

N-Arachidonoyl-serotonin

TRPV1 antagonist

Unnatural fatty acid-neurotransmitter conjugate. Potent TRPV1 antagonist (IC_{50} =40nM). Inhibits fatty acid amide hydrolase (FAAH) (IC_{50} =5.6mM). Enhances levels of anandamide and 2-AG *in vivo*. Displays analgesic activity in both acute and chronic peripheral pain models but shows little efficacy in an inflammatory bowel disease model.

Citations: 4

[View Online »](#)

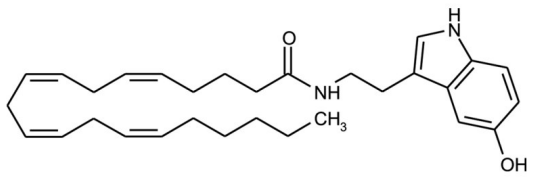
Ordering Information

[Order Online »](#)

BML-VR106-0010	10mg
----------------	------

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.
Handling	Protect from air.
Long Term Storage	-20°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Pale yellow oil.
CAS	187947-37-1
Couple Target	FAAH, TRP channel
Couple Type	Blocker, Inhibitor
Formula	$C_{30}H_{42}N_2O_2$
MW	462.7
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (>25mg/ml), 100% ethanol (15mg/ml) or dimethyl formamide (30mg/ml); also soluble in PBS, pH 7.2 (290µg/ml).

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-fr@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