

FAHD2A Polyclonal Antibody

(Catalog # A66294)

Background

May have hydrolase activity.

Description

FAHD2A Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human

Isotype

IgG

Uniprot ID

Q96GK7

Purification

Affinity purified

Immunogen

Recombinant Human Fumarylacetoacetate hydrolase domain-containing protein 2A protein (4-314AA)

Storage

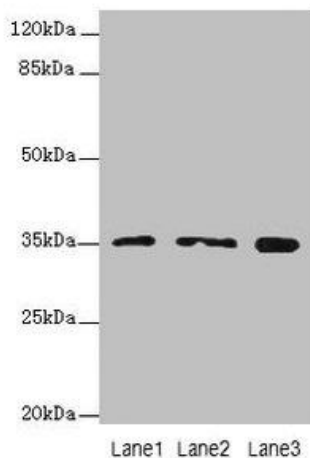
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Fumarylacetoacetate hydrolase domain-containing protein 2A, FAHD2A, CGI-105

Application

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



Western blot

All lanes: FAHD2A Polyclonal Antibody at 0.5µg/ml

Lane 1: K562 whole cell lysate

Lane 2: U87 whole cell lysate

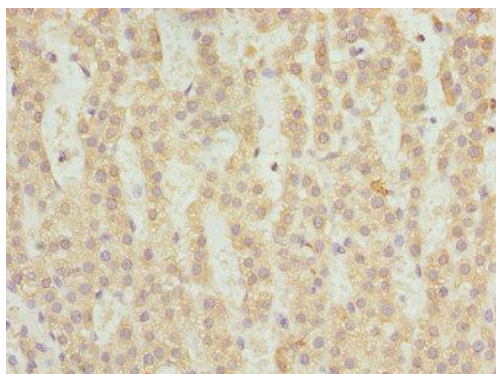
Lane 3: A549 whole cell lysate

Secondary

Goat polyclonal to rabbit IgG at 1/10000 dilution

Predicted band size: 35 kDa

Observed band size: 35 kDa



Immunohistochemistry of paraffin-embedded human adrenal gland tissue using FAHD2A Polyclonal Antibody at dilution of 1:100