

# Ampicillin . sodium salt

Antibiotic

$\beta$ -Lactam antibiotic. Inhibits bacterial cell-wall synthesis (peptidoglycan cross-linking). Inactivates transpeptidases on the inner surface of the bacterial cell membrane. Effective against Gram-negative and Gram-positive bacteria. Is inactivated by  $\beta$ -lactamase.

Citations: 8

[View Online »](#)

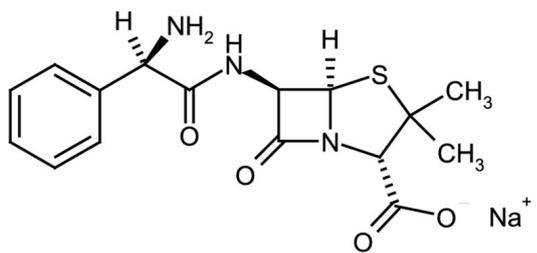
Ordering Information

[Order Online »](#)

ALX-380-268-G005	5g
------------------	----

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	D-(-)-α-Aminobenzylpenicillin . Na
Appearance	White to light yellow powder.
CAS	69-52-3
Couple Target	Transpeptidase
Couple Type	Inhibitor
Formula	$C_{16}H_{18}N_3O_4S \cdot Na$
MI	14: 586
MW	348.4 . 23.0
Purity	Assay (anhydrous basis): 845-988 µg/mg.
RTECS	XH8400000
Solubility	Soluble in water (50mg/ml).
Source	Semisynthetic from penicillin.

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)