

IDH1 Polyclonal Antibody

(Catalog # A53345)

Description

IDH1 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Mouse

Isotype

IgG

Uniprot ID

O88844

Purification

Protein G purified

Immunogen

Recombinant Mouse Isocitrate dehydrogenase [NADP] cytoplasmic protein (1-240AA)

Storage

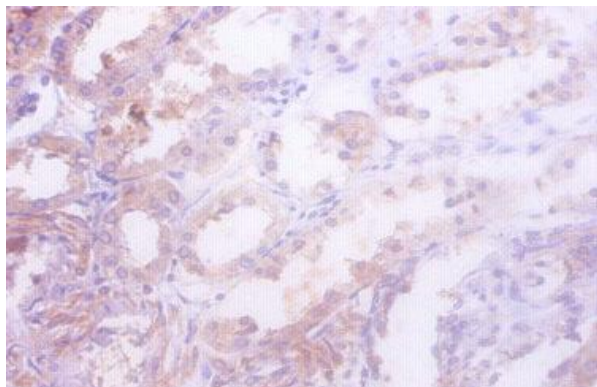
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

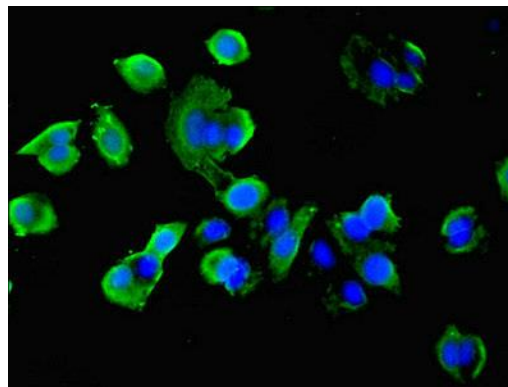
Cytosolic NADP-isocitrate dehydrogenase IDP NADP(+)-specific ICDH Oxalosuccinate decarboxylase Idh1

Application

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



Immunohistochemistry of paraffin-embedded human kidney using IDH1 Polyclonal Antibody at dilution of 1:50



Immunofluorescent analysis of MCF-7 cells using IDH1 Polyclonal Antibody at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG (H+L)