

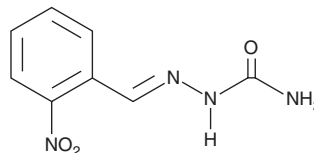
PRODUCT INFORMATION



2-Nitrobenzaldehyde semicarbazone

Item No. 42892

CAS Registry No.: 16004-43-6
Formal Name: 2-[(2-nitrophenyl)methylene]-hydrazinecarboxamide
Synonyms: o-Nitrobenzaldehyde semicarbazone, NPSCA, 2-NP-SCA
MF: C₈H₈N₄O₃
FW: 208.2
Purity: ≥95%
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

2-Nitrobenzaldehyde semicarbazone is supplied as a solid. A stock solution may be made by dissolving the 2-nitrobenzaldehyde semicarbazone in the solvent of choice, which should be purged with an inert gas. 2-Nitrobenzaldehyde semicarbazone is soluble (≥10 mg/ml) in DMSO.

Description

2-Nitrobenzaldehyde semicarbazone is an acid hydrolysis and thermal degradation product of nitrofurazone (Item No. 21656).¹ It also can be formed by thermal degradation of azodicarbonamide in azodicarbonamide-treated flour.²

References

1. Kulikovskii, A.V., Gorlov, I.F., Slozhenkina, M.I., *et al.* Determination of nitrofuran metabolites in muscular tissue by high-performance liquid chromatography with mass spectrometric detection. *J. Anal. Chem.* **74**, 906-912 (2019).
2. Becalski, A., Lau, B.P., Lewis, D., *et al.* Semicarbazide formation in azodicarbonamide-treated flour: A model study. *J. Agric. Food Chem.* **52**(18), 5730-5734 (2004).

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, 03/07/2025

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