

DDX31 Polyclonal Antibody

(Catalog # A58746)

Background

Probable ATP-dependent RNA helicase. Plays a role in ribosome biogenesis and TP53/p53 regulation through its interaction with NPM1 (PubMed:23019224).

Description

DDX31 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

Q9H8H2

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Probable ATP-dependent RNA helicase DDX31 protein (560-851AA)

Storage

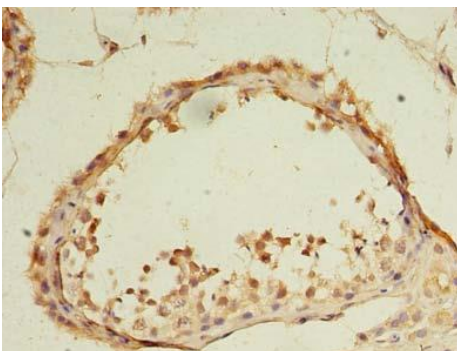
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

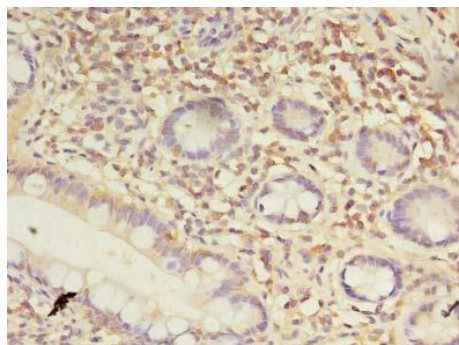
Probable ATP-dependent RNA helicase DDX31 (EC:3.6.4.13), DDX31, DEAD box protein 31, Helicain

Application

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



Immunohistochemistry of paraffin-embedded human testis tissue using DDX31 Polyclonal Antibody at dilution of 1:100



Immunohistochemistry of paraffin-embedded human small intestine tissue using DDX31 Polyclonal Antibody at dilution of 1:100