
EIF3I Polyclonal Antibody

(Catalog # A51665)

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. Unconjugated. Raised in: Rabbit.

Formulation

0.03% Proclin 300, 50% Glycerol, 0.01M PBS, PH 7.4

Specificity

Human, Rat

Isotype

IgG

Uniprot ID

Q13347

Purification

Protein G purified

Immunogen

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

Storage

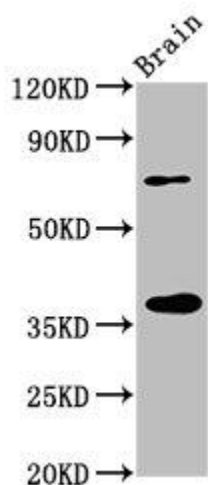
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Eukaryotic translation initiation factor 3 subunit I

Application

ELISA, WB, IF; Recommended dilution: WB:1:2000-1:10000, IF:1:200-1:500



Western Blot

Positive WB detected in: Rat brain tissue

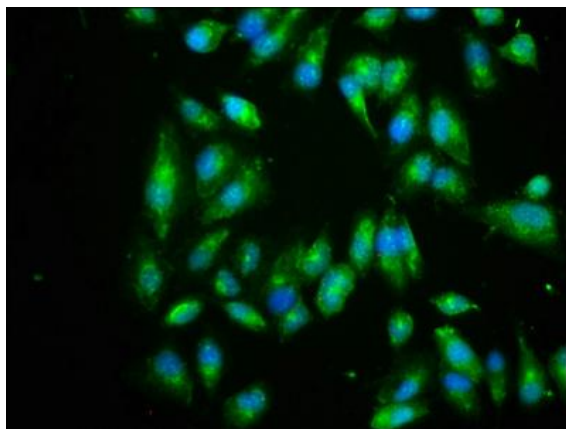
All lanes: EIF3I antibody at 3.2ug/ml

Secondary

Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 37 kDa

Observed band size: 37 kDa



Immunofluorescence staining of HeLa cells with EIF3I Antibody at 1:266, counter-stained with DAPI. The cells were fixed in 4% formaldehyde, permeabilized using 0.2% Triton X-100 and blocked in 10% normal Goat Serum. The cells were then incubated with the antibody overnight at 4°C. The secondary antibody was Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L).