

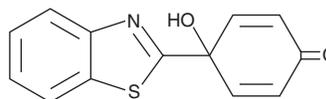
PRODUCT INFORMATION



PMX-464

Item No. 15343

CAS Registry No.: 485842-97-5
Formal Name: 4-(2-benzothiazolyl)-4-hydroxy-2,5-cyclohexadien-1-one
Synonym: AW-464
MF: C₁₃H₉NO₂S
FW: 243.3
Purity: ≥98%
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

PMX-464 is supplied as a solid. A stock solution may be made by dissolving the PMX-464 in the solvent of choice, which should be purged with an inert gas. PMX-464 is slightly soluble (0.1-1 mg/ml) in acetonitrile and DMSO.

Description

PMX-464 is a quinol inhibitor of thioredoxin reductase (TrxR; mean IC₅₀ = 6.5 μM for the rat enzyme).¹ It inhibits cell growth in the NCI-60 panel of cancer cell lines (mean GI₅₀ = 0.23 μM). PMX-464 (30 μM) inhibits IL-1β-induced NF-κB DNA binding in A549 cells.²

References

1. Chew, E.-H., Lu, J., Bradshaw, T.D., *et al.* Thioredoxin reductase inhibition by antitumor quinols: A quinol pharmacophore effect correlating to antiproliferative activity. *FASEB J.* **22(6)**, 2072-2083 (2008).
2. Callister, M.E., Pinhu, L., Catley, M.C., *et al.* PMX464, a thiol-reactive quinol and putative thioredoxin inhibitor, inhibits NF-kappaB-dependent proinflammatory activation of alveolar epithelial cells. *Br. J. Pharmacol.* **155(5)**, 661-672 (2008).

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, 06/09/2025

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