

FAH Polyclonal Antibody

(Catalog # A62538)

Background

cytosol, extracellular exosome, fumarylacetoacetase activity, cellular nitrogen compound metabolic process, L-phenylalanine catabolic process, small molecule metabolic process, tyrosine catabolic process

Description

FAH Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Liquid. 0.03% Proclin 300, 50% Glycerol, 0.01M PBS, PH 7.4.

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

P16930

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Fumarylacetoacetase protein (2-250AA)

Storage

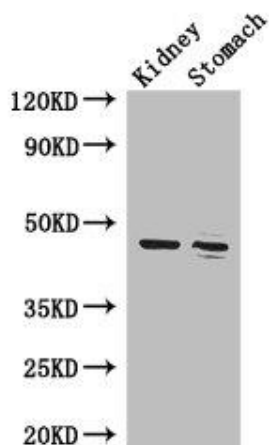
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Fumarylacetoacetase, Beta-diketonase, Fumarylacetoacetate hydrolase, FAH

Application

ELISA, WB, IHC, IF; Recommended dilution: WB:1:2000-1:5000, IHC:1:20-1:200, IF:1:50-1:200



Western Blot

Positive WB detected in: Mouse kidney tissue, Mouse stomach tissue

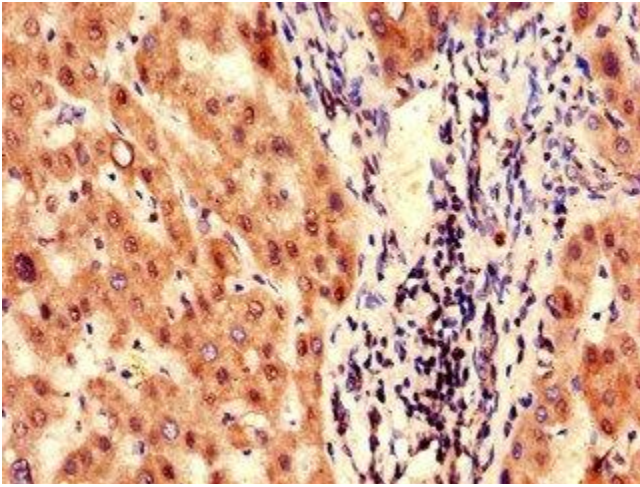
All lanes: FAH Polyclonal Antibody at 3.4ug/ml

Secondary

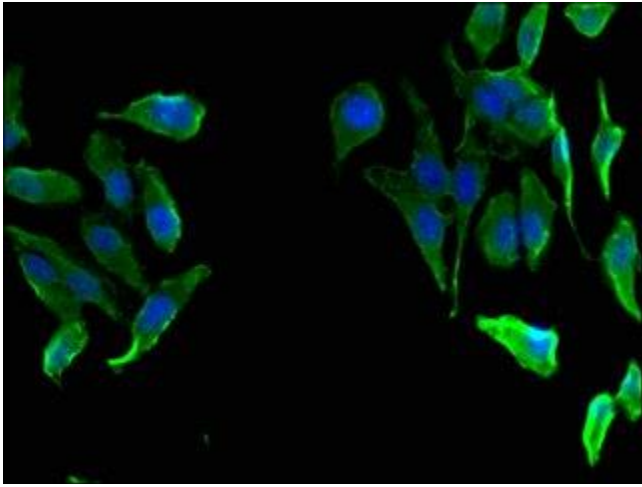
Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 47, 39 kDa

Observed band size: 47 kDa



Immunohistochemistry of paraffin-embedded human liver tissue at dilution of 1:100



Immunofluorescent analysis of HeLa cells at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)