

# Hygromycin B (sterile-filtered solution in PBS)

## Protein synthesis inhibitor

Antibiotic. Inhibits growth of prokaryotic and eukaryotic microorganisms and mammalian cells. Inhibits protein synthesis at the translocation step on the 70S ribosome and causes misreading of the mRNA. Penetrates cells that have been permeabilized by virus infection, hence it can act as an effective antiviral agent.

Citations: 17

[View Online »](#)

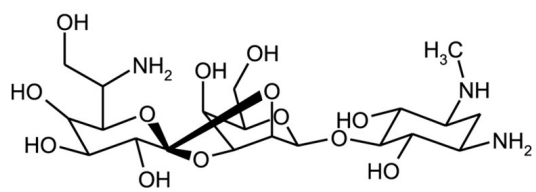
## Ordering Information

[Order Online »](#)

ALX-380-306-G001	1g
------------------	----

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Working hygromycin solutions ( $\leq 2\text{mg/ml}$ ) are stable for up to 1 month when stored at $+4^{\circ}\text{C}$ .
Handling	Protect from light.
Long Term Storage	$+4^{\circ}\text{C}$
Shipping	Ambient Temperature – Dangerous Good

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	O-6-Amino-6-deoxy-L-glycero-D-galacto-heptopyranosylidene-(1-2-3)-O- $\beta$ -D-talopyranosyl(1-5)-2-deoxy-N3-methyl-D-streptamine
Appearance	Liquid. In PBS.
CAS	31282-04-9
Formula	$\text{C}_{20}\text{H}_{37}\text{N}_3\text{O}_{13}$
MI	14: 4852
MW	527.5
Purity	$\geq 90\%$ (HPLC)
Purity Detail	Sterile filtered.
RTECS	WK2130000
Source	Isolated from <i>Streptomyces hygroscopicus</i> .

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-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)

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