

HELSS

Phospholipase inhibitor

HELSS is a potent, irreversible, mechanism-based inhibitor of Ca^{2+} -independent PLA_2 (70% inhibition at 100 nM). It possesses a 1000-fold selectivity for Ca^{2+} -independent PLA_2 vs. Ca^{2+} -dependent PLA_2 . It has been used to identify the PLA_2 involved in the release of arachidonate induced by vasopressin stimulation of A-10 smooth muscle cells ($\text{IC}_{50}=1 \mu\text{M}$). It has also been shown to inhibit the Mg^{2+} -dependent phosphatidate phosphohydrolase from P388D1 macrophages ($\text{IC}_{50}\sim 8 \mu\text{M}$).

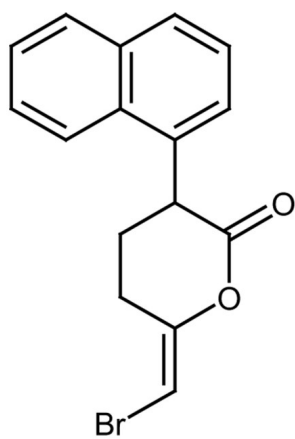
Ordering Information

[Order Online »](#)

BML-ST340-0005	5mg
BML-ST340-0025	25mg

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 under at atmosphere of argon or nitrogen at -20°C for up to 1 month.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Haloenol lactone suicide substrate
Appearance	White solid.
CAS	88070-98-8
Couple Target	Phospholipase, PLA
Couple Type	Inhibitor
Formula	$C_{16}H_{13}BrO_2$
MW	317.2
MeltingPoint	104-106°C
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (25mg/ml).

Last modified: May 29, 2024



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-
fr@enzolifesciences.com](mailto:info-fr@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)