₃ in the solvent of choice, which should be purged with an inert gas. Hydroxy ipronidazole-d₃ is slightly soluble in chloroform and methanol.

Description

Hydroxy ipronidazole is a metabolite of the nitroimidazole antiprotozoal agent ipronidazole (Item No. 35082).¹

Reference

1. Fraselle, S., Derop, V., Degroodt, J.-M., et al. Validation of a method for the detection and confirmation of nitroimidazoles and the corresponding hydroxy metabolites in pig plasma by high performance liquid chromatography-tandem mass spectrometry. *Anal. Chim. Acta* **586**(1-2), 383-393 (2007).

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, 07/10/2023

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