
Phospho-ACLY-S455 Polyclonal Antibody

(Catalog # A72329)

Background

ATP citrate lyase is the primary enzyme responsible for the synthesis of cytosolic acetyl-CoA in many tissues. The enzyme is a tetramer (relative molecular weight approximately 440,000) of apparently identical subunits. It catalyzes the formation of acetyl-CoA and oxaloacetate from citrate and CoA with a concomitant hydrolysis of ATP to ADP and phosphate. The product, acetyl-CoA, serves several important biosynthetic pathways, including lipogenesis and cholesterol synthesis. In nervous tissue, ATP citrate-lyase may be involved in the biosynthesis of acetylcholine. Multiple transcript variants encoding distinct isoforms have been identified for this gene.

Description

Phospho-ACLY-S455 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Mouse, Rat

Isotype

IgG

Uniprot ID

P53396

Purification

Affinity Purified

Immunogen

A synthetic phosphorylated peptide around S455 of human ACLY (NP_001087.2).

Storage

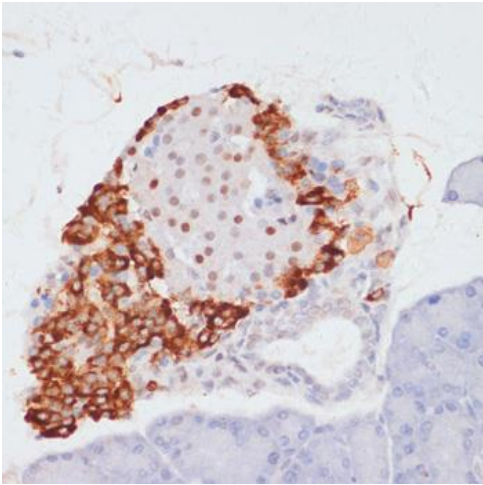
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

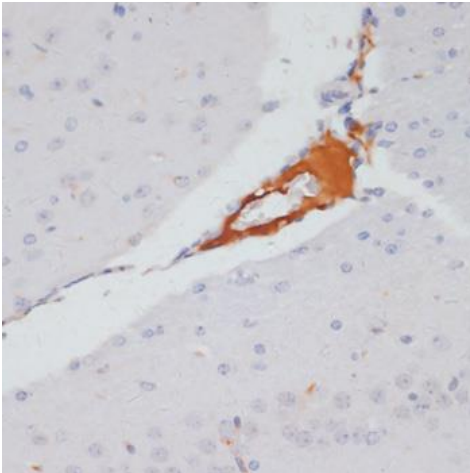
ACLY; ACL; ATPCL; CLATP; ATP citrate lyase

Application

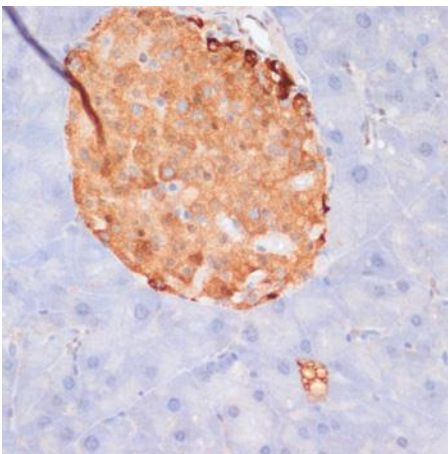
WB, IHC, IP; Recommended dilution: WB: 1:500-1:2000, IHC: 1:50-1:100, IP:1:50-1:100



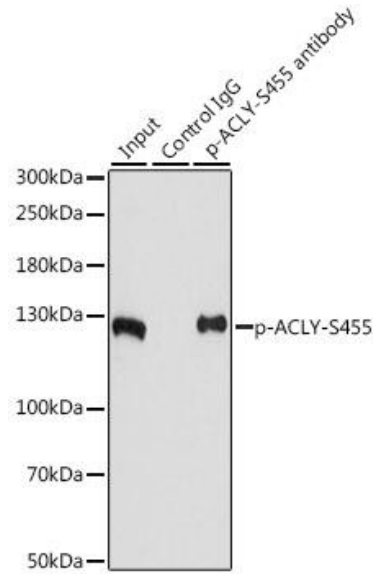
Immunohistochemistry of paraffin-embedded rat pancreas using Phospho-ACLY-S455 Polyclonal Antibody at dilution of 1:100 (40x lens).



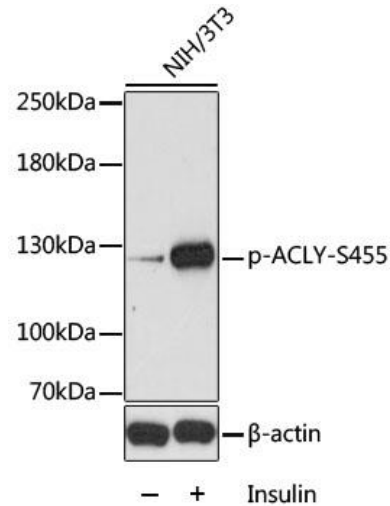
Immunohistochemistry of paraffin-embedded mouse brain using Phospho-ACLY-S455 Polyclonal Antibody at dilution of 1:100 (40x lens).



Immunohistochemistry of paraffin-embedded mouse pancreas using Phospho-ACLY-S455 Polyclonal Antibody at dilution of 1:100 (40x lens).



Immunoprecipitation analysis of 200ug extracts of NIH/3T3 cells, using 3 ug Phospho-ACLY-S455 Polyclonal Antibody. Western blot was performed from the immunoprecipitate using Phospho-ACLY-S455 Polyclonal Antibody at a dilution of 1:1000. NIH/3T3 cells were treated by Insulin (100 nM) at 37°C for 10 minutes after serum-starvation overnight.



Western blot analysis of extracts of NIH/3T3 cells, using Phospho-ACLY-S455 antibody at 1:2000 dilution. NIH/3T3 cells were treated by Insulin (100nM) for 10 minutes after serum-starvation overnight. Secondary antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution. Lysates/proteins: 25ug per lane. Blocking buffer: 3% BSA. Exposure time: 60s.