

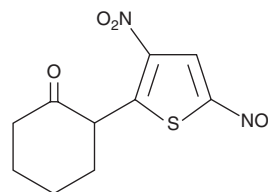
PRODUCT INFORMATION



NSC 245214

Item No. 9002543

CAS Registry No.: 31554-45-7
Formal Name: 2-(3,5-dinitro-2-thienyl)-cyclohexanone
MF: C₁₀H₁₀N₂O₅S
FW: 270.3
Purity: ≥95%
Supplied as: A crystalline solid
UV/Vis.: λ_{max}: 260, 302 nm
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

NSC 245214 is supplied as a crystalline solid. A stock solution may be made by dissolving the NSC 245214 in the solvent of choice, which should be purged with an inert gas. NSC 245214 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of NSC 245214 in these solvents is approximately 0.25, 10, and 15 mg/ml, respectively.

NSC 245214 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, NSC 245214 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. NSC 245214 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

NSC 245214 is a dinitrothiophene derivative.¹

Reference

1. Hellberg, L.H., Adams, C.C., Milligan, R.J., *et al.* Reaction of enamines of some bicyclic and monocyclic ketones with 2,4-dinitrochlorobenzene and 2-chloro-3,5-dinitrothiophene. *Org. Prep. Proced. Int.* **3(2)**, 77-81 (1971).

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, 06/29/2021

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