
SPAM1 Polyclonal Antibody

(Catalog # A50496)

Background

Involved in sperm-egg adhesion. Upon fertilization sperm must first penetrate a layer of cumulus cells that surrounds the egg before reaching the zona pellucida. The cumulus cells are embedded in a matrix containing hyaluronic acid which is formed prior to ovulation. This protein aids in penetrating the layer of cumulus cells by digesting hyaluronic acid.

Description

SPAM1 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

P38567

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Hyaluronidase PH-20 protein (36-490AA)

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

Hyal-PH20, Hyaluronoglucosaminidase PH-20, Sperm adhesion molecule 1, Sperm surface protein PH-20, HYAL3, PH20, SPAM1

Application

ELISA; Recommended dilutions: ELISA 1:500 to 1:2000
