

# 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

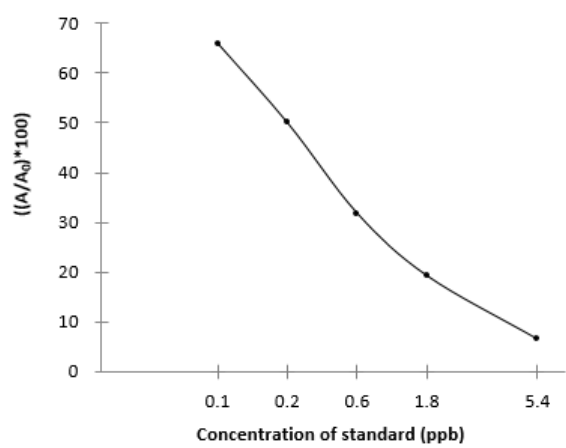
[Order Online »](#)

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 [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, Neomycin, Kanamycin < 0.01%

**Sensitivity** 0.1 ppb (ng/ml)

**Specificity** Gentamicin – 100%

**Wavelength** 450 nm/630 nm

**Last modified: June 18, 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-  
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)