

Tobramycin

Antibiotic

Aminoglycoside antibiotic with similar structure to the kanamycins and the gentamycins. Used in the treatment of respiratory infection and cystic fibrosis.

Citations: 3

[View Online »](#)

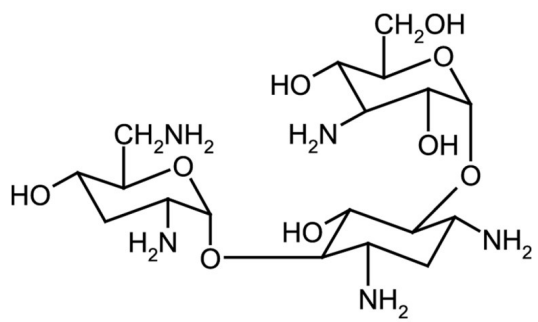
Ordering Information

[Order Online »](#)

ALX-380-018-M250	250mg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Activity	≥900µg/mg
Alternative Name	Deoxykanamycin B, Distobram, Gernebcin, Obramycin, Tobradistin
Appearance	White to off-white powder.
CAS	32986-56-4
Formula	$C_{18}H_{37}N_5O_9$
Formulation	Free base.
MI	14: 9490
MW	467.5
RTECS	WK2100000
Solubility	Soluble in water; very slightly soluble in 100% ethanol.
Source	Isolated from <i>Streptomyces tenebarius</i> .

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