
TRPA1 (extracellular) Polyclonal Antibody (Cat. No A-0656)

Background

TRPA1 (Transient receptor potential cation channel subfamily A member 1) is expressed in peripheral sensory neurons, where it is suggested to be involved in the detection of pain and possibly cold perception. The structure of TRPA1 is highly related to both the protein ankyrin and transmembrane proteins. The specific function of this protein has not yet been determined; however, studies indicate the function may involve a role in signal transduction and growth control. Diseases related to this gene include episodic pain syndrome, familial and familial episodic pain syndrome with predominantly upper body involvement.

Description

Rabbit polyclonal antibody to TRPA1 (extracellular)

Formulation

Lyophilized from PBS, pH 7.4, 1% BSA and 0.05% sodium azide.

Immunogen

Synthetic peptide corresponding to aa 747-760 of human TRPA1.

Specificity

Mouse, Rat

Purification

Serum

Storage

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

Alternative Names

Transient receptor potential cation channel subfamily A member 1, Ankyrin-like protein with transmembrane domains 1

Application

IHC, WB

Ordering Information**Products**

TRPA1 (extracellular) Polyclonal Antibody

Size

50 µl

Cat. No.

A-0656-050

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