

HKDC1 Polyclonal Antibody

(Catalog # A64866)

Background

cytosol, fructokinase activity, glucokinase activity, mannokinase activity, cellular glucose homeostasis, glucose 6-phosphate metabolic process, glycolytic process

Description

HKDC1 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

Q2TB90

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Putative hexokinase HKDC1 protein (129-427AA)

Storage

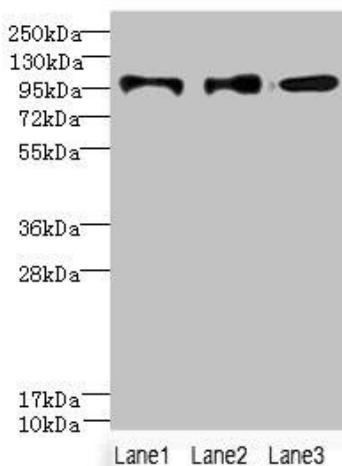
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Putative hexokinase HKDC1, Hexokinase domain-containing protein 1, HKDC1

Application

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



Western blot

All lanes: HKDC1 Polyclonal Antibody at 4µg/ml

Lane 1: U87 whole cell lysate

Lane 2: Caco-2 whole cell lysate

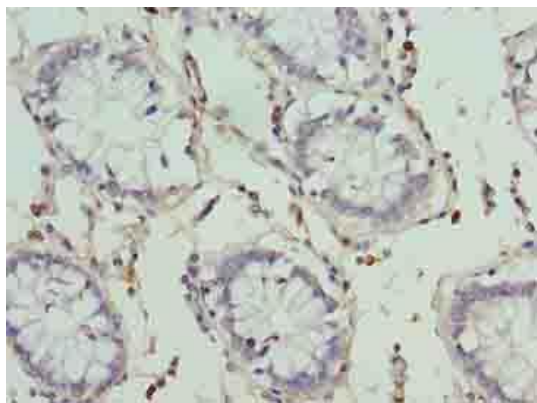
Lane 3: PC-3 whole cell lysate

Secondary

Goat polyclonal to Rabbit IgG at 1/10000 dilution

Predicted band size: 103, 91, 83 kDa

Observed band size: 103 kDa



Immunohistochemistry of paraffin-embedded human colon cancer using HKDC1 Polyclonal Antibody at dilution of 1:100