
ESM1 Polyclonal Antibody

(Catalog #A58962)

Background

Involved in angiogenesis; promotes angiogenic sprouting. May have potent implications in lung endothelial cell-leukocyte interactions.

Description

ESM1 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

Q9NQ30

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Endothelial cell-specific molecule 1 protein (20-184AA)

Storage

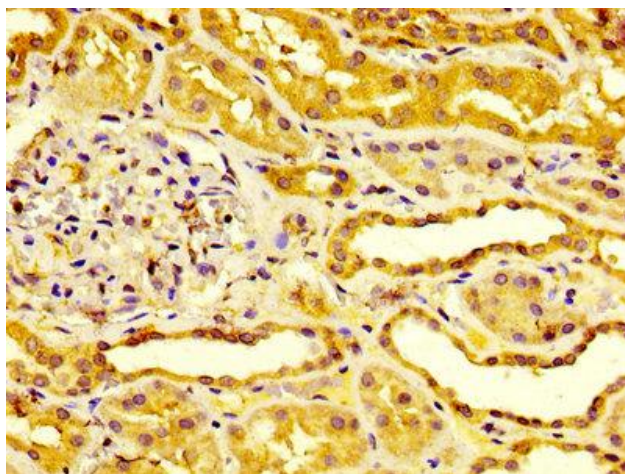
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Endothelial cell-specific molecule 1, ESM-1, ESM1

Application

ELISA, IHC; Recommended dilution: IHC:1:200-1:500



IHC image of ESM1 Polyclonal Antibody diluted at 1:400 and staining in paraffin-embedded human kidney tissue performed on a Leica Bond™ system. After dewaxing and hydration, antigen retrieval was mediated by high pressure in a citrate buffer (pH 6.0). Section was blocked with 10% normal goat serum 30min at RT. Then primary antibody (1% BSA) was incubated at 4°C overnight. The primary is detected by a biotinylated secondary antibody and visualized using an HRP conjugated SP system.