

CGGBP1 Polyclonal Antibody

(Catalog #A72863)

Description

CGGBP1 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Buffer: PBS with 0.01% thimerosal, 50% glycerol, pH 7.3.

Specificity

Human

Isotype

IgG

Uniprot ID

Q9UFW8

Purification

Affinity Purification

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 1-167 of human CGGBP1 (NP_003654.3).

Storage

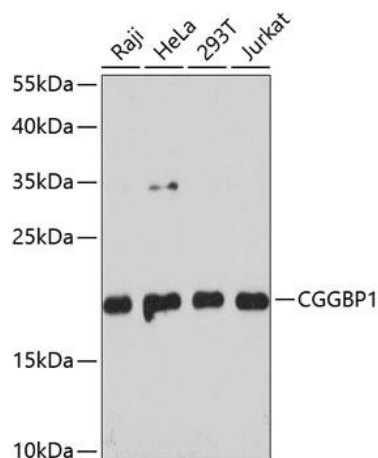
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

CGGBP1; CGGBP; p20-CGGBP; CGG triplet repeat-binding protein 1

Application

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



Western blot analysis of extracts of various cell lines, using CGGBP1 Polyclonal Antibody at 1:3000 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: 60s.