

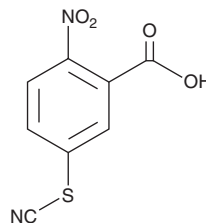
PRODUCT INFORMATION



2-Nitro-5-thiocyanatobenzoic Acid

Item No. 18236

CAS Registry No.: 30211-77-9
Formal Name: 2-nitro-5-thiocyanato-benzoic acid
Synonym: NTCB
MF: C₈H₄N₂O₄S
FW: 224.2
Purity: ≥98%
UV/Vis.: λ_{max}: 280 nm
Supplied as: A crystalline 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-Nitro-5-thiocyanatobenzoic acid is supplied as a crystalline solid. A stock solution may be made by dissolving the 2-nitro-5-thiocyanatobenzoic acid in the solvent of choice, which should be purged with an inert gas. 2-Nitro-5-thiocyanatobenzoic acid is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 2-nitro-5-thiocyanatobenzoic acid in ethanol and DMF is approximately 25 mg/ml and approximately 11 mg/ml in DMSO.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of 2-nitro-5-thiocyanatobenzoic acid can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of 2-nitro-5-thiocyanatobenzoic acid in PBS, pH 7.2, is approximately 0.3 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

2-Nitro-5-thiocyanatobenzoic acid is used to cyanilate proteins and to specifically cleave the amino-terminal peptide bond of cysteine residues.^{1,2}

References

1. Tang, H.-Y. and Speicher, D.W. Identification of alternative products and optimization of 2-nitro-5-thiocyanatobenzoic acid cyanilation and cleavage at cysteine residues. *Anal. Biochem.* **334**, 48-61 (2004).
2. Degani, Y. and Patchornik, A. Cyanilation of sulfhydryl groups by 2-nitro-5-thiocyanobenzoic acid. High-yield modification and cleavage of peptides at cysteine residues. *Biochem.* **13(1)**, 1-11 (1974).

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, 11/21/2022

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