
TRPV2 Polyclonal Antibody, FITC Conjugated

(Catalog # A69357)

Background

is not activated by vanilloids and acidic pH.

Description

TRPV2 Polyclonal Antibody, FITC Conjugated. FITC. Raised in: Rabbit.

Formulation

Liquid. 0.03% Proclin 300, 50% Glycerol, 0.01M PBS, pH 7.4.

Specificity

Human

Isotype

IgG

Uniprot ID

Q9Y5S1

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Transient receptor potential cation channel subfamily V member 2 protein (1-117AA)

Storage

Shipped at 4°C. Upon receipt, store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Transient receptor potential cation channel subfamily V member 2, TrpV2, Osm-9-like TRP channel 2, OTRPC2, Vanilloid receptor-like protein 1, VRL-1, TRPV2, VRL

Application

ELISA; Recommended dilution: ELISA 1:100-1:500
