

SPCS1 Polyclonal Antibody

(Catalog # A61118)

Background

Component of the microsomal signal peptidase complex which removes signal peptides from nascent proteins as they are translocated into the lumen of the endoplasmic reticulum.

Description

SPCS1 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

Q9Y6A9

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Signal peptidase complex subunit 1 protein (70-102AA)

Storage

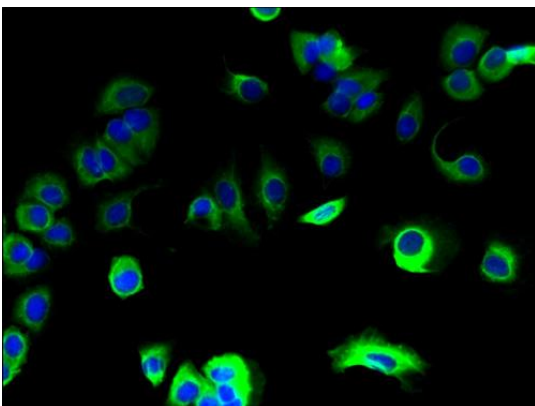
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Signal peptidase complex subunit 1 (EC:3.4.-.-), SPCS1, SPC12, HSPC033, Microsomal signal peptidase 12 kDa subunit, SPase 12 kDa subunit

Application

ELISA, IF; Recommended dilution: IF:1:50-1:200



Immunofluorescent analysis of MCF-7 cells using SPCS1 Polyclonal Antibody at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)