

Streptomycin ELISA kit

Competitive immunoassay for the detection of streptomycin in cell culture supernatants and tissue samples

The Streptomycin EIA kit is a colorimetric competitive enzyme immunoassay kit with results in 75 minutes. Absorbance is read at 450 or 630 nm. Very low reactivity with other antibiotics provides confidence in assay results. Sensitive quantification of 0.1 ng/mL Streptomycin. Reagents and protocol for sample extraction is included in the kit.

Streptomycin is an antibiotic medication used to treat a number of bacterial infections. Streptomycin is an aminoglycoside. It works by blocking the ability of 30S ribosomal subunits to make proteins, which results in bacterial death. Streptomycin, in combination with penicillin, is used in a standard antibiotic cocktail to prevent bacterial infection in cell culture. When purifying protein from a biological extract, streptomycin sulfate is sometimes added as a means of removing nucleic acids.

Ordering Information

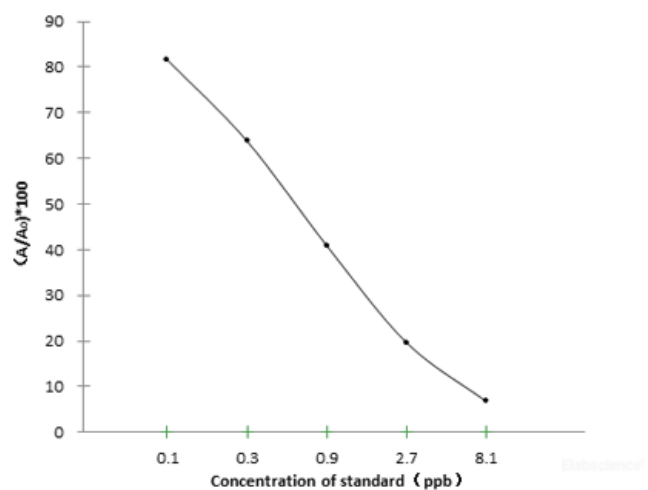
[Order Online »](#)

ENZ-KIT219-0001	96 wells
-----------------	----------

Manuals, SDS & CofA

[View Online »](#)

- Sensitive measurement, detecting as little as 0.1 ppb (ng/mL) streptomycin
- Cost effective and fast, with results in 75 minutes
- Easy-to-use liquid color-coded reagents and pre-coated plates minimize error



Standard curve of Streptomycin 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 Streptomycin in cell culture supernatants, tissue (muscle), milk, milk powder, honey.
Assay Time	75 min
Compatibility	This product is compatible with the Absorbance 96 Plate Reader .
Contents	Microtiter plate, Conjugate, Antibody Working Solution, Substrate Reagents, Wash Buffer Concentrate, Standard, Stop solution, Reconstitution Buffer
Crossreactivity	Clarithromycin – 6.5%, Gentamicin – 2.5%
Sensitivity	0.1 ppb (ng/ml)
Species Reactivity	Species independent
Specificity	Streptomycin – 100%, Dihydrostreptomycin – 100%
Wavelength	450 nm/630nm



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