

Carbenicillin . disodium salt

Antibiotic

Analog to ampicillin (Prod. No. ALX-380-268). Inhibits bacterial cell-wall synthesis (peptidoglycan cross-linking) by inactivating transpeptidases on the inner surface of the bacterial cell membrane. Effective against Gram-positive and Gram-negative bacteria and *Pseudomonas*.

Citations: 4

[View Online »](#)

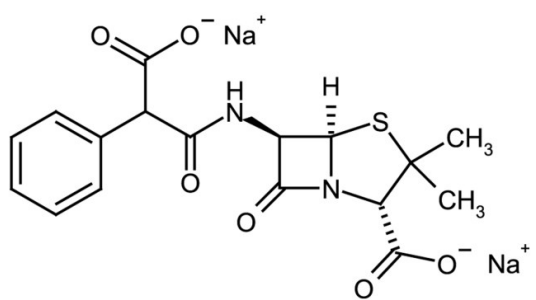
Ordering Information

[Order Online »](#)

ALX-380-270-G001	1g
------------------	----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Hygroscopic.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	α -Carboxybenzylpenicillin . 2Na
Appearance	White to off-white powder.
CAS	4800-94-6
Couple Target	Transpeptidase
Couple Type	Inhibitor
Formula	$C_{17}H_{16}N_2O_6S \cdot 2Na$
MI	14: 1792
MW	376.4 . 46.0
Purity	≥92% (Assay)
RTECS	ON9105000
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)