

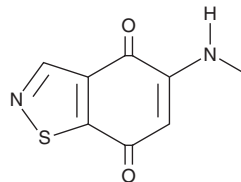
PRODUCT INFORMATION



Pronqodine A

Item No. 43836

CAS Registry No.: 1424195-39-0
Formal Name: 5-(methylamino)-1,2-benzisothiazole-4,7-dione
MF: C₈H₆N₂O₂S
FW: 194.2
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Bacterium/*Streptomyces* sp. MK832-95F2



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

Laboratory Procedures

Pronqodine A is supplied as a solid. A stock solution may be made by dissolving the pronqodine A in the solvent of choice, which should be purged with an inert gas. Pronqodine A is soluble in methanol and DMSO. Pronqodine A is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

Pronqodine A is a quinone that has been found in *Streptomyces* and has enzyme inhibitory and anti-inflammatory activities.^{1,2} It inhibits NAD(P)H:quinone acceptor oxidoreductase 1 (NQO1; IC₅₀ = 131.5 nM).² Pronqodine A inhibits bradykinin-induced prostaglandin E₂ (PGE₂; Item No. 14010), 6-keto PGF_{1α} (Item No. 15210), and PGD₂ (Item No. 12010) production in SW 982 human synovial sarcoma cells (IC₅₀s = 9, 19, and 7 nM, respectively).¹ It also induces the production of reactive oxygen species (ROS) in SW 982 cells.

References

1. Nakae, K., Adachi, H., Sawa, R., *et al.* NAD(P)H quinone oxidoreductase 1 (NQO1)-bioactivated pronqodine A regulates prostaglandin release from human synovial sarcoma cells. *J. Nat. Prod.* **76(4)**, 510-515 (2013).
2. Blunt, C.E., Torcuk, C., Liu, Y., *et al.* Synthesis and intracellular redox cycling of natural quinones and their analogues and identification of indoleamine-2,3-dioxygenase (IDO) as potential target for anticancer activity. *Angew. Chem. Int. Ed. Engl.* **54(30)**, 8740-8745 (2015).

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, 07/09/2025

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