
VKORC1L1 Polyclonal Antibody

(Catalog #A66186)

Background

Involved in vitamin K metabolism. Can reduce inactive vitamin K 2,3-epoxide to active vitamin K (in vitro), and may contribute to vitamin K-mediated protection against oxidative stress. Plays a role in vitamin K-dependent gamma-carboxylation of Glu residues in target proteins.

Description

VKORC1L1 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

Q8N0U8

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Vitamin K epoxide reductase complex subunit 1-like protein 1 protein (38-91AA)

Storage

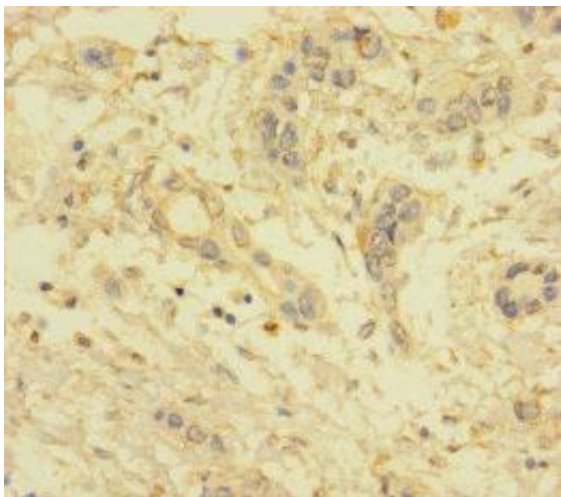
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

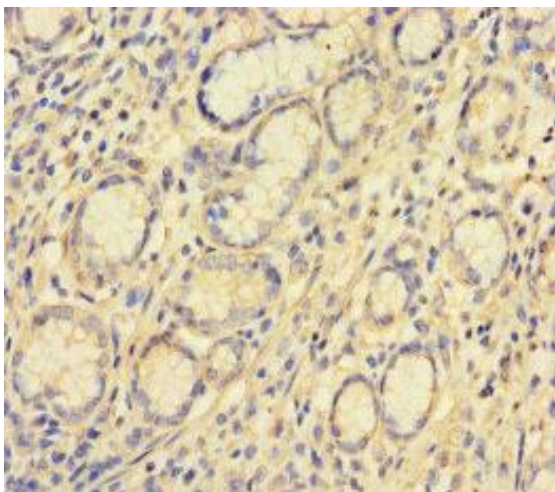
Vitamin K epoxide reductase complex subunit 1-like protein 1, VKORC1L1

Application

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



Immunohistochemistry of paraffin-embedded human pancreatic cancer using VKORC1L1 Polyclonal Antibody at dilution of 1:100



Immunohistochemistry of paraffin-embedded human stomach tissue using VKORC1L1 Polyclonal Antibody at dilution of 1:100