

GPI Polyclonal Antibody

(Catalog # A53853)

Background

Besides its role as a glycolytic enzyme, mammalian GPI can function as a tumor-secreted cytokine and an angiogenic factor (AMF) that stimulates endothelial cell motility. GPI is also a neurotrophic factor (Neuroleukin) for spinal and sensory neurons.

Description

GPI Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

P06744

Purification

Protein G purified

Immunogen

Recombinant Human Glucose-6-phosphate isomerase protein (2-554AA)

Storage

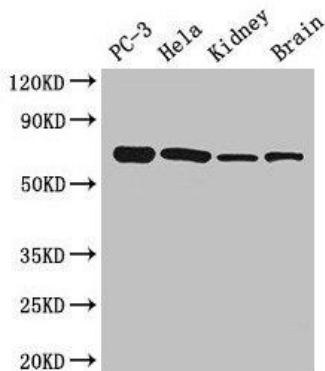
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Autocrine motility factor, Phosphoglucose isomerase, Phosphohexose isomerase, Sperm antigen 36

Application

ELISA, WB, IHC, IF, IP; Recommended dilution: WB:1:500-1:5000, IHC:1:20-1:200, IF:1:50-1:200, IP:1:200-1:2000



Western Blot

Positive WB detected in: PC-3 whole cell lysate, HeLa whole cell lysate, Mouse kidney tissue, Mouse brain tissue

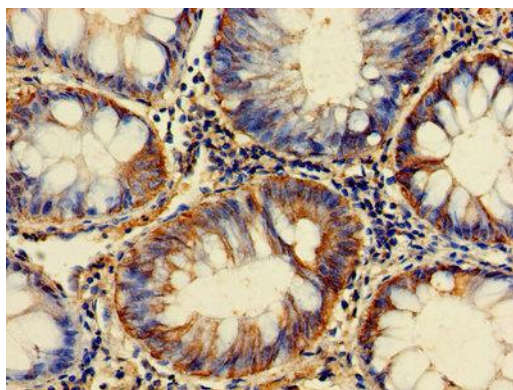
All lanes: GPI antibody at 2.7ug/ml

Secondary

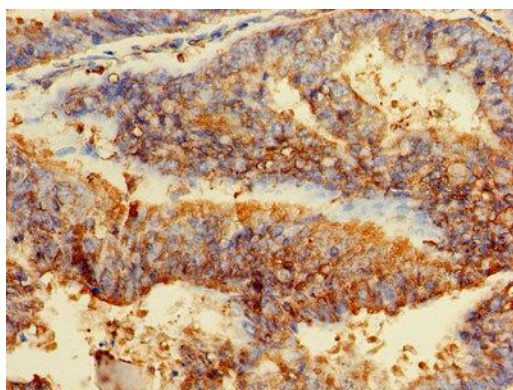
Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 64, 65 kDa

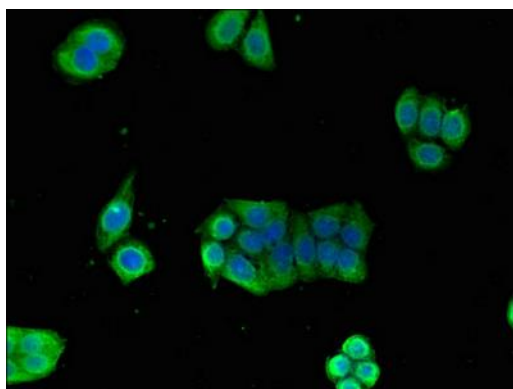
Observed band size: 64 kDa



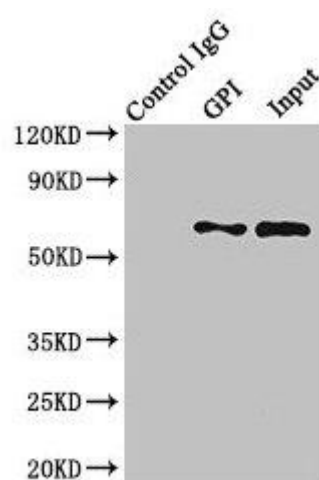
Immunohistochemistry of paraffin-embedded human colon cancer using GPI Antibody at dilution of 1:100



Immunohistochemistry of paraffin-embedded human endometrial cancer using GPI Antibody at dilution of 1:100



Immunofluorescent analysis of PC-3 cells using GPI Antibody at dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)
Immunofluorescent analysis of PC-3 cells using GPI Antibody at dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)



Immunoprecipitating GPI in HeLa whole cell lysate
Lane 1: Rabbit control IgG (1ug) instead of GPI Antibody in HeLa whole cell lysate. For western blotting, an HRP-conjugated Protein G antibody was used as the secondary antibody (1/2000)
Lane 2: GPI Antibody (8ug) + HeLa whole cell lysate (500ug)
Lane 3: HeLa whole cell lysate (10ug)