

---

## TESC Polyclonal Antibody

(Catalog # A57293)

---

### Background

Functions as an integral cofactor in cell pH regulation by controlling plasma membrane-type Na<sup>+</sup>/H<sup>+</sup> exchange activity. Promotes the maturation, transport, cell surface stability and exchange activity of SLC9A1/NHE1 at the plasma membrane. Promotes the induction of hematopoietic stem cell differentiation toward megakaryocytic lineage. Essential for the coupling of ERK cascade activation with the expression of ETS family genes in megakaryocytic differentiation. Also involved in granulocytic differentiation in a ERK-dependent manner. Inhibits the phosphatase activity of calcineurin.

### Description

TESC Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

### Formulation

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

### Specificity

Human, Mouse

### Isotype

IgG

### Uniprot ID

Q96BS2

### Purification

Protein G purified

### Immunogen

Recombinant Human Calcineurin B homologous protein 3 protein (2-214AA)

### Storage

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

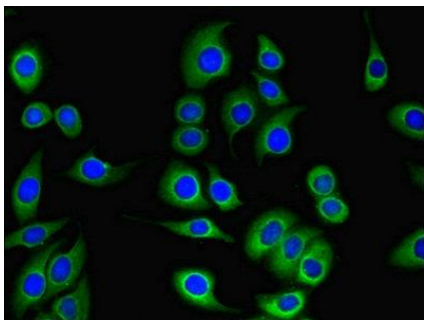
### Alternative Names

Tescalcin, CHP3, TESC

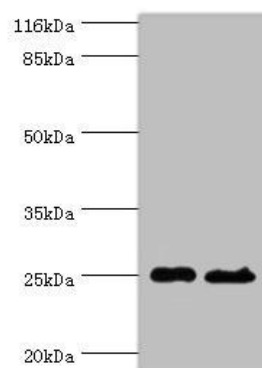
### Application

ELISA, WB, IF; Recommended dilution: WB:1:1000-1:5000, IF:1:50-1:200

---



Immunofluorescent analysis of A549 cells using Polyclonal Antibody at dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)



#### Western blot

All lanes: Polyclonal Antibody at 2 $\mu$ g/ml

Lane 1: Hela whole cell lysate

Lane 2: Mouse brain tissue

Secondary

Goat polyclonal to rabbit at 1/10000 dilution

Predicted band size: 25, 19, 22 kDa

Observed band size: 25 kDa