

(±)11,12-Epoxyeicosa-5Z,8Z,14Z-trienoic acid

Inhibits Na⁺K⁺-ATPase

Biosynthesis *via* liver and renal cytochrome P-450. Induces glucagon release from pancreatic islets. Vasodilator in intestinal microcirculation. Inhibits Na⁺K⁺-ATPase. Urinary excretion in humans is increased during pregnancy-induced hypertension.

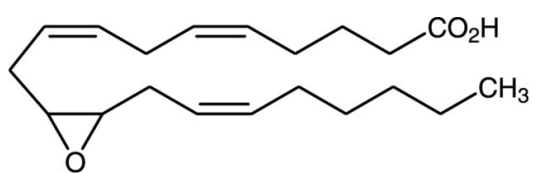
Ordering Information

[Order Online »](#)

BML-EE11-0050	50µg
---------------	------

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

CAS 123931-40-8

Couple Target ATPase

Couple Type Inhibitor

Formulation Supplied at 50 µg/ml in ethanol.

Purity ≥98%

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-
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)