

PRODUCT INFORMATION



Nicarbazine

Item No. 35110

CAS Registry No.: 330-95-0

Formal Name: N,N'-bis(4-nitrophenyl)-urea,
compd. with 4,6-dimethyl-2(1H)-
pyrimidinone

Synonyms: MK 75, NCZ, NSC 7171, Nicarbazine

MF: $C_{13}H_{10}N_4O_5 \cdot C_6H_8N_2O$

FW: 426.4

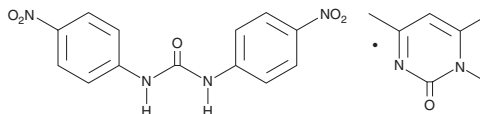
Purity: $\geq 95\%$

UV/Vis.: λ_{max} : 222, 341 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

Nicarbazine is supplied as a solid. A stock solution may be made by dissolving the nicarbazine in the solvent of choice, which should be purged with an inert gas. Nicarbazine is slightly soluble in DMSO.

Description

Nicarbazine is a coccidiostat.¹ It is composed of an equimolar complex of 4,4'-dinitrocarbanilide (DNC) and 2-hydroxy-4,6-dimethylpyrimidine (HDP). Nicarbazine (0.01-0.04% in the diet) prevents mortality of, and suppresses oocyst production in, chickens infected with the parasites *E. acervulina*, *E. necatrix*, *E. tenella*, and *P. gallinaceum*. Dietary administration of nicarbazine (100 ppm) decreases the production, weight, and pigmentation of eggs produced by broiler breeder chickens.² Formulations containing nicarbazine have been used in the prevention of coccidiosis in chickens.

References

1. Ott, W.H., Kuna, S., Porter, C.C., *et al.* Biological studies on nicarbazine, a new anticoccidial agent. *Poult. Sci.* **35**(6), 1355-1367 (1956).
2. Hughes, B.L., Jones, J.E., and Toler, J.E. Effects of exposing broiler breeders to nicarbazine contaminated feed. *Poult. Sci.* **70**(3), 476-482 (1991).

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