
ALPI Polyclonal Antibody, FITC Conjugated

(Catalog # A52184)

Background

Alkaline phosphatase, intestinal also known as ALPI is a type of alkaline phosphatase that in humans is encoded by the ALPI gene. There are at least four distinct but related alkaline phosphatases: intestinal, placental, placental-like, and liver/bone/kidney (tissue non-specific). The intestinal alkaline phosphatase gene encodes a digestive brush-border enzyme. This enzyme is a component of the gut mucosal defense system and is thought to function in the detoxification of lipopolysaccharide, and in the prevention of bacterial translocation in the gut.

Description

ALPI 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

P09923

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Intestinal-type alkaline phosphatase protein (26-500AA)

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

IAP, Intestinal alkaline phosphatase, ALPI

Application

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