

# Palmitylethanolamide

## Endocannabinoid

Endogenous cannabinoid. Weak ligand of CB<sub>1</sub> (K<sub>i</sub>=23.8μM) and CB<sub>2</sub> (K<sub>i</sub>=13.9μM) receptor. Inhibits fatty acid amide hydrolase (FAAH) (IC<sub>50</sub>=5.1μM). Immunosuppressant, anti-inflammatory, anti-nociceptive and anti-convulsant *in vivo*. The exact mode of action has not yet been revealed. It has been suggested that PEA: i) binds to a yet to be discovered cannabinoid receptor similar to CB<sub>2</sub>; ii) administered *in vivo* elicits the synthesis of endogenous agonists of CB<sub>2</sub>; iii) acts as an "entourage" compound by enhancing the activity and/or by influencing the turnover of endogenous agonists of CB<sub>2</sub>, possibly but not uniquely, by inhibiting their degradation.

Citations: 10

[View Online »](#)

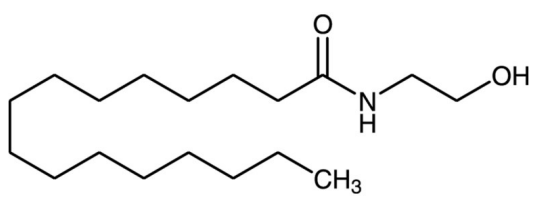
## Ordering Information

[Order Online »](#)

BML-FA018-0050	50mg
----------------	------

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored, as supplied, at 0-4°C. Stock solutions are stable for up to 3 months at -20°C.
Long Term Storage	+4°C
Shipping	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	PEA, Palmidrol
Appearance	White solid.
CAS	544-31-0
Couple Target	Cannabinoid receptor, FAAH
Couple Type	Inhibitor, Ligand
Formula	$C_{18}H_{37}NO_2$
MW	299.5
Purity	≥98% (TLC)
Solubility	Soluble in DMSO or 100% ethanol.



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)