

XAGE1A Polyclonal Antibody

(Catalog # A61802)

Description

XAGE1A 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

Q9HD64

Purification

>95%, Protein G purified

Immunogen

Recombinant Human X antigen family member 1 protein (1-81AA)

Storage

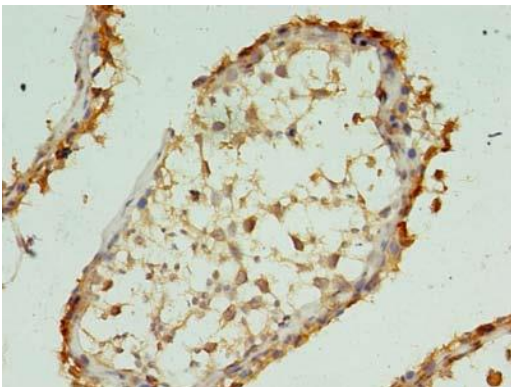
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

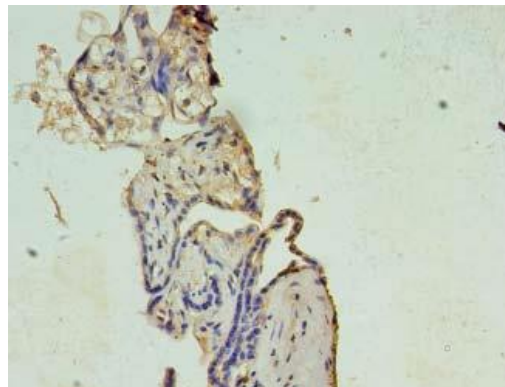
X antigen family member 1, XAGE-1, XAGE1A, GAGED2, XAGE1, AND, XAGE1B, AND, XAGE1C, AND, XAGE1D, AND, XAGE1E, Cancer/testis antigen 12.1, CT12.1, G antigen family D member 2

Application

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



Immunohistochemistry of paraffin-embedded human testis using XAGE1A Polyclonal Antibody at dilution of 1:100



Immunohistochemistry of paraffin-embedded human placenta using XAGE1A Polyclonal Antibody at dilution of 1:100