
EIF3I Polyclonal Antibody, FITC Conjugated

(Catalog # A51667)

Background

EIF3S2 is the largest of the EIFs. It consists of at least 10 nonidentical subunits in mammals. In *S. cerevisiae* the p39 subunit contains WD repeats; these are thought to mediate protein-protein interactions. The p39 protein appears to be essential for maintaining the integrity of the yeast EIF3 complex. The mammalian EIF3-p36 subunit is homologous to yeast p39.

Description

EIF3I 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

Q13347

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Eukaryotic translation initiation factor 3 subunit I protein (1-323AA)

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

EIF3S2, TRIP1

Application

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