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 Grampositive *Staphylococcus*. The antibiotic is collected from the culture of the Micromonospora by perforating the cell wall of the bacterium.

Ordering Information

Order Online »

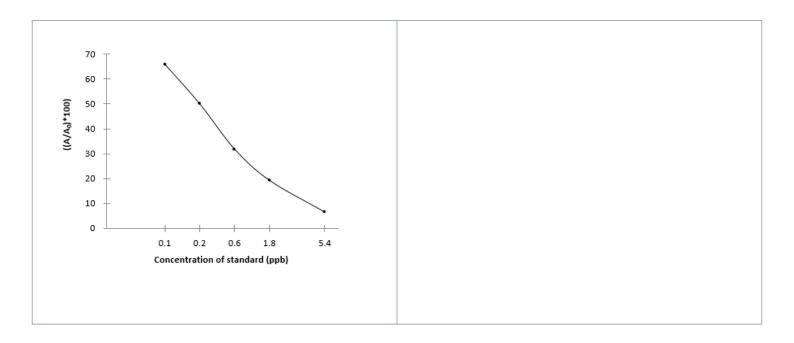
ENZ-KIT218-0001

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 NotesFor 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)

Species Reactivity Species independent

Specificity Gentamicin – 100%

Wavelength 450 nm/630 nm



European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland Phone (UK customers): 0845 601 1488 Phone: +44 1392 825900 info-

uk@enzolifesciences.com