

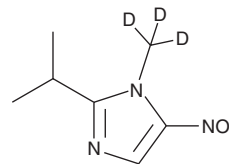
PRODUCT INFORMATION



Ipronidazole-d₃

Item No. 37194

CAS Registry No.: 1015855-83-0
Formal Name: 1-(methyl-d₃)-2-(1-methylethyl)-5-nitro-1H-imidazole
MF: C₇H₈D₃N₃O₂
FW: 172.2
Chemical Purity: ≥98% (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

Ipronidazole-d₃ is intended for use as an internal standard for the quantification of ipronidazole (Item No. 35082) 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).

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

Description

Ipronidazole is a nitroimidazole antiprotozoal agent.¹ Dietary administration of ipronidazole (0.00312-0.0094% w/w) reduces *H. meleagridis*-induced mortality in young turkeys. It also induces 100% disease clearance in bulls infected with *T. foetus* when administered at a dose of 60 g/animal.² Formulations containing ipronidazole have been used in the treatment of veterinary histomoniasis in turkeys and swine dysentery.

References

1. Bowen, T.E., Sullivan, T.W., and Grace, O.D. Effect of cupric sulfate on the prophylactic efficacy of 2-acetylamino-5-nitrothiazole, nifursol and ipronidazole against histomoniasis in turkeys. *Poult. Sci.* **50**(6), 1668-1672 (1971).
2. Skirrow, S.Z. and Bondurant, R.H. Treatment of bovine trichomoniasis with ipronidazole. *Aust. Vet. J.* **65**(5), 156 (1988).

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, 12/21/2022

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