

Amh Polyclonal Antibody

(Catalog # A54922)

Background

This glycoprotein, produced by the Sertoli cells of the testis, causes regression of the Muellerian duct. It is also able to inhibit the growth of tumors derived from tissues of Muellerian duct origin.

Description

Amh Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

P27106

Purification

Protein G purified

Immunogen

Recombinant Mouse Muellerian-inhibiting factor protein (450-552AA)

Storage

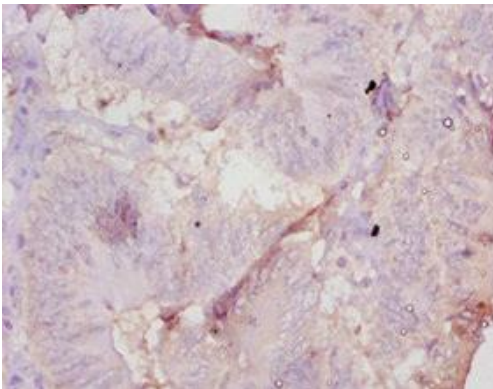
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

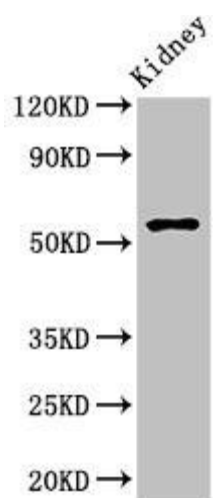
Anti-Muellerian hormone, AMH, MIS, Amh

Application

ELISA, WB, IHC; Recommended dilution: WB:1:500-1:5000, IHC:1:10-1:100



Immunohistochemistry of paraffin-embedded human colon cancer using Amh Polyclonal Antibody at dilution 1:10.



Western Blot

Positive WB detected in: Mouse kidney tissue

All lanes: Amh Polyclonal Antibody at 2.7ug/ml

Secondary

Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 60 KDa

Observed band size: 60 KDa