

7-HC-arachidonate

Phospholipase substrate

PLA₂ fluorogenic substrate for use in assays with phospholipid and Ca²⁺.
For use with purified enzyme, as this substrate is nonspecific. Ideal for drug discovery, HTS, and similar applications. Ex.: 360 nm, Em.: 460 nm

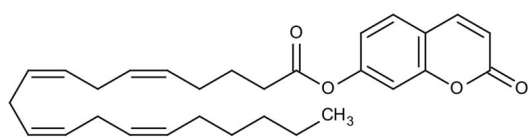
Ordering Information

[Order Online »](#)

BML-ST502-0010	10mg
BML-ST502-0050	50mg

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. Hexane should be evaporated by nitrogen or argon.
Handling	Protect from oxygen. Protect from air.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Colorless oil.
CAS	161180-11-6
Couple Target	Phospholipase, PLA
Couple Type	Substrate
Formula	$C_{29}H_{36}O_4$
MW	448.6
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (50mg/ml) or hexane. Use only de-gassed exane. Do not use ethanol as it may degrade product.



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)