

Product datasheet

Neomycin ELISA Kit ab285276

[1 Image](#)

Overview

Product name	Neomycin ELISA Kit
Detection method	Colorimetric
Sample type	Milk, Tissue sections
Assay type	Competitive
Sensitivity	= 0.1 ng/ml
Range	0.1 ng/ml - 8.1 ng/ml
Assay duration	Multiple steps standard assay
Product overview	<p>Neomycin is an aminoglycoside antibiotic found in many topical medications such as creams, ointments, and eyedrops. It is typically used as a topical preparation, such as Neosporin. It can also be given orally, where it is usually combined with other antibiotics. Neomycin, like other aminoglycosides, has been shown to be ototoxic causing tinnitus, hearing loss, and vestibular problems in a small number of patients. Standard techniques/instruments (HPLC or GC-MS) are utilized to detect neomycin. However, these techniques are complex, expensive, laborious, and time-consuming. Immunoassay techniques, such as ELISAs are commonly preferred as a simple, reliable and rapid method for the quantification of neomycin in various samples. The Neomycin ELISA Kit (ab285276, E4611) is a competitive-based ELISA that can be used for the determination of this antibiotic in tissue, honey, milk and milk powder.</p>
Notes	<p>This product is manufactured by BioVision, an Abcam company and was previously called E4611 Neomycin ELISA Kit. E4611-100 is the same size as the 96 test size of ab285276.</p>
Platform	Microplate (12 x 8 well strips)

Properties

Storage instructions Store at +4°C. Please refer to protocols.

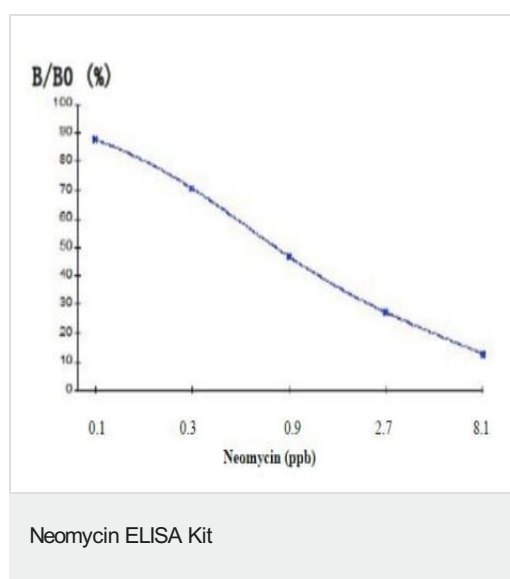
Components	96 tests	96 tests
20X Wash buffer	1 x 15ml	1 x 15ml
Antibody Working Solution	1 x 7ml	1 x 7ml

Components	96 tests	96 tests
Redissolving Solution (2X)	2 x 50ml	2 x 50ml
Enzyme Conjugate	1 x 7ml	1 x 7ml
Micro ELISA Plate	1 unit	1 unit
Standard (S0 / S6)	7 x 1ml	7 x 1ml
Stop Solution	1 x 7ml	1 x 7ml
Substrate A	1 x 7ml	1 x 7ml
Substrate B	1 x 7ml	1 x 7ml

Relevance

Neomycin is an aminoglycoside broad spectrum antibiotic effective against both gram positive and gram negative bacteria. It interferes with protein synthesis in sensitive bacterial cells such as species of Proteus and Staphylococcus.

Images



Typical Standard Curve: These standard curves are for demonstration only. A standard curve must be run with each assay.

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you

- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors