

---

## ASPH Polyclonal Antibody, Biotin Conjugated

(Catalog # A54362)

---

### Background

Isoform 1: specifically hydroxylates an Asp or Asn residue in certain epidermal growth factor-like (EGF) domains of a number of proteins. Ref. 13 Isoform 8: membrane-bound  $\text{Ca}^{2+}$ -sensing protein, which is a structural component of the ER-plasma membrane junctions. Isoform 8 regulates the activity of  $\text{Ca}^{+2}$  released-activated  $\text{Ca}^{+2}$  (CRAC) channels in T-cells.

### Description

ASPH Polyclonal Antibody, Biotin Conjugated. Biotin. Raised in: Rabbit.

### Formulation

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

### Specificity

Human

### Isotype

IgG

### Uniprot ID

Q12797

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Human Aspartyl/asparaginyl beta-hydroxylase protein (75-270AA)

### 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

Aspartate beta-hydroxylase Peptide-aspartate beta-dioxygenase ASPH

### Application

ELISA; Recommended dilution: ELISA 1:500-1:1000

---