

WDR37 Polyclonal Antibody

(Catalog # A61762)

Description

WDR37 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

0.03% Proclin 300, 50% Glycerol, 0.01M PBS, PH 7.4.

Specificity

Human

Isotype

IgG

Uniprot ID

Q9Y2I8

Purification

Protein G purified

Immunogen

Recombinant Human WD repeat-containing protein 37 protein (1-264AA)

Storage

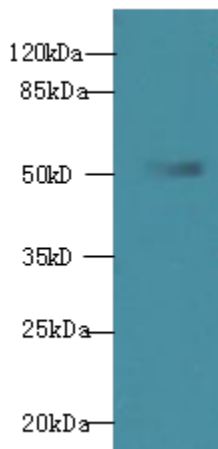
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

WD repeat-containing protein 37, WDR37, KIAA0982

Application

ELISA, WB, IHC; Recommended dilution: WB:1:500-1:2000, IHC:1:20-1:200



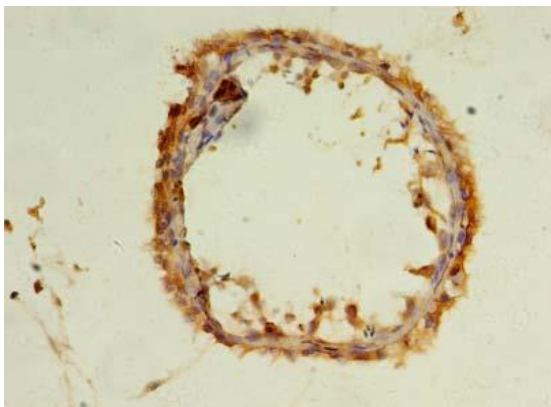
Western blot

All lanes: WDR37 Polyclonal Antibody at 7ug/ml+ HepG-2 whole cell lysate

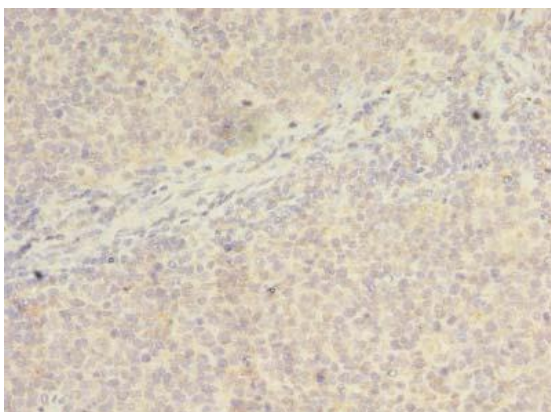
Goat polyclonal to rabbit at 1/10000 dilution

Predicted band size: 55 kDa

Observed band size: 55 kDa



Immunohistochemistry of paraffin-embedded human testis tissue using WDR37 Polyclonal Antibody at dilution of 1:100



Immunohistochemistry of paraffin-embedded human tonsil tissue using WDR37 Polyclonal Antibody at dilution of 1:100