

# 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: 32

[View Online »](#)

## Ordering Information

[Order Online »](#)

ADI-900-018A

96 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

# 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 type includes tissue.

**Assay Time** 3 hours

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

**Contents** GxR IgG Microtiter plate, Conjugate, Antibody, Assay buffer, 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