

ZNF689 Polyclonal Antibody

(Catalog # A64008)

Background

May be involved in transcriptional regulation.

Description

ZNF689 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

Q96CS4

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Zinc finger protein 689 protein (1-300AA)

Storage

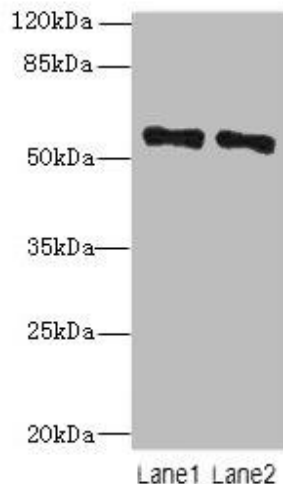
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Zinc finger protein 689, ZNF689

Application

ELISA, WB; Recommended dilution: WB:1:1000-1:5000



Western blot

All lanes: ZNF689 Polyclonal Antibody at 5ug/ml

Lane 1: Hela whole cell lysate

Lane 2: HepG-2 whole cell lysate

Secondary

Goat polyclonal to Rabbit IgG at 1/10000 dilution

Predicted band size: 57 kDa

Observed band size: 57 kDa