
ACTA1 Polyclonal Antibody, FITC Conjugated

(Catalog #A54418)

Background

Actins are highly conserved proteins that are involved in various types of cell motility and are ubiquitously expressed in all eukaryotic cells.

Description

ACTA1 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

P68133

Purification

Protein G purified

Immunogen

Recombinant Human Actin, alpha skeletal muscle protein (3-377AA)

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

Alpha-actin-1 ACTA1 ACTA

Application

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