

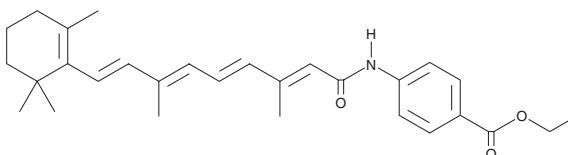
PRODUCT INFORMATION



Viaminate

Item No. 40118

CAS Registry No.: 53839-71-7
Formal Name: N-[4-(ethoxycarbonyl)phenyl]-retinamide
MF: C₂₉H₃₇NO₃
FW: 447.6
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

Viaminate is supplied as a solid. A stock solution may be made by dissolving the viaminate in the solvent of choice, which should be purged with an inert gas. Viaminate is soluble in DMSO. It is slightly soluble in chloroform and methanol

Description

Viaminate is a derivative of the retinoid vitamin A (Item No. 20241).^{1,2} It reverses keratinization of vitamin A-deficient primary hamster tracheal epithelial cells when used at a concentration of 10 nM.¹ Viaminate (1.7-15 mg/kg, p.o.) reduces epidermal keratinization, oil accumulation, ear thickness and edema, and ear skin levels of Il-6, Il-1 β , and Tnf- α in a rat model of acne induced by *P. acnes*.²

References

1. Jaegar, E.P., Jurs, P.C., and Stouch, T.R. Structure-activity relationship studies of retinoid cancer inhibition. *Eur. J. Med. Chem.* **28(4)**, 275-290 (1993).
2. Cao, J., Xu, M., Zhu, L., et al. Viaminate ameliorates *Propionibacterium acnes*-induced acne via inhibition of the TLR2/NF- κ B and MAPK pathways in rats. *Naunyn Schmiedebergs Arch Pharmacol.* **396(7)**, 1487-1500 (2023).

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, 03/05/2024

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