

12-

Methoxydodecanoic acid

Myristate analog

A myristate analog which has similar chain length but reduced hydrophobicity compared to myristate. Inhibits replication of HIV-1. Its specific mechanism of action is not clear but it is believed to either inhibit the activity of N-myristoyltransferase or modulate the function of target acylated proteins.

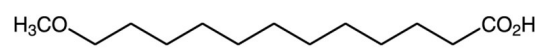
Ordering Information

[Order Online »](#)

BML-G234-0010	10mg
BML-G234-0050	50mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at room temperature for up to 1 year. Store solutions at -20°C for up to 6 months.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance	White to off-white solid.
CAS	92169-28-3
Formula	$C_{13}H_{26}O_3$
MW	230.3
MeltingPoint	45-46°C
Purity	≥98% (TLC)
Solubility	Soluble in ethanol (25 mg/mL) and DMSO (25 mg/mL).

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com