

Substance P ELISA kit

Highly sensitive Substance P ELISA kit enabling detection of Substance P in plasma, serum, saliva, urine, and supernatants in just 3 hours.

The Substance P ELISA kit is a colorimetric competitive enzyme immunoassay kit with results in 3 hours. Absorbance is read at 405 nm. This kit utilized low sample volume and provides the ability to assay up to 38 samples in duplicate per 96 well kit.

Substance P is an undecapeptide hormone (RPKPQQFFGLM) that mediates a number of biological activities including nausea, anxiety, pain, and vasodilation. It is a member of the tachykinin family of neuropeptides, and is widely distributed in the peripheral and central nervous systems of vertebrates.

Citations: 31

[View Online »](#)

Ordering Information

[Order Online »](#)

ADI-901-018A

5x96 wells

Manuals, SDS & CofA

[View Online »](#)

- Highly sensitive assay, measuring as little as 5.3 pg/ml of Substance P
- Widely cited in numerous peer reviewed publications
- High throughput format, assay up to 38 samples in duplicate per 96 well kit
- Fully quantitative results that surpass semi-quantitative Western blot analysis
- Save with bulk kit package!

Handling & Storage

Use/Stability Store components at temperatures indicated on the component labeling. The Antibody, Conjugate and Standard are stored at -20°C and the remaining components are stored at 4°C. For convenience upon receipt, the unopened kit containing all components can be stored at -20°C and are stable at that temperature until the kit's expiration date. After first thaw, store components at temperature indicated on the labeling as indicated above. Freeze-thaw cycles should be limited for optimal kit performance. Please see Reagent Preparation section in the product manual for more detailed component handling information.

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Colorimetric detection, ELISA

Application Notes For the quantitative determination of Substance P in culture supernatants, plasma, serum, saliva, and urine from any species. Cited sample types include tissue.

Assay Time 3 hours

Compatibility This product is compatible with the [Absorbance 96 Plate Reader](#).

Contents GxR Microtiter plate, Conjugate, Antibody, Assay buffer concentrate, Wash buffer concentrate, Standard, pNpp Substrate, Stop solution

Sensitivity 5.3 pg/ml (range 9.76 – 10,000 pg/ml)

Species Reactivity Species independent

UniProt ID P20366

Wavelength 405 nm



Last modified: August 18, 2024

ENZO LIFE SCIENCES

Phone: 800.942.0430

info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG

Phone: +41 61 926 8989

info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg

Phone: +32 3 466 0420

info-be@enzolifesciences.com

France
Phone: +33 472 440 655

info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526

info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):

0845 601 1488
Phone: +44 1392 825900

info-uk@enzolifesciences.com