

AM-404

Anandamide transport inhibitor

Analog of anandamide (Prod. No. BML-FA017). Potentiates the activity of endogenous anandamide by blocking its re-uptake into presynaptic membranes. Activates TRPV1 ($EC_{50}=0.04\mu\text{M}$) at concentrations lower than those required to inhibit anandamide transport into the cell (neuronal (C6 glioma cells): $IC_{50}=10\mu\text{M}$; non-neuronal (rat RBL-2H3 cells): $IC_{50}=11\mu\text{M}$). Low affinity to FAAH ($IC_{50}=5.9\mu\text{M}$), to CB_1 receptor ($IC_{50}=1.76\mu\text{M}$) and to CB_2 receptor ($IC_{50}>1\mu\text{M}$).

Citations: 12

[View Online »](#)

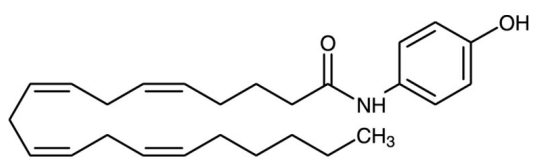
Ordering Information

[Order Online »](#)

BML-CR106-0050	50mg
----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for up to one year after receipt when stored at -20°C. Solutions are stable for up to 3 months when stored at -20°C under an inert atmosphere of nitrogen or argon.
Handling	Protect from air. Keep under inert gas.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	N-(4-Hydroxyphenyl)-5Z,8Z,11Z,14Z-eicosatetraenamide, N-(4-Hydroxyphenyl)arachidonoylamide
Appearance	Waxy solid.
CAS	183718-77-6
Couple Target	Anandamide transporter, TRP channel
Couple Type	Activator, Inhibitor
Formula	$C_{26}H_{37}NO_2$
MW	395.6
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (>25mg/ml) or 100% ethanol (>25mg/ml).



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