

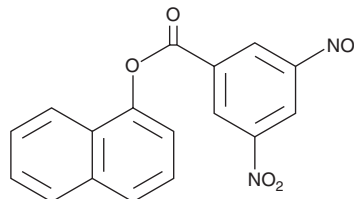
PRODUCT INFORMATION



1-Naphthyl 3,5-dinitrobenzoate

Item No. 11729

CAS Registry No.: 93261-39-3
Formal Name: 1-(3,5-dinitrobenzoate) 1-naphthalenol
Synonyms: 1-(3,5-Dinitrobenzoyloxy)naphthalene,
1-Naphthalenol 3,5-dinitrobenzoate
MF: C₁₇H₁₀N₂O₆
FW: 338.3
Purity: ≥98%
UV/Vis.: λ_{max}: 221 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

1-Naphthyl 3,5-dinitrobenzoate is supplied as a crystalline solid. A stock solution may be made by dissolving the 1-naphthyl 3,5-dinitrobenzoate in the solvent of choice, which should be purged with an inert gas. 1-Naphthyl 3,5-dinitrobenzoate is soluble in organic solvents such as methanol and DMSO. The solubility of 1-naphthyl 3,5-dinitrobenzoate in these solvents is approximately 0.1 and 0.5 mg/ml, respectively, with gentle warming required.

Description

1-Naphthyl 3,5-dinitrobenzoate has been identified as a dual 5-lipoxygenase and microsomal prostaglandin E synthase-1 inhibitor with IC₅₀ values of 1 and 3.6 μM, respectively.¹ The scaffold of this compound is being explored for anti-inflammatory drug design.¹

Reference

1. Wu, Y., He, C., Gao, Y., *et al.* Dynamic modeling of human 5-lipoxygenase-inhibitor interactions helps to discover novel inhibitors. *J. Med. Chem.* **55**, 2597-2605 (2012).

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, 12/13/2022

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