

EIF3G Polyclonal Antibody

(Catalog # A70579)

Description

EIF3G Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Buffer: PBS with 0.02% sodium azide, 50% glycerol, pH7.3.

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

O75821

Purification

Affinity Purification

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 1-320 of human EIF3G (NP_003746.2).

Storage

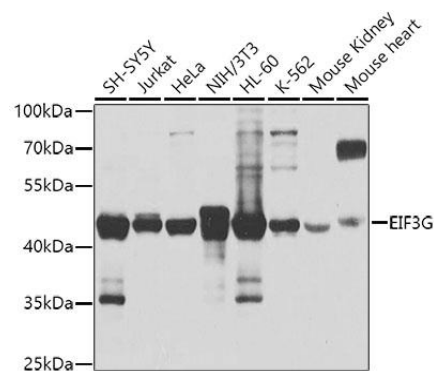
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

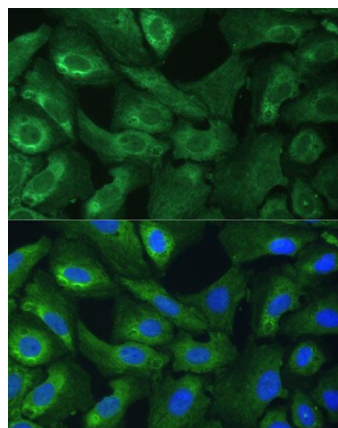
EIF3G; EIF3-P42; EIF3S4; eIF3-delta; eIF3-p44; eukaryotic translation initiation factor 3 subunit G

Application

WB, IF; Recommended dilution: WB 1:200 - 1:2000, IF 1:50 - 1:200



WB analysis of extracts of various cell lines, using EIF3G Polyclonal Antibody at 1:1000 dilution.
Secondary Ab: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution.
Lysates/proteins: 25ug per lane.
Blocking buffer: 3% nonfat dry milk in TBST.
Exposure time: 1s.



Immunofluorescence analysis of U-2 OS cells using EIF3G Polyclonal Antibody at dilution of 1:100 (40x lens). Blue: DAPI for nuclear staining.