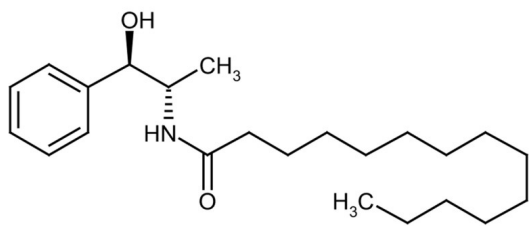


# L-erythro MAPP

Negative control for D-erythro-MAPP

Useful as a negative control for D-erythro-MAPP. It does not raise endogenous ceramide levels or inhibit the growth of HL-60 cells.

**Manuals, SDS & CofA** [View Online »](#)



## Handling & Storage

<b>Use/Stability</b>	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.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	(1S,2R-L-erythro-2-N-Myristoylamino)-1-phenyl-1-propanol
<b>Appearance</b>	White to off-white solid.
<b>CAS</b>	143492-39-1
<b>Formula</b>	$C_{23}H_{39}NO_2$
<b>MW</b>	361.6
<b>Purity</b>	≥98% (TLC)
<b>Solubility</b>	Soluble in DMSO (25 mg/ml) or ethanol (25 mg/ml).

Last modified: May 29, 2024



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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