

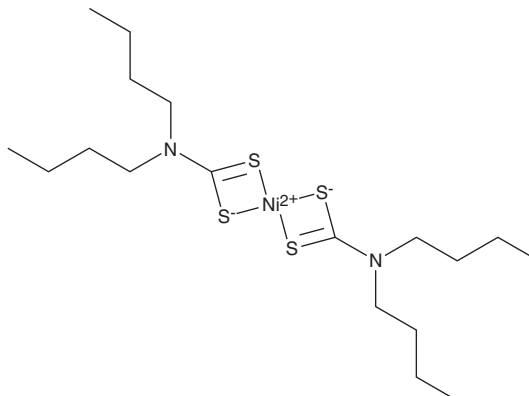
PRODUCT INFORMATION



NBC

Item No. 42583

CAS Registry No.: 13927-77-0
Formal Name: (SP-4-1)-bis(N,N-dibutylcarbamodithioato- κ S, κ S')-nickel
Synonyms: Nickel Dibutyldithiocarbamate, Nickel(II) bis(dibutyldithiocarbamate), NiDBC, NiDBDTC
MF: $C_{18}H_{36}N_2NiS_4$
FW: 467.4
Purity: $\geq 98\%$
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

NBC is supplied as a solid. A stock solution may be made by dissolving the NBC in the solvent of choice, which should be purged with an inert gas. NBC is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in ethanol.

Description

NBC is an antioxidant and antiozonant.¹ Topical application of NBC (35% in a powder) protects against ozone-induced bronzing of *P. vulgaris* leaves. It is not mutagenic to *S. typhimurium* in the Ames test.²

References

1. Dass, H.C. and Weaver, G.M. Modification of ozone damage to *Phaseolus vulgaris* by antioxidants, thiols and sulfhydryl reagents. *Can. J. Plant Sci.* **48**(6), 569-574 (1968).
2. Hedenstedt, A., Rannug, U., Ramel, C., et al. Mutagenicity and metabolism studies on 12 thiuram and dithiocarbamate compounds used as accelerators in the Swedish rubber industry. *Mutat. Res.* **68**(4), 313-325 (1979).

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, 02/19/2025

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