

ACTL7B Polyclonal Antibody

(Catalog # A68234)

Background

Actin cytoskeleton, structural constituent of cytoskeleton

Description

ACTL7B 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

Q9Y614

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Actin-like protein 7B protein (1-415AA)

Storage

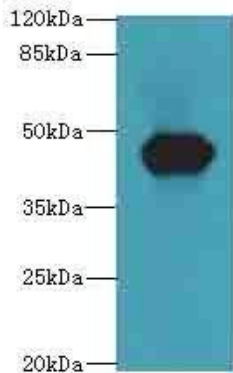
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

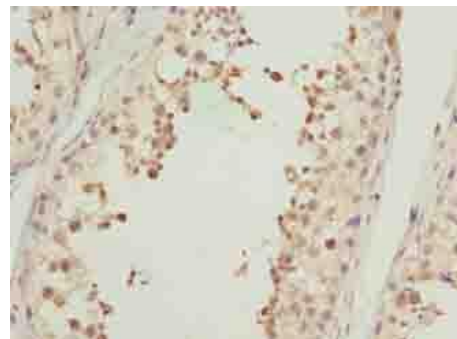
Actin-like protein 7B, Actin-like-7-beta, ACTL7B

Application

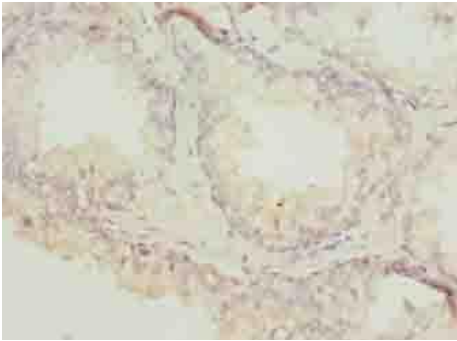
ELISA, WB, IHC; Recommended dilution: WB:1:1000-1:5000, IHC:1:20-1:200



Western blot
All lanes: ACTL7B Polyclonal Antibody at 0.2ug/ml+Mouse gonadal tissue
Goat polyclonal to rabbit at 1/10000 dilution
Predicted band size: 45 kDa
Observed band size: 45 kDa



Immunohistochemistry of paraffin-embedded human testis tissue using ACTL7B Polyclonal Antibody at dilution of 1:100



Immunohistochemistry of paraffin-embedded human prostate cancer using ACTL7B Polyclonal Antibody at dilution of 1:100