

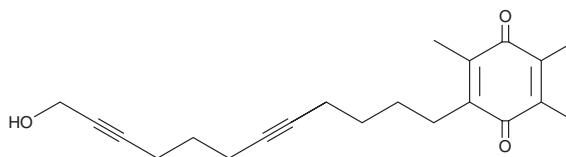
PRODUCT INFORMATION



AA-861

Item No. 13263

CAS Registry No.: 80809-81-0
Formal Name: 2-(12-hydroxy-5,10-dodecadiyn-1-yl)-3,5,6-trimethyl-2,5-cyclohexadiene-1,4-dione
Synonyms: A-61589, Docebenone
MF: C₂₁H₂₆O₃
FW: 326.4
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

AA-861 is supplied as a solid. A stock solution may be made by dissolving the AA-861 in the solvent of choice, which should be purged with an inert gas. AA-861 is soluble in the organic solvent DMSO (sonicated) at a concentration of approximately 250 mg/ml.

Description

AA-861 is an inhibitor of 5-lipoxygenase (5-LO; IC₅₀ = 0.8 μM).¹ It is greater than 100-fold selective for 5-LO over 12-LO and cyclooxygenase (COX). AA-861 inhibits the production of 5-HETE and leukotriene B₄ (LTB₄; Item Nos. 20110 | 10007240) induced by A23187 (Item No. 22030) with IC₅₀ values of 0.03 and 0.08 μM, respectively, in isolated rat peritoneal macrophages.² It reduces paw edema, pleurisy, and cell infiltration induced by carrageenan in rats when administered at a dose of 80 mg/kg. AA-861 (20 mg/kg) reduces ovalbumin-induced bronchoconstriction in an ovalbumin-sensitized guinea pig model of experimental allergic asthma.

References

1. Yoshimoto, T., Yokoyama, C., Ochi, K., *et al.* 2,3,5-Trimethyl-6-(12-hydroxy-5,10-dodecadiynyl)-1,4-benzoquinone (AA861), a selective inhibitor of the 5-lipoxygenase reaction and the biosynthesis of slow-reacting substance of anaphylaxis. *Biochim. Biophys. Acta* **713**(2), 470-473 (1982).
2. Ashida, Y., Saijo, T., Kuriki, H., *et al.* Pharmacological profile of AA-861, a 5-lipoxygenase inhibitor. *Prostaglandins* **26**(6), 955-972 (1983).

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, 07/14/2023

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