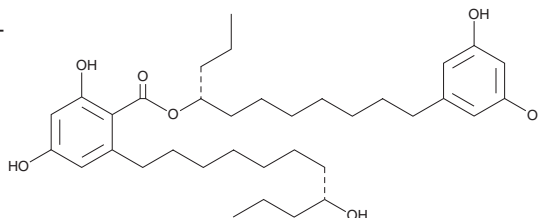


PRODUCT INFORMATION



Integracin B Item No. 28088

CAS Registry No.: 224186-05-4
Formal Name: 2,4-dihydroxy-6-[(8R)-8-hydroxyundecyl]-benzoic acid, (1R)-8-(3,5-dihydroxyphenyl)-1-propyloctyl ester
MF: $C_{35}H_{54}O_7$
FW: 586.8
Purity: $\geq 95\%$
Supplied as: An oil
Storage: -20°C
Stability: ≥ 4 years
Item Origin: Fungus/*Cytospora* sp.



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

Laboratory Procedures

Integracin B is supplied as an oil. A stock solution may be made by dissolving the integracin B in the solvent of choice, which should be purged with an inert gas. Integracin B is soluble in methanol and DMSO.

Description

Integracin B is a fungal metabolite and an inhibitor of HIV-1 integrase ($IC_{50} = 6.1 \mu\text{M}$ in an HIV-1 integrase coupled assay).¹ It also inhibits HIV-1 integrase in a strand transfer assay ($IC_{50} = 17 \mu\text{M}$).

References

1. Singh, S.B., Zink, D.L., Bills, G.F., et al. Discovery, structure and HIV-1 integrase inhibitory activities of integracins, novel dimeric alkyl aromatics from *Cytospora* sp. *Tetrahedron Lett.* **43**(9), 1617-1620 (2002)

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, 08/11/2023

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