

# **EMILIN1 Polyclonal Antibody**

(Catalog # A58934)

## **Background**

May be responsible for anchoring smooth muscle cells to elastic fibers, and may be involved not only in the formation of the elastic fiber, but also in the processes that regulate vessel assembly. Has cell adhesive capacity.

## **Description**

EMILIN1 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**

Q9Y6C2

#### **Purification**

>95%, Protein G purified

#### **Immunogen**

Recombinant Human EMILIN-1 protein (1-347AA)

### Storage

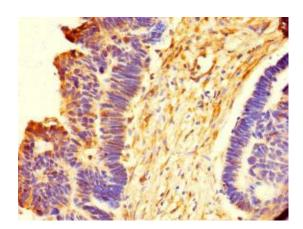
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

## **Alternative Names**

EMILIN-1, EMILIN1, EMI, Elastin microfibril interface-located protein 1, Elastin microfibril interfacer 1

#### **Application**

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



Immunohistochemistry of paraffin-embedded human ovarian tissue using EMILIN1 Polyclonal Antibody at dilution of 1:100