

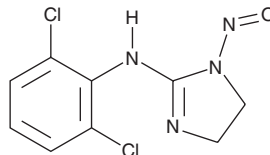
PRODUCT INFORMATION



N-Nitroso Clonidine

Item No. 34736

CAS Registry No.: 148950-49-6
Formal Name: N-(2,6-dichlorophenyl)-4,5-dihydro-1-nitroso-H-imidazol-2-amine
MF: $C_9H_8Cl_2N_4O$
FW: 259.1
Purity: $\geq 90\%$
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

N-Nitroso clonidine is supplied as a solid. A stock solution may be made by dissolving the N-nitroso clonidine in the solvent of choice, which should be purged with an inert gas. N-Nitroso clonidine is soluble in DMSO.

Description

N-Nitroso clonidine is a nitrosamine.¹ It is formed from clonidine (Item No. 15949) in acidic aqueous solutions.

Reference

1. Iley, J., Norberto, F., Rosa, E., *et al.* N-nitroso compounds. Part 2. The synthesis of N-nitrosoclonidine and its decomposition in aqueous acidic media. *J. Chem. Soc. Perk. T. 2* 591-594 (1993).

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, 12/08/2022

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