

Epoxide hydrolase inhibitor

Novel inhibitor of soluble epoxide hydrolase (IC_{50} =18nM and 69nM for the mouse and human enzymes respectively). Produces a hypotensive effect in angiotensin-infused rats and prevents and reverses cardiac hypertrophy in mouse models of cardiac hypertrophy.

Citations: 3

[View Online »](#)

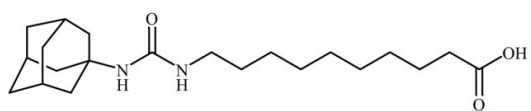
Ordering Information

[Order Online »](#)

BML-EI392-0050	50mg
----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store solutions at -20°C for up to 3 months.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	12-(3-Adamantan-1-yl-ureido)-dodecanoic acid
Appearance	White solid.
CAS	479413-70-2
Couple Target	Epoxide hydrolase
Couple Type	Inhibitor
Formula	$C_{21}H_{36}N_2O_3$
MW	392.5753
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (10mg/ml) or 100% ethanol (4mg/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)