

MPG Polyclonal Antibody

(Catalog #A73371)

Description

MPG Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

P29372

Purification

Affinity Purified

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 1-293 of human MPG (NP_001015052.1).

Storage

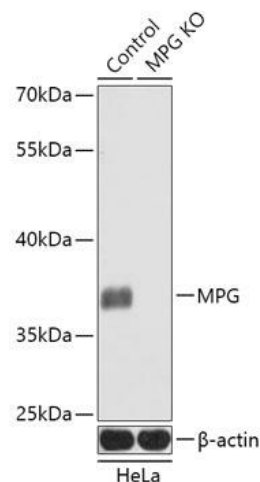
Shipped at 4°C. Store at -20°C. Avoid freeze / thaw cycles.

Alternative Names

MPG; AAG; ADPG; APNG; CRA36.1; MDG; Mid1; PIG11; PIG16; anpg

Application

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



Western blot analysis of extracts from normal (control) and MPG knockout (KO) HeLa cells, using MPG 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: 60s.