
SFTPC Polyclonal Antibody, FITC Conjugated

(Catalog # A60876)

Background

Pulmonary surfactant associated proteins promote alveolar stability by lowering the surface tension at the air-liquid interface in the peripheral air spaces.

Description

SFTPC Polyclonal Antibody, FITC Conjugated. FITC. Raised in: Rabbit.

Formulation

Liquid. 0.03% Proclin 300. 50% Glycerol, 0.01M PBS, pH 7.4.

Specificity

Human

Isotype

IgG

Uniprot ID

P11686

Purification

Affinity purified

Immunogen

Recombinant Human Pulmonary surfactant-associated protein C protein (24-58AA)

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

Pulmonary surfactant-associated protein C, SP-C, SFTPC, SFTP2, Pulmonary surfactant-associated proteolipid SPL(Val), SP5

Application

ELISA; Recommended dilution: ELISA 1:100-1:500
