
S100A2 Polyclonal Antibody, FITC Conjugated

(Catalog # A63368)

Background

May function as calcium sensor and modulator, contributing to cellular calcium signaling. May function by interacting with other proteins, such as TPR-containing proteins, and indirectly play a role in many physiological processes. May also play a role in suppressing tumor cell growth.

Description

S100A2 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

P29034

Purification

Affinity purified

Immunogen

Recombinant Human Protein S100-A2 protein (1-98AA)

Storage

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

Alternative Names

Protein S100-A2, CAN19, Protein S-100L, S100 calcium-binding protein A2, S100A2, S100L

Application

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