

Neomycin ELISA kit

Simple, fast, and sensitive method for detection of Neomycin in cell culture and tissue

The Neomycin ELISA kit is a colorimetric competitive enzyme immunoassay kit with results in less than 1 hour. Absorbance is read at 450 or 630 nm. Very low reactivity with other antibiotics provides confidence in assay results. Sensitive quantification of 0.5 ng/mL neomycin. Reagents and protocol for sample extraction is included in the kit.

Neomycin is an antibiotic that displays bactericidal activity against gram-negative bacteria and is partially effective against gram positive bacteria. Neomycin belongs to the aminoglycoside class of antibiotics that contain two or more amino sugars connected by glycosidic bonds. All aminoglycosides are potentially toxic compounds. Neomycin is included, in small quantities, as a preservative in some vaccines.

Ordering Information

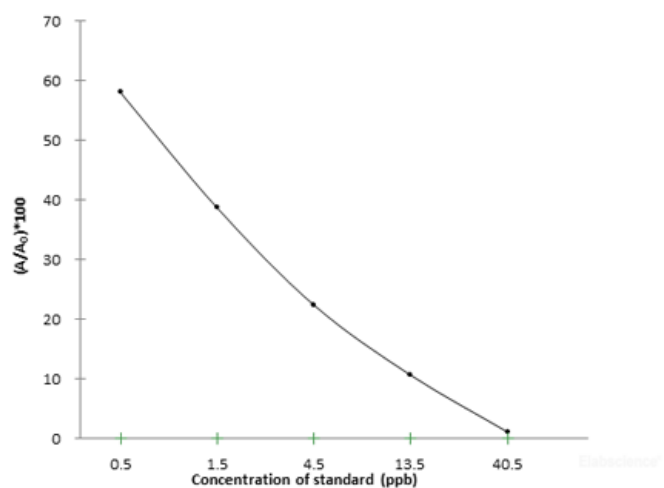
[Order Online »](#)

ENZ-KIT217-0001	96 wells
-----------------	----------

Manuals, SDS & CofA

[View Online »](#)

- Highly specific ELISA which does not detect streptomycin, kanamycin or gentamicin, providing confidence in assay results
- Sensitive measurement, detecting as little as 0.5 ppb (ng/mL) neomycin
- Cost effective and fast, with results in less than 1h
- Easy-to-use liquid color-coded reagents and pre-coated plates minimize error



Standard curve of Neomycin ELISA Kit

Handling & Storage

Use/Stability	Store all components at 2-8°C
Handling	Do not freeze
Long Term Storage	+4°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application	Colorimetric detection, ELISA
Application Notes	For the quantitative determination of neomycin in cell culture supernatants, muscle, milk, egg.
Assay Time	45 min
Compatibility	This product is compatible with the Absorbance 96 Plate Reader .
Contents	Microtiter plate, Conjugate, Antibody Working Solution, Substrate Reagents, Concentrated Sample Solution, Wash Buffer Concentrate, Standard, Stop solution. Additional: Concentrated Tissue Extraction Solution
Crossreactivity	Streptomycin, Kanamycin, Gentamicin < 0.1%
Sensitivity	0.5 ppb (ng/ml)
Species Reactivity	Species independent
Wavelength	450nm / 630nm

Last modified: June 13, 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)