

Fatty Acid Synthase Polyclonal Antibody

(Catalog # A73225)

Background

The enzyme encoded by this gene is a multifunctional protein. Its main function is to catalyze the synthesis of palmitate from acetyl-CoA and malonyl-CoA, in the presence of NADPH, into long-chain saturated fatty acids. In some cancer cell lines, this protein has been found to be fused with estrogen receptor-alpha (ER-alpha), in which the N-terminus of FAS is fused in-frame with the C-terminus of ER-alpha.

Description

Fatty Acid Synthase Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Buffer: PBS with 0.02% sodium azide, 50% glycerol, pH7.3.

Specificity

Human

Isotype

IgG

Uniprot ID

P49327

Purification

Affinity Purified

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 2212-2511 of human Fatty Acid Synthase (NP_004095.4).

Storage

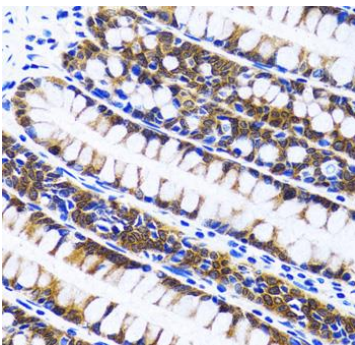
Shipped at 4°C. Store at -20°C. Avoid freeze / thaw cycles.

Alternative Names

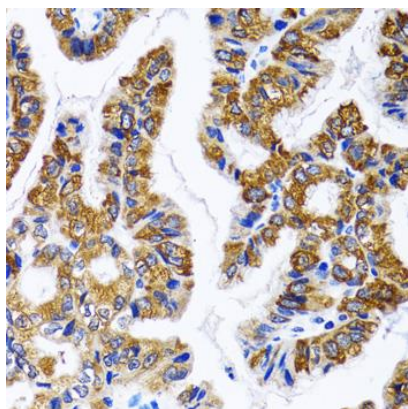
FAS; OA-519; SDR27X1; FASN

Application

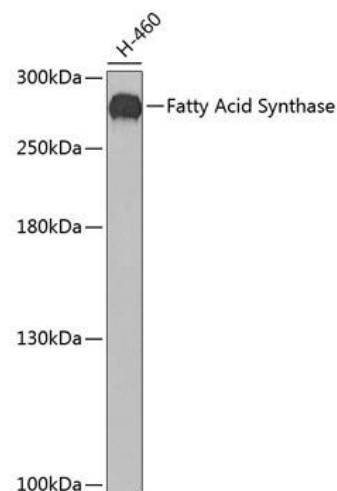
WB, IHC, IF, IP; Recommended dilution: WB 1:500 - 1:2000, IHC 1:50 - 1:100, IF 1:50 - 1:200, IP 1:50 - 1:100



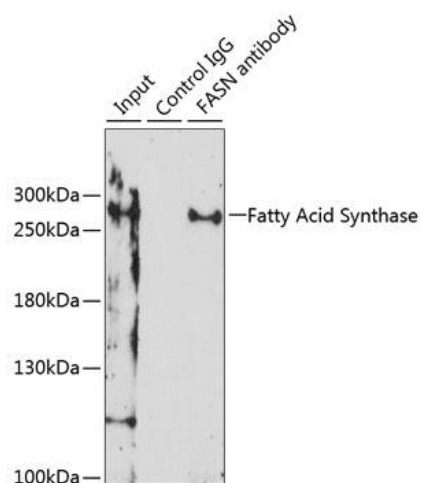
Immunohistochemistry of paraffin-embedded human colon using Fatty Acid Synthase Polyclonal Antibody at dilution of 1:100 (40x lens).



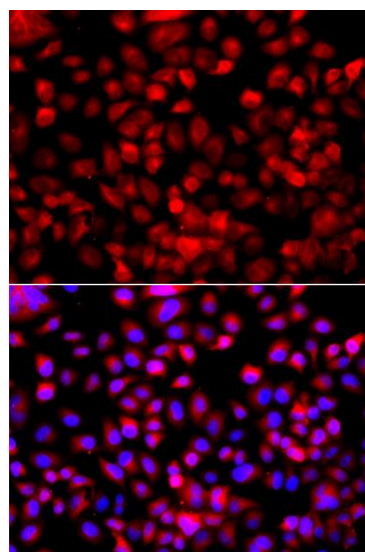
Immunohistochemistry of paraffin-embedded human gastric cancer using Fatty Acid Synthase Polyclonal Antibody at dilution of 1:100 (40x lens).



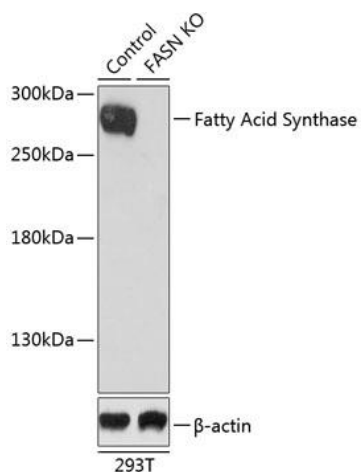
WB of extracts of H460 cells, using Fatty Acid Synthase Polyclonal Antibody at 1:200 dilution. Secondary Antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution. Lysates/proteins: 25ug per lane. Blocking buffer: 3% nonfat dry milk in TBST. Exposure time: 90s.



Immunoprecipitation analysis of 200ug extracts of H460 cells, using 3 ug Fatty Acid Synthase Polyclonal Antibody. Western blot was performed from the immunoprecipitate using Fatty Acid Synthase Polyclonal Antibody at a dilution of 1:1000.



Immunofluorescence analysis of A549 cells using Fatty Acid Synthase Polyclonal Antibody. Blue: DAPI for nuclear staining.



WB of extracts from normal (control) and Fatty Acid Synthase knockout (KO) 293T cells, using Fatty Acid Synthase 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: 90s.