

Arachidonoyl-AMC

FAAH substrate

Fluorogenic substrate for fatty acid amide hydrolase (FAAH). This substrate is the basis for a simple, selective and sensitive assay for FAAH which is amenable to high-throughput screening.

Citations: 1

[View Online »](#)

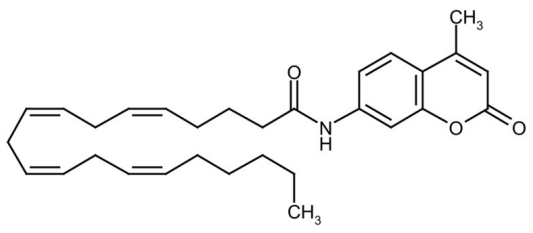
Ordering Information

[Order Online »](#)

BML-ST504-0005	5mg
BML-ST504-0025	25mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Off-white solid.
Couple Target	FAAH
Couple Type	Substrate
Formula	$C_{30}H_{39}O_3$
MW	461.6
Purity	≥99%
Solubility	Soluble in DMSO (>25mg/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-be@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