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.

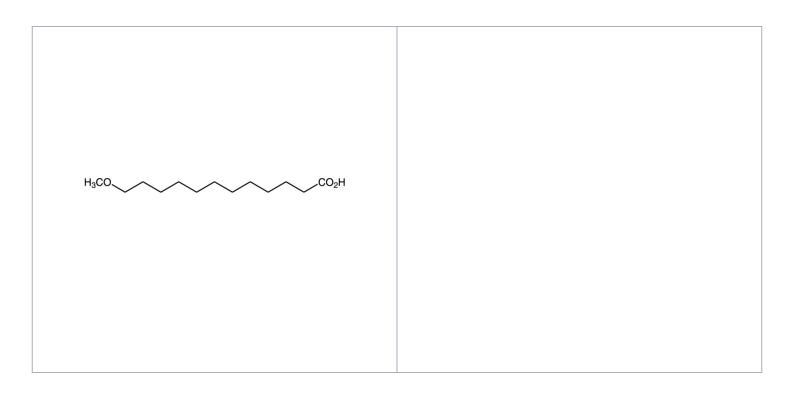
Ordering Information

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

