
Substance P Polyclonal Antibody

(Catalog No. A-0638)

Background

Substance P (SP) is an undecapeptide (a peptide composed of a chain of 11 amino acid residues) member of the tachykinin neuropeptide family. Tachykinins are active peptides which excite neurons, evoke behavioral responses, are potent vasodilators and secretagogues, and contract (directly or indirectly) many smooth muscles.

Description

Rabbit polyclonal antibody to Substance P

Formulation

Liquid. Antiserum containing 0.1% sodium azide.

Specificity

Rat, Human, Pig, Guinea pig

Uniprot ID

P20366

Purification

Serum

Immunogen

Synthetic substance P (H-RPKPQQFFGLM-NH₂)

Storage

Store at -20°C. Avoid multiple freeze/thaw cycles.

Application

IHC

Ordering Information

Products	Size	Cat. No.
Substance P Polyclonal Antibody	100 µl	A-0638-100

This product is for research purposes only. Not intended for use in diagnostic procedures.