

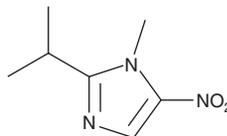
PRODUCT INFORMATION



Ipronidazole

Item No. 35082

CAS Registry No.: 14885-29-1
Formal Name: 1-methyl-2-(1-methylethyl)-5-nitro-1H-imidazole
Synonyms: NSC 109212, Ro 07-1554
MF: C₇H₁₁N₃O₂
FW: 169.2
Purity: ≥90%
UV/Vis.: λ_{max}: 310 nm
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 is supplied as a solid. A stock solution may be made by dissolving the ipronidazole in the solvent of choice, which should be purged with an inert gas. Ipronidazole is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of ipronidazole in these solvents is approximately 12, 11, and 14 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of ipronidazole can be prepared by directly dissolving the solid in aqueous buffers. The solubility of ipronidazole in PBS (pH 7.2) is approximately 0.16 mg/ml. We do not recommend storing the aqueous solution for more than one day.

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-5nitrothiazole, 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, 11/11/2022

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