

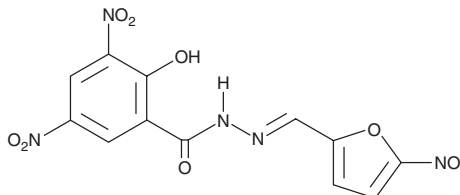
PRODUCT INFORMATION



Nifursol

Item No. 35139

CAS Registry No.: 16915-70-1
Formal Name: 2-hydroxy-3,5-dinitro-benzoic acid, 2-[(5-nitro-2-furanyl)methylene]hydrazide
Synonyms: Histomon, Salfuride
MF: C₁₂H₇N₅O₉
FW: 365.2
Purity: ≥98%
UV/Vis.: λ_{max}: 369 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

Nifursol is supplied as a solid. A stock solution may be made by dissolving the nifursol in the solvent of choice, which should be purged with an inert gas. Nifursol is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of nifursol in these solvents is approximately 5 mg/ml.

Description

Nifursol is an antiprotozoal.¹ Dietary administration of nifursol (37.5-100 ppm) reduces morbidity and mortality in a turkey model of histomoniasis. Formulations containing nifursol have previously been used in the prevention of histomoniasis in poultry.

Reference

1. Sullivan, T.W., Mitchell, R.J., and Grace, O.D. Prophylactic efficacy of nifursol against different levels of exposure to histomoniasis in turkeys 4 to 9 weeks of age. *Poult. Sci.* **51(6)**, 1956-1959 (1972).

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, 12/12/2022

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