
Somatostatin 14 Polyclonal Antibody

(Catalog No. A-0636)

Background

Somatostatin is a peptide hormone that has two active forms produced by alternative cleavage of a single preproprotein: one of 14 amino acids and the other of 28 amino acids. Expressed throughout the body, this hormone hinders the release of numerous secondary hormones by binding to high-affinity G-protein-coupled somatostatin receptors. Somatostatin interacts with the pituitary growth hormone, the thyroid stimulating hormone, and nearly all of the gastrointestinal tract hormones, making it an essential regulator of the endocrine system. Somatostatin affects neurotransmission and cell proliferation, and is also responsible for inhibiting insulin and glucagon secretion.

Description

Rabbit polyclonal antibody to somatostatin-14 (SOM-14)

Formulation

Liquid. Neat serum containing 0.01M sodium azide

Immunogen

Synthetic peptide corresponding to aa 15-28 of human somatostatin-14

Specificity

Species Independent. Recognizes the C-terminus of somatostatin-14 in a wide range of mammalian and non-mammalian species, including mouse and human

Purification

Serum

Storage

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

Application

IHC, IMMUNOEM

Ordering Information

Products	Size	Cat. No.
Somatostatin 14 Polyclonal Antibody	100 µl	A-0636-100

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