

ZBTB43 Polyclonal Antibody

(Catalog # A61838)

Background

May be involved in transcriptional regulation.

Description

ZBTB43 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

O43298

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Zinc finger and BTB domain-containing protein 43 protein (165-467AA)

Storage

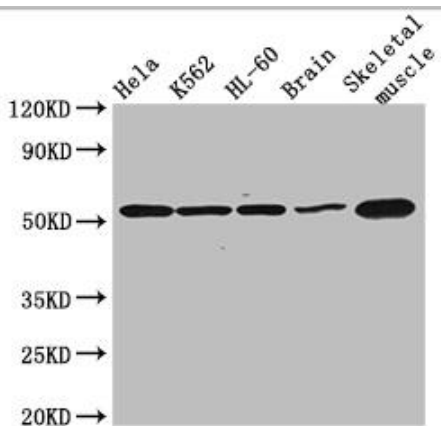
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Zinc finger and BTB domain-containing protein 43, ZBTB43, KIAA0414, ZBTB22B, ZNF297B, Zinc finger and BTB domain-containing protein 22B, Zinc finger protein 297B, ZnF-x

Application

ELISA, WB, IHC; Recommended dilution: WB:1:500-1:5000, IHC:1:20-1:200



Western Blot

Positive WB detected in: HeLa whole cell lysate, K562 whole cell lysate, HL60 whole cell lysate, Mouse brain tissue, Mouse skeletal muscle tissue

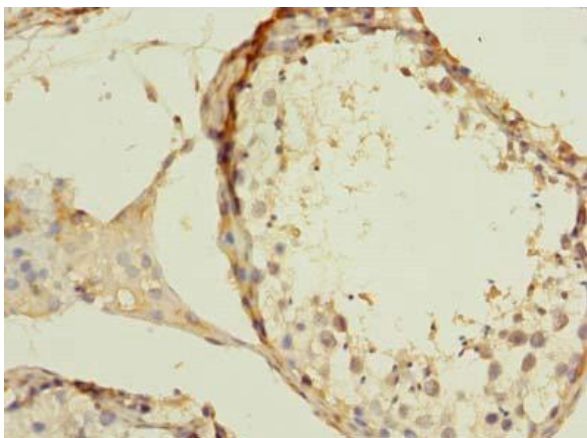
All lanes: ZBTB43 Polyclonal Antibody at 3ug/ml

Secondary

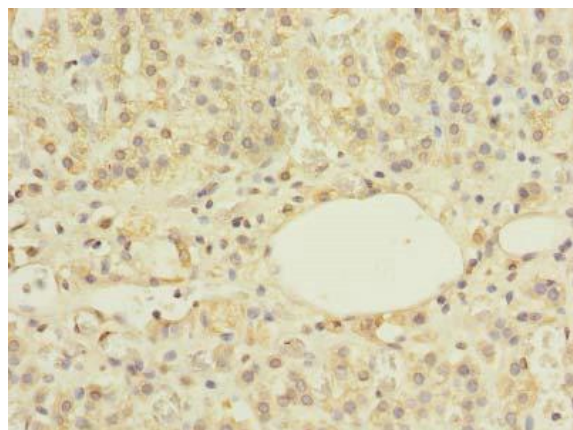
Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 53 kDa

Observed band size: 53 kDa



Immunohistochemistry of paraffin-embedded human testis at dilution of 1:100



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