

SPACA1 Polyclonal Antibody

(Catalog # A67290)

Background

May be involved in sperm-egg fusion.

Description

SPACA1 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

Q9HBV2

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Sperm acrosome membrane-associated protein 1 protein (30-221AA)

Storage

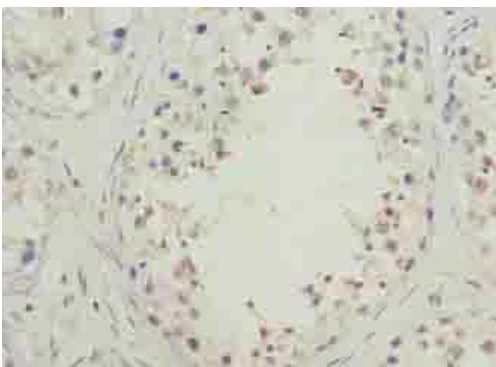
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

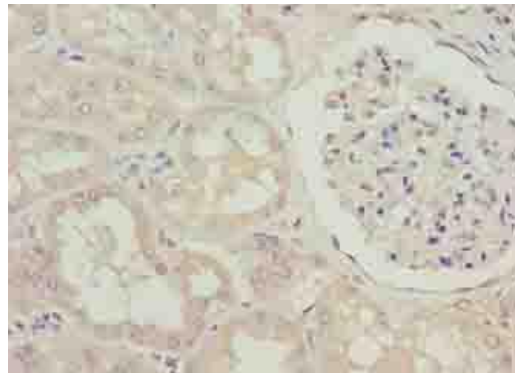
Sperm acrosome membrane-associated protein 1, Sperm acrosomal membrane-associated protein 32, SPACA1, SAMP32

Application

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



Immunohistochemistry of paraffin-embedded human testis tissue using SPACA1 Polyclonal Antibody at dilution of 1:100



Immunohistochemistry of paraffin-embedded human kidney tissue using SPACA1 Polyclonal Antibody at dilution of 1:100