
SLPI Polyclonal Antibody

(Catalog #A51962)

Background

Acid-stable proteinase inhibitor with strong affinities for trypsin, chymotrypsin, elastase, and cathepsin G. May prevent elastase-mediated damage to oral and possibly other mucosal tissues.

Description

SLPI 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

P03973

Purification

>95%, Protein G purified

Immunogen

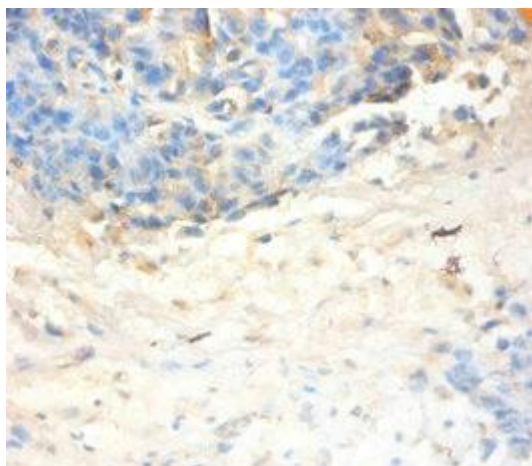
Recombinant Human Antileukoproteinase protein (26-132AA)

Storage

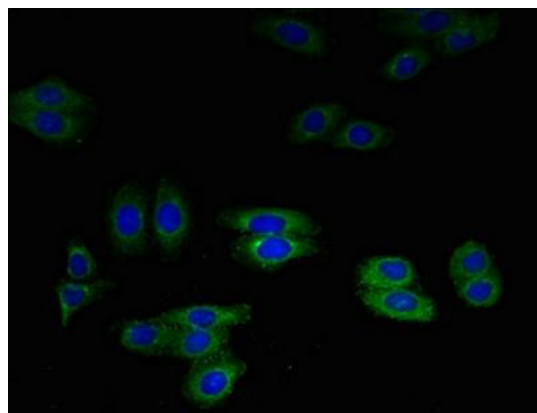
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Application

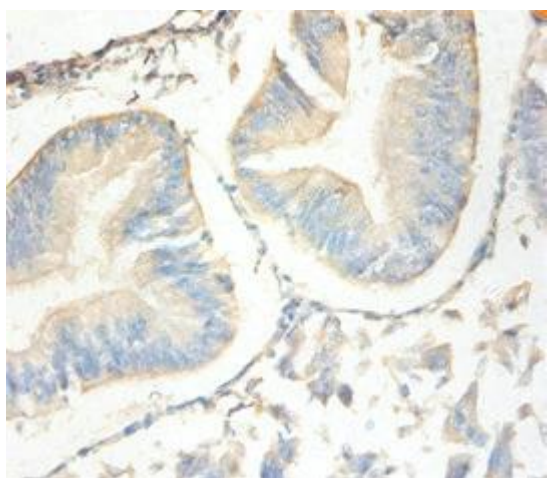
ELISA, IHC, IF; Recommended dilution: IHC:1:20-1:200, IF:1:50-1:200



Immunohistochemistry of paraffin-embedded human ovarian cancer using SLPI Antibody at dilution of 1:100



Immunofluorescent analysis of HepG2 cells using SLPI Antibody at dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)



Immunohistochemistry of paraffin-embedded human endometrial cancer using SLPI Antibody at dilution of 1:100