

# N-Arachidonyl-GABA

Ultra-pure

One member of the class of endocannabinoids related to anandamide (arachidonoyl aminoacids) isolated from bovine brain. Displays analgesic activity.

Citations: 2

[View Online »](#)

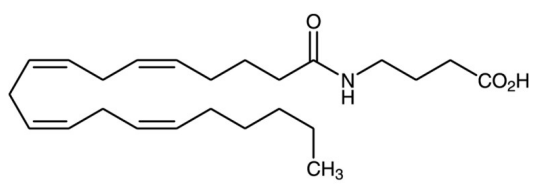
## Ordering Information

[Order Online »](#)

BML-FA035-0005	5mg
BML-FA035-0025	25mg

Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

<b>Use/Stability</b>	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored, as supplied, at -80°C. Stock solutions are stable for up to 3 months at -80°C when stored under inert atmosphere.
<b>Handling</b>	Protect from air. Keep under inert gas.
<b>Long Term Storage</b>	-80°C
<b>Shipping</b>	Dry Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	NA-GABA, 4-(1-oxo-5Z,8Z,11Z,14Z-Eicosatetraenyl)amino-butanoic acid, N-Arachidonoyl-γ-aminobutyric acid
<b>Appearance</b>	Pale yellow viscous oil.
<b>CAS</b>	128201-89-8
<b>Formula</b>	$C_{24}H_{39}NO_3$
<b>MW</b>	389.6
<b>Purity</b>	≥98% (TLC)
<b>Solubility</b>	Soluble in DMSO (25mg/ml) or ethanol (25mg/ml).



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)