

# Gentamicin ELISA kit

Competitive ELISA for rapid detection of gentamicin in cell culture supernatants and tissue samples

Gentamicin ELISA kit rapidly and accurately analysis Gentamicin residues in samples with results in 45 minutes. Detection limit: Culture Supernatants - 0.12ppb; Muscle - 5ppb; Raw milk/Finished milk - 15ppb

Gentamicin is active against a wide range of bacterial infections, mostly Gram-negative bacteria including *Pseudomonas*, *Proteus*, *Escherichia coli*, *Klebsiella pneumoniae*, *Enterobacter aerogenes*, *Serratia*, and the Gram-positive *Staphylococcus*. The antibiotic is collected from the culture of the Micromonospora by perforating the cell wall of the bacterium.

## Ordering Information

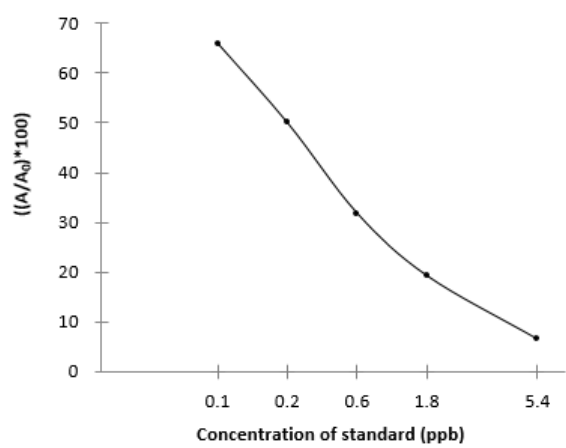
[Order Online »](#)

ENZ-KIT218-0001	96 wells
-----------------	----------

## Manuals, SDS & CofA

[View Online »](#)

- Specific - antibody specifically binds to gentamicin
- Sensitive - measure as little as 0.1ng/ml
- Rapid - analyze samples in 45 minutes



## 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	ELISA
Application Notes	For the quantitative and sensitive analysis of gentamicin in cell culture supernatants, tissue (muscle), milk.
Assay Time	45 minutes
Compatibility	This product is compatible with the <a href="#">Absorbance 96 Plate Reader</a> .
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, Neomycin, Kanamycin < 0.01%
Sensitivity	0.1 ppb (ng/ml)
Specificity	Gentamicin – 100%
Wavelength	450 nm/630 nm



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

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

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