
GSS Polyclonal Antibody

(Catalog # A73107)

Background

Glutathione is important for a variety of biological functions, including protection of cells from oxidative damage by free radicals, detoxification of xenobiotics, and membrane transport. The protein encoded by this gene functions as a homodimer to catalyze the second step of glutathione biosynthesis, which is the ATP-dependent conversion of gamma-L-glutamyl-L-cysteine to glutathione. Defects in this gene are a cause of glutathione synthetase deficiency.

Description

GSS Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Buffer: PBS with 0.01% thiomersal, 50% glycerol, pH7.3

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

P48637

Purification

Affinity Purification

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 373-474 of human GSS (NP_000169.1).

Storage

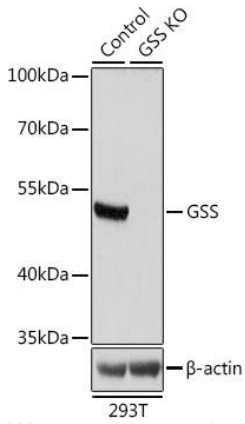
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

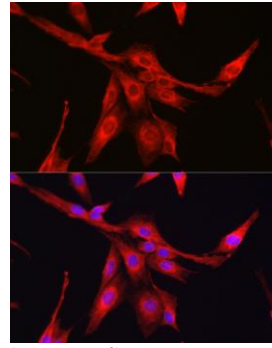
GSS; GSHS; HEL-S-64p; HEL-S-88n; glutathione synthetase

Application

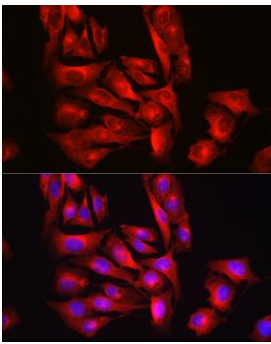
WB, IF/ICC; Recommended Dilutions: WB 1:500 - 1:2000, IF/ICC 1:50 - 1:200



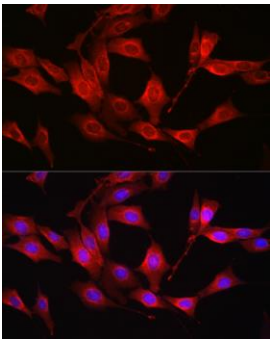
Western blot analysis of extracts from normal (control) and GSS knockout (KO) 293T cells, using GSS Polyclonal Antibody at 1:1000 dilution.
 Secondary Antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution.
 Lysates/proteins: 25ug per lane.
 Blocking buffer: 3% nonfat dry milk in TBST.
 Exposure time: 5s.



Immunofluorescence analysis of PC-12 cells using [KO Validated] Glutathione Synthetase (GSS) Rabbit pAb at dilution of 1:100 (40x lens). Blue: DAPI for nuclear staining.



Immunofluorescence analysis of U2OS cells using [KO Validated] Glutathione Synthetase (GSS) Rabbit pAb at dilution of 1:100 (40x lens). Blue: DAPI for nuclear staining.



Immunofluorescence analysis of NIH/3T3 cells using [KO Validated] Glutathione Synthetase (GSS) Rabbit pAb at dilution of 1:100 (40x lens). Blue: DAPI for nuclear staining.