

C2 Dihydroceramide

Negative control for C2 ceramide

C₂ Dihydroceramide is inactive and may be used as a negative control for C₂ ceramide.

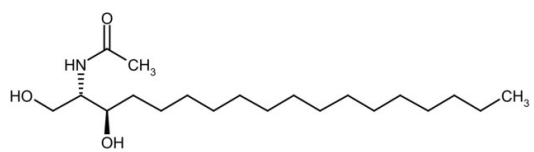
Ordering Information

[Order Online »](#)

BML-SL101-0005	5mg
----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at -20°C for up to 1 year. Store solutions at -20°C for up to 3 months.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	White solid.
CAS	13031-64-6
Formula	$C_{20}H_{41}NO_3$
MW	343.5
Purity	≥98% (TLC)
Solubility	Soluble in ethanol (5mg/ml) or DMSO (5mg/ml: dissolve in hot DMSO, then cool to RT).

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