
AMH Polyclonal Antibody, FITC Conjugated

(Catalog # A56140)

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, FITC Conjugated. FITC. Raised in: Rabbit.

Formulation

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

Specificity

Human

Isotype

IgG

Uniprot ID

P03971

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Muellerian-inhibiting factor protein (26-560AA)

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, Muellerian-inhibiting substance, MIF, AMH

Application

ELISA; Recommended dilution: ELISA 1:100-1:500
