

---

## MAGEA9 Polyclonal Antibody

(Catalog #A59730)

---

### Background

Not known, though may play a role in embryonal development and tumor transformation or aspects of tumor progression.

### Description

MAGEA9 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

P43362

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Human Melanoma-associated antigen 9 protein (1-315AA)

### Storage

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

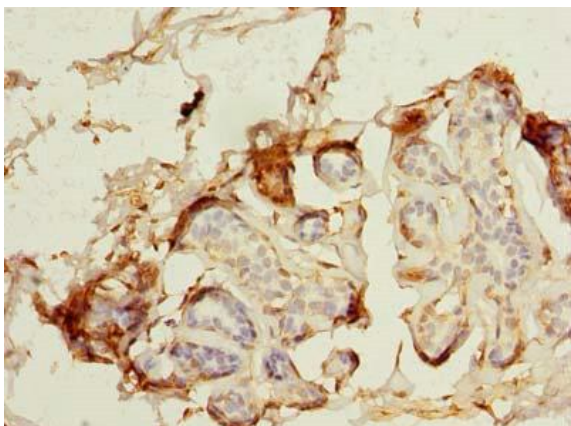
### Alternative Names

Melanoma-associated antigen 9, MAGEA9, MAGE9, MAGEA9A, AND, MAGEA9B, Cancer/testis antigen 1.9, CT1.9, MAGE-9 antigen

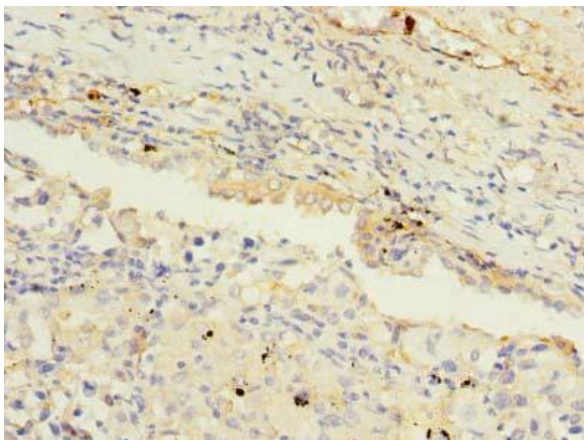
### Application

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

---



Immunohistochemistry of paraffin-embedded human breast cancer at dilution of 1:100



Immunohistochemistry of paraffin-embedded human lung cancer at dilution of 1:100