

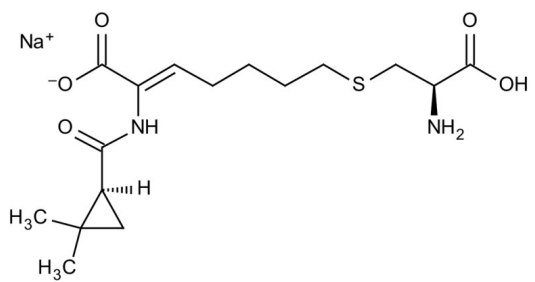
Cilastatin Sodium

DPP inhibitor

Cilastatin is a brush border dipeptidase inhibitor. It inhibits the conversion of leukotriene D4 to leukotriene E4. Nephroprotective in cyclosporine A-induced nephrotoxicity.

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 [R-[R*,S*-(Z)]]-7-[(2-Amino-2-carboxyethyl)thio]-2- [[(2,2-dimethylcyclopropyl)carbonyl]amino]-2- heptenoic acid monosodium salt

Appearance Off-white solid.

CAS 81129-83-1

Couple Target DPP

Couple Type Inhibitor

Formula C₁₆H₂₅N₂O₅S . Na

MW 380.5

Purity ≥98%

Solubility Soluble in water (25mg/ml) or DMSO (10mg/ml at 37°C).

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com