

ZNF821 Polyclonal Antibody

(Catalog # A64036)

Background

May be involved in transcriptional regulation.

Description

ZNF821 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

O75541

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Zinc finger protein 821 protein (1-370AA)

Storage

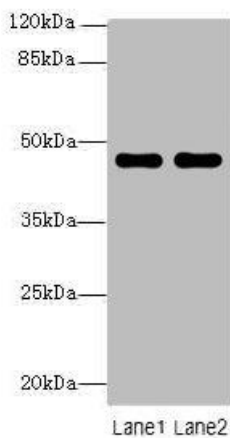
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Zinc finger protein 821, ZNF821

Application

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



Western blot

All lanes: ZNF821 Polyclonal Antibody at 8ug/ml

Lane 1: MCF7 whole cell lysate

Lane 2: HCT116 whole cell lysate

Secondary

Goat polyclonal to Rabbit IgG at 1/10000 dilution

Predicted band size: 47, 43 kDa

Observed band size: 47 kDa



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



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