

# N-Oleylethanolamide

TRPV1 activator

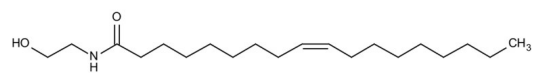
Activates TRPV1. Endogenous anandamide-like lipid which regulates appetite but does not bind to cannabinoid receptors. In rats, food intake is inhibited by oral oleylethanolamide. Induces satiety by activation of PPAR $\alpha$ . Inhibits ceramidase.

Citations: 10

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

Appearance	White solid.
CAS	111-58-0
Couple Target	Ceramidase, PPAR, TRP channel
Couple Type	Activator, Inhibitor, Ligand
Formula	$C_{20}H_{39}NO_2$
MW	325.5
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (>25mg/ml), methanol or 100% ethanol (>25mg/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-be@enzolifesciences.com](mailto:info-be@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)