

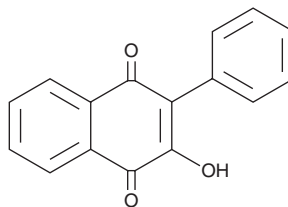
PRODUCT INFORMATION



NSC 102533

Item No. 42435

CAS Registry No.: 1150-59-0
Formal Name: 2-hydroxy-3-phenyl-1,4-naphthalenedione
Synonym: AN-04
MF: C₁₆H₁₀O₃
FW: 250.3
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

NSC 102533 is supplied as a solid. A stock solution may be made by dissolving the NSC 102533 in the solvent of choice, which should be purged with an inert gas. NSC 102533 is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in ethanol.

Description

NSC 102533 is an antagonist of the purinergic P2X₇ receptor.¹ It inhibits ATP-induced dye uptake in HEK293 cells expressing human purinergic P2X₇ receptors (IC₅₀ = 22.4 nM). NSC 102533 also inhibits common lancehead (*B. atrox*) venom proteolysis in a concentration-dependent manner.² It inhibits ATP-induced IL-1β secretion in human THP-1 monocytes and primary mouse peritoneal macrophages (IC₅₀s = 9.8 and 11.2 nM, respectively).¹ Intradermal administration of NSC 102533 (3 mg/kg) inhibits common lancehead (*B. atrox*) and jararaca (*B. jararaca*) venom-induced skin hemorrhage in mice.² It inhibits carrageenan-induced paw edema in mice when administered at doses ranging from 0.01 to 1 mg/kg.¹

References

1. Faria, R.X., Oliveira, F.H., Salles, J.P., *et al.* 1,4-Naphthoquinones potently inhibiting P2X₇ receptor activity. *Eur. J. Med. Chem.* **143**, 1361-1372 (2018).
2. Strauch, M.A., Tomaz, M.A., Monteiro-Machado, M., *et al.* Lapachol and synthetic derivatives: *In vitro* and *in vivo* activities against *Bothrops* snake venoms. *PLoS One* 14(1), (2019).

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, 12/12/2024

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