

PAH Polyclonal Antibody

(Catalog # A63090)

Background

cytosol, extracellular exosome, phenylalanine 4-monooxygenase activity, catecholamine biosynthetic process, cellular amino acid biosynthetic process, cellular nitrogen compound metabolic process, L-phenylalanine catabolic process, neurotransmitter biosynthetic process, small molecule metabolic process

Description

PAH Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

P00439

Purification

Protein G purified

Immunogen

Recombinant Human Phenylalanine-4-hydroxylase protein (2-452AA)

Storage

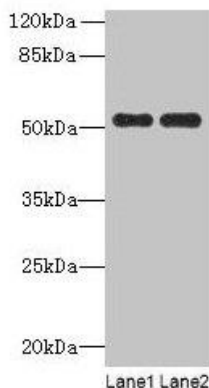
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Phenylalanine-4-hydroxylase, PAH, Phe-4-monooxygenase, PAH

Application

ELISA, WB, IHC, IF; Recommended dilution: WB:1:1000-1:5000, IHC:1:20-1:200, IF:1:50-1:200



Western blot

All lanes: PAH antibody at 2.5ug/ml

Lane 1: Mouse kidney tissue

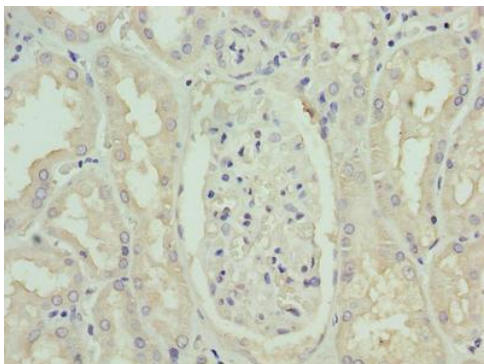
Lane 2: Mouse liver tissue

Secondary

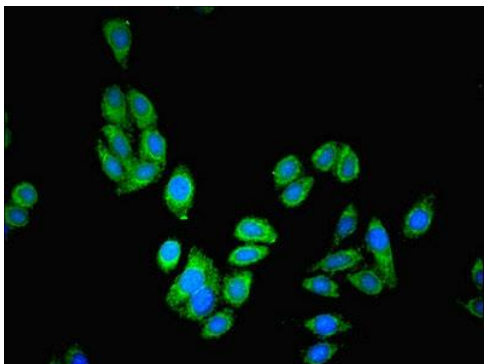
Goat polyclonal to rabbit IgG at 1/10000 dilution

Predicted band size: 52 kDa

Observed band size: 52 kDa



Immunohistochemistry of paraffin-embedded human kidney tissue using PAH Antibody at dilution of 1:100



Immunofluorescent analysis of HepG2 cells using PAH Antibody at dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)