

Zoledronic acid

Acid sphingomyelinase inhibitor

Inhibitor of osteoclastic bone resorption. Prevents osteoporosis and skeletal fractures, particularly in patients with multiple myeloma or prostate cancer. Also used to treat hypercalcemia. Inhibition of acid sphingomyelinase (aSMase) with an IC_{50} 5 μ m.

Citations: 13

[View Online »](#)

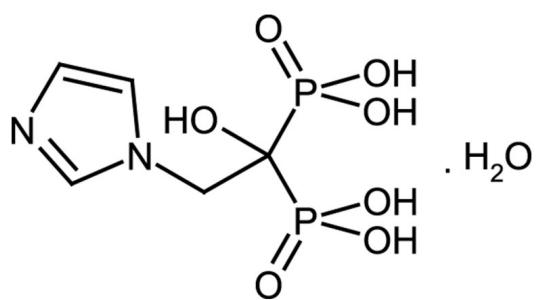
Ordering Information

[Order Online »](#)

ALX-430-153-M025	25mg
ALX-430-153-M100	100mg

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	1-Hydroxy-2-(1H-imidazol-1-yl)ethylidene]bisphosphonic acid, CGP 42446
Appearance	White to off-white powder.
CAS	165800-06-6
Couple Target	Sphingomyelinase
Couple Type	Inhibitor
Formula	$C_5H_{10}N_2O_7P_2 \cdot H_2O$
MI	14: 10187
MW	272.1 . 18.0
Purity	≥98% (HPLC)
Solubility	Soluble in 0.1N NaOH. Sparingly soluble in water.
Source	Synthetic.

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-fr@enzolifesciences.com

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

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

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