

PRODUCT INFORMATION



PDDHV

Item No. 20445

CAS Registry No.: 179469-40-0

Formal Name: 4-hydroxy-3-methoxy-benzeneacetic acid, [(1aR,1bS,4aR,7aS,7bS,8R,9R,9aS)-1a,1b,4,4a,5,7a,7b,8,9,9a-decahydro-4a,7b-dihydroxy-1,1,6,8-tetramethyl-5-oxo-9,9a-bis[(1-oxodecyl)oxy]-1H-cyclopropa[3,4]benz[1,2-e]azulen-3-yl)methyl ester

Synonym: Phorbol 12,13-didecanoate 20-homovanillate

MF: $C_{49}H_{72}O_{11}$

FW: 837.1

Purity: $\geq 95\%$

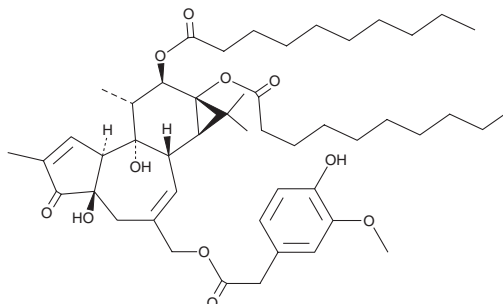
Supplied as: A solid

Storage: -20°C

Stability: ≥ 4 years

Special Conditions: Protect from light

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



Laboratory Procedures

PDDHV is supplied as a solid. A stock solution may be made by dissolving the PDDHV in the solvent of choice, which should be purged with an inert gas. PDDHV is soluble in the organic solvent methanol.

Description

PDDHV is a resiniferatoxin-type phorboid vanilloid with capsaicin-like selectivity for the cloned rat transient receptor potential cation channel subfamily V member 1 (TRPV1), formerly known as vanilloid receptor 1.¹ It induces Ca^{2+} -uptake by rat dorsal root ganglion neurons with an EC_{50} value of 70 nM.¹

Reference

1. Szallasi, A., Szabó, T., Bíró, T., *et al.* Resiniferatoxin-type phorboid vanilloids display capsaicin-like selectivity at native vanilloid receptors on rat DRG neurons and at the cloned vanilloid receptor VR1. *Br. J. Pharmacol.* **128**(2), 428-434 (1999).

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, 12/19/2022

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