

TADA2B Polyclonal Antibody

(Catalog #A65754)

Background

Coactivates PAX5-dependent transcription together with either SMARCA4 or GCN5L2.

Description

TADA2B 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

Q86TJ2

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Transcriptional adapter 2-beta protein (29-328AA)

Storage

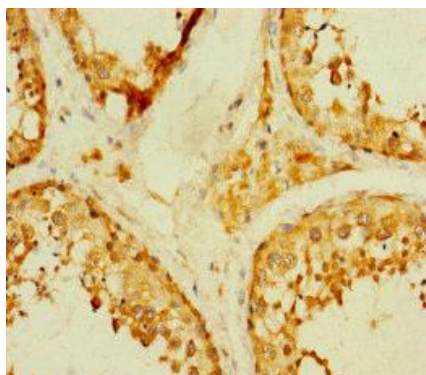
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

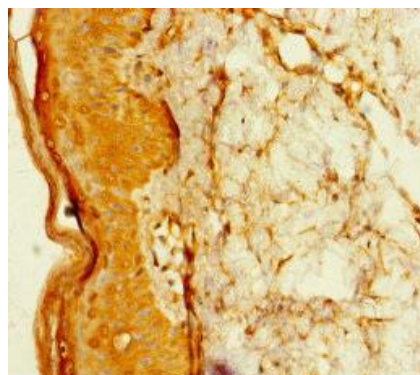
Transcriptional adapter 2-beta, ADA2-like protein beta, ADA2-beta, TADA2B, ADA2B

Application

ELISA, IHC; Recommended dilution: IHC:1:20-1:200



Immunohistochemistry of paraffin-embedded human testis tissue using TADA2B Polyclonal Antibody at dilution of 1:100



Immunohistochemistry of paraffin-embedded human skin tissue using TADA2B Polyclonal Antibody at dilution of 1:100