

EIF3I Polyclonal Antibody

(Catalog # A70747)

Description

EIF3I Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human

Isotype

IgG

Uniprot ID

Q13347

Purification

Affinity Purification

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 1-325 of human EIF3I (NP_003748.1).

Storage

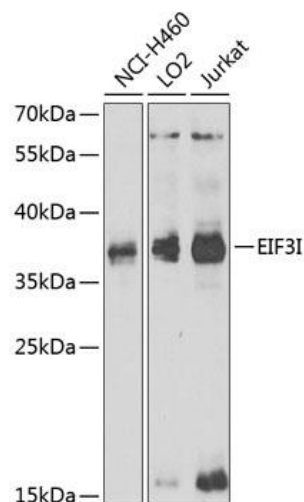
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

EIF3I; EIF3S2; PRO2242; TRIP-1; TRIP1; eIF3-beta; eIF3-p36; eukaryotic translation initiation factor 3 subunit I

Application

WB; Recommended dilution: WB 1:500 - 1:2000



Western blot analysis of extracts of various cell lines, using EIF3I Polyclonal Antibody at 1:1000 dilution. Secondary antibody: 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: 90s.