

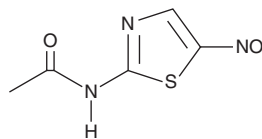
PRODUCT INFORMATION



Nithiamide

Item No. 36006

CAS Registry No.: 140-40-9
Formal Name: N-(5-nitro-2-thiazolyl)-acetamide
Synonyms: 2-Acetylamino-5-nitrothiazole, Aminitrozole, CL 5,279, NSC 31539, NSC 45914
MF: C₅H₅N₃O₃S
FW: 187.2
Purity: ≥95%
UV/Vis.: λ_{max}: 235, 339 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

Nithiamide is supplied as a solid. A stock solution may be made by dissolving the nithiamide in the solvent of choice, which should be purged with an inert gas. Nithiamide is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of nithiamide in these solvents is approximately 10 and 3 mg/ml, respectively.

Description

Nithiamide is an antiprotozoal agent.¹ It is active against *G. intestinalis* and *T. vaginalis* (IC₅₀s = 0.49 and 0.022 μM, respectively). Nithiamide (0.05% in the diet) reduces mortality in a turkey model of histomoniasis.²

References

1. Navarrete-Vázquez, G., Chávez-Silva, F., Colín-Lozano, B., *et al.* Synthesis of nitro(benzo)thiazole acetamides and in vitro antiprotozoal effect against amitochondriate parasites *Giardia intestinalis* and *Trichomonas vaginalis*. *Bioorg. Med. Chem.* **23(9)**, 2204-2210 (2015).
2. Flowers, A.I., Hall, C.F., and Grumbles, L.C. Chemotherapy of histomoniasis of turkeys I. The value of 1:2-dimethyl-5-nitroimidazole in prevention and treatment. *Avian Dis.* **9(3)**, 394-400 (1965).

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, 12/15/2022

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