
ZNF410 Polyclonal Antibody

(Catalog #A63888)

Background

Transcription factor that activates transcription of matrix-remodeling genes such as MMP1 during fibroblast senescence.

Description

ZNF410 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

Q86VK4

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Zinc finger protein 410 protein (1-300AA)

Storage

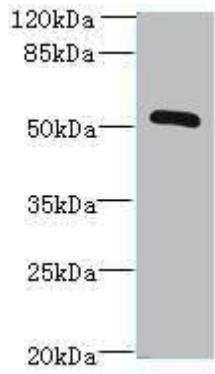
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Zinc finger protein 410, Another partner for ARF 1, Zinc finger protein APA-1, ZNF410, APA1

Application

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



Western blot

All lanes: ZNF410 Polyclonal Antibody at 3ug/ml+Mouse brain tissue

Goat polyclonal to rabbit at 1/10000 dilution

Predicted band size: 53, 30, 48, 45, 56 kDa

Observed band size: 52 kDa