
Fatty Acid Synthase Polyclonal Antibody

(Catalog #A70604)

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 0.09% Sodium azide, 50% glycerol, pH 7.3.

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

P49327

Purification

Affinity Purified

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 2221-2511 of human Fatty Acid Synthase (FASN) (NP_004095.4)

Storage

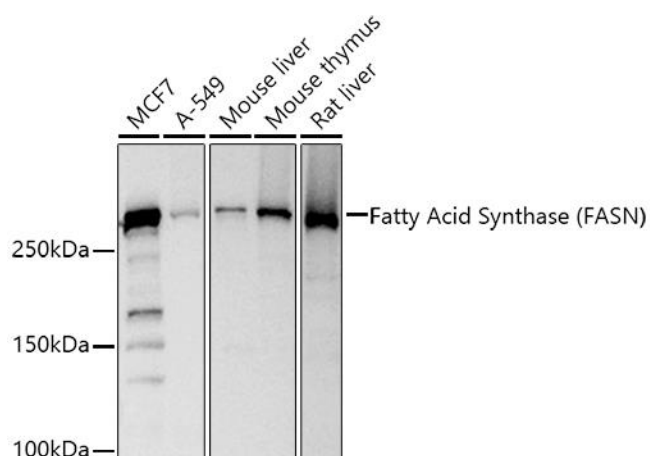
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

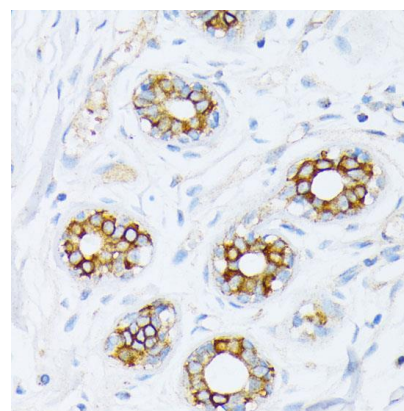
FASN; FAS; OA-519; SDR27X1; fatty acid synthase

Application

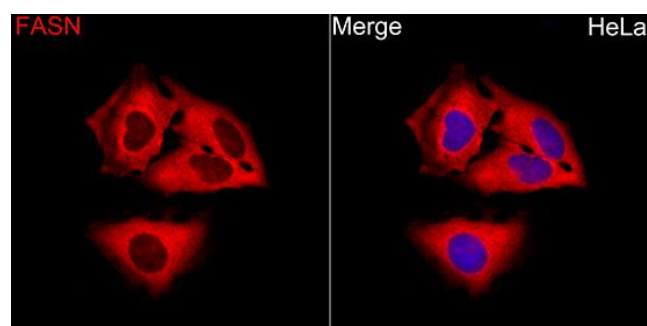
ELISA, WB, IHC, IF; Recommended dilution: WB 1:500 - 1:1000, IHC 1:50 - 1:200, IF 1:50 - 1:200



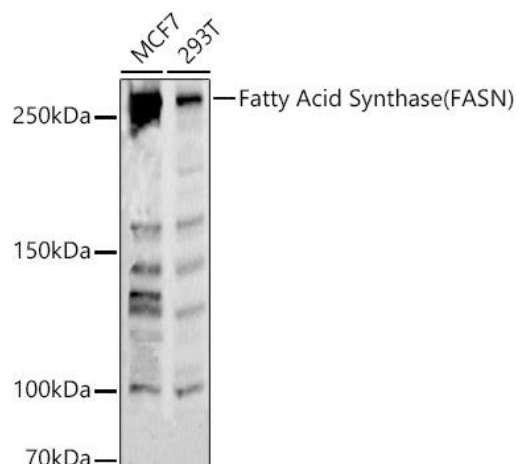
Western blot analysis of various lysates using Fatty Acid Synthase Polyclonal Antibody at 1:1000 dilution.
 Secondary antibody: HRP-conjugated Goat anti-Rabbit IgG (H+L) at 1:10000 dilution.
 Lysates/proteins: 25µg per lane.
 Blocking buffer: 3% nonfat dry milk in TBST.



Immunohistochemistry of paraffin-embedded human breast cancer using Fatty Acid Synthase Polyclonal Antibody at dilution of 1:200 (40x lens).



Immunofluorescence analysis of HeLa cells using Fatty Acid Synthase Polyclonal Antibody at a dilution of 1:200 (40x lens).
 Secondary antibody: Cy3 Goat Anti-Rabbit IgG (H+L) at 1:500 dilution. Blue: DAPI for nuclear staining.



Western blot analysis of various lysates, using Fatty Acid Synthase Polyclonal Antibody at 1:1000 dilution.
 Secondary antibody: HRP-conjugated Goat anti-Rabbit IgG (H+L) at 1:10000 dilution.
 Lysates/proteins: 25µg per lane.
 Blocking buffer: 3% nonfat dry milk in TBST.