

PHGDH Polyclonal Antibody

(Catalog # A53941)

Background

Catalyzes the reversible oxidation of 3-phospho-D-glycerate to 3-phosphonooxypyruvate, the first step of the phosphorylated L-serine biosynthesis pathway. Also catalyzes the reversible oxidation of 2-hydroxyglutarate to 2-oxoglutarate and the reversible oxidation of (S)-malate to oxaloacetate.

Description

PHGDH 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

O43175

Purification

>95%, Protein G purified

Immunogen

Recombinant Human D-3-phosphoglycerate dehydrogenase protein (2-251AA)

Storage

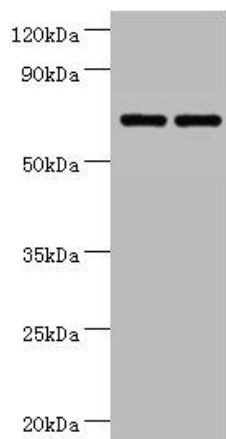
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

PGDH3, PHGDH, 3-PGDH, D-3-phosphoglycerate dehydrogenase

Application

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



Western blot

All lanes: PHGDH Polyclonal Antibody at 2ug/ml

Lane 1: EC109 whole cell lysate

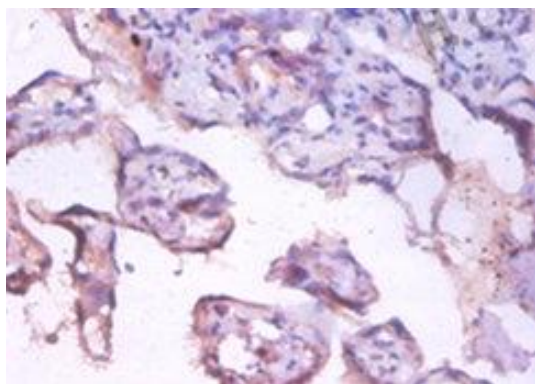
Lane 2: 293T whole cell lysate

Secondary

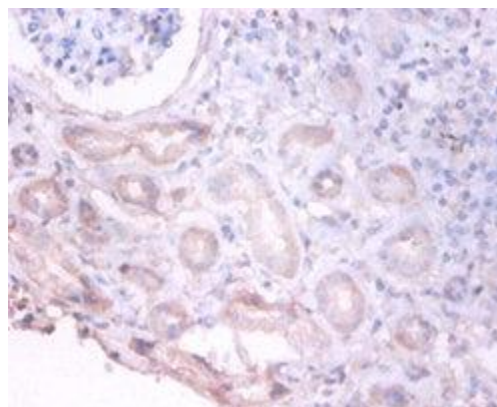
Goat polyclonal to rabbit at 1/10000 dilution

Predicted band size: 59kDa

Observed band size: 59kDa



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



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