

G2E3 Polyclonal Antibody

(Catalog # A73171)

Description

G2E3 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

Q7L622

Purification

Affinity Purification

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 1-180 of human G2E3 (NP_060239.2).

Storage

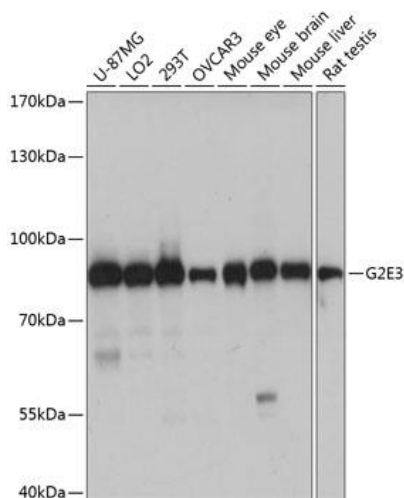
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

G2E3; KIAA1333; PHF7B; G2/M phase-specific E3 ubiquitin-protein ligase

Application

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



Western blot analysis of extracts of various cell lines, using G2E3 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: 1s.