

Acivicin

Irreversible inhibitor of γ -glutamyl transpeptidase (GGT) $ID_{50} = 0.54$ mM. Inhibits the enzymatic conversion of leukotriene C_4 to leukotriene D_4 . Displays antitumor and anti-metastatic activity. Blocks *H. Pylori* GGT-induced T-cell inhibition. Inhibits bioactivation/release of nitric oxide from GSNO.

Citations: 11

[View Online »](#)

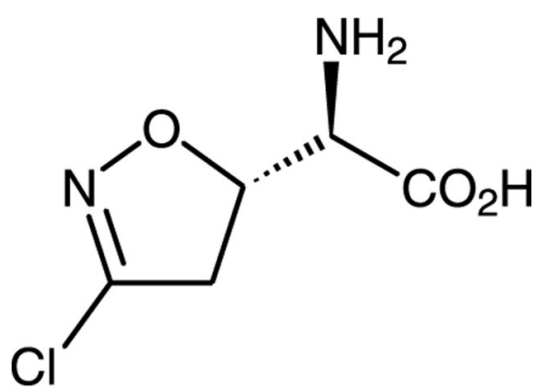
Ordering Information

[Order Online »](#)

BML-EI113-0010	10mg
----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name U-42126, AT-125

Appearance White solid.

CAS 42228-92-2

Couple Target gamma-Glutamyl transpeptidase

Couple Type Inhibitor

Formula $C_5H_7ClN_2O_3$

MW 178.6

Purity ≥98% (HPLC)

Solubility Soluble in water (max. 18mg/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)