
ARMC10 Polyclonal Antibody

(Catalog #A58146)

Background

May play a role in cell survival and cell growth. May suppress the transcriptional activity of p53/TP53.

Description

ARMC10 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

Q8N2F6

Purification

Protein G purified

Immunogen

Recombinant Human Armadillo repeat-containing protein 10 protein (28-160AA)

Storage

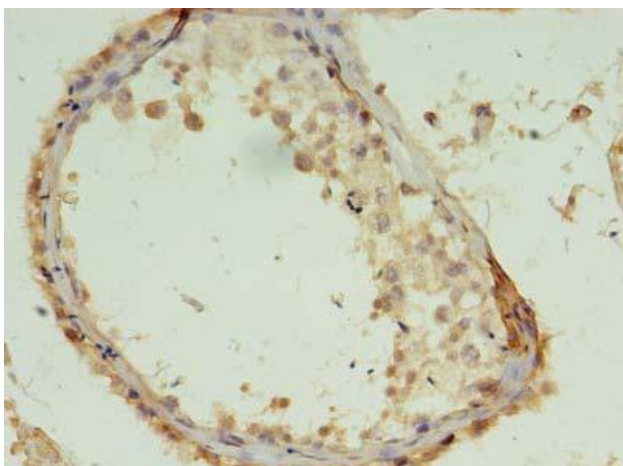
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

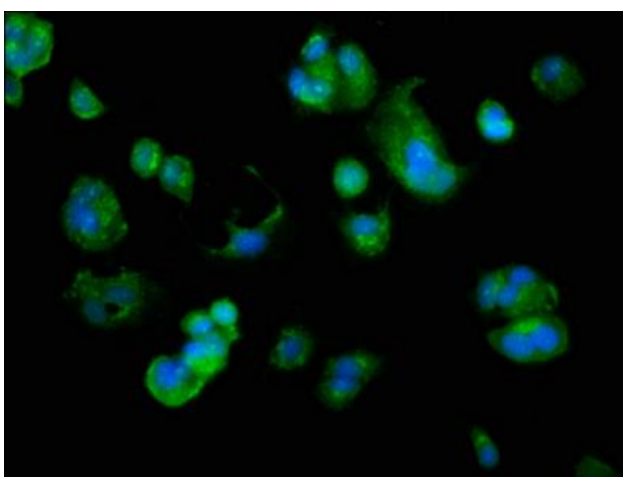
Armadillo repeat-containing protein 10, ARMC10, SVH, PSEC0198, Splicing variant involved in hepatocarcinogenesis protein

Application

ELISA, IHC, IF; Recommended dilution: IHC:1:20-1:200, IF:1:50-1:200



Immunohistochemistry of paraffin-embedded human testis tissue using ARMC10 Polyclonal Antibody at dilution of 1:100



Immunofluorescence staining of MCF-7 cells with ARMC10 Polyclonal Antibody at 1:166, counter-stained with DAPI. The cells were fixed in 4% formaldehyde, permeabilized using 0.2% Triton X-100 and blocked in 10% normal Goat Serum. The cells were then incubated with the antibody overnight at 4°C. The secondary antibody was Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L).