

EIF5A2 Polyclonal Antibody

(Catalog # A72823)

Description

EIF5A2 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

Q9GZV4

Purification

Affinity Purification

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 1-153 of human EIF5A2 (NP_065123.1).

Storage

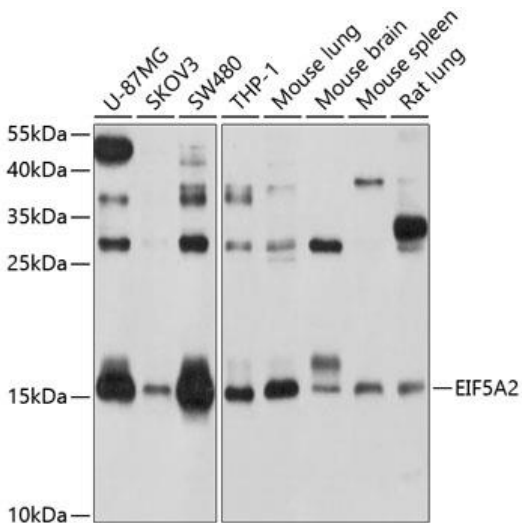
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

EIF5A2; EIF-5A2; eIF5AII; eukaryotic translation initiation factor 5A2

Application

WB; Recommended dilution: WB 1:200 - 1:2000



Western blot analysis of extracts of various cell lines, using EIF5A2 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: 30s.