

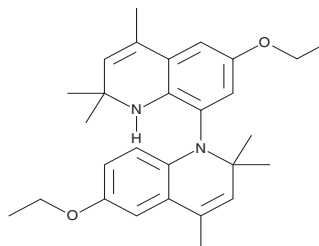
PRODUCT INFORMATION



Ethoxyquin Dimer

Item No. 29823

CAS Registry No.: 74681-77-9
Formal Name: 4,4'-hexamethyl-2'-dihydro-2,2,2',2'6,6'-diethoxy-1'1(2H),8'-biquinoline
Synonym: EQDM
MF: C₂₈H₃₆N₂O₂
FW: 432.6
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

Ethoxyquin dimer is supplied as a solid. A stock solution may be made by dissolving the ethoxyquin dimer in the solvent of choice, which should be purged with an inert gas. Ethoxyquin dimer is slightly soluble in chloroform and ethyl acetate.

Description

Ethoxyquin dimer is an antioxidant and metabolite of ethoxyquin (Item No. 23094).¹ It prevents oxidation of polyunsaturated fatty acids in fish meal and fish oil. Dietary administration of ethoxyquin dimer (0.1, 0.3, and 0.5% w/w) induces microvesicular steatosis and hepatocyte necrosis, as well as increases liver levels of oxidized glutathione and total lipids in mice.²

References

1. Koning, A.J. A new method for measuring efficacies of antioxidants in fish meal. *Int. J. Food Prop.* **1(3)**, 255-261 (2009).
2. Bernhard, A., Rasinger, J.D., Wisløff, H., *et al.* Subchronic dietary exposure to ethoxyquin dimer induces microvesicular steatosis in male BALB/c mice. *Food Chem. Toxicol.* **118**, 608-625 (2018).

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, 12/09/2022

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