

---

## MEIS3 Polyclonal Antibody

(Catalog # A68623)

---

### Background

Transcriptional regulator which directly modulates PDPK1 expression, thus promoting survival of pancreatic beta-cells. Also regulates expression of NDFIP1, BNIP3, and CCNG1.

### Description

MEIS3 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

Q99687

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Human Homeobox protein Meis3 protein (178-264AA)

### Storage

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

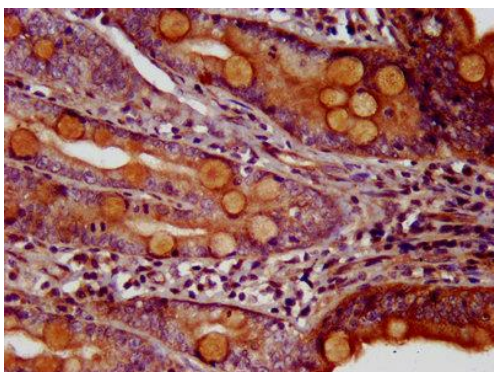
### Alternative Names

Homeobox protein Meis3, Meis1-related protein 2, MEIS3, MRG2

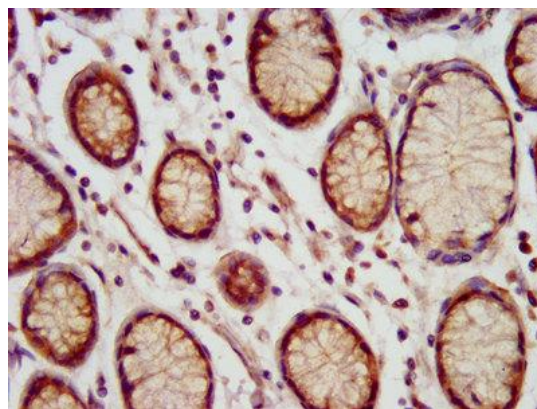
### Application

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

---



IHC image of MEIS3 Polyclonal Antibody diluted at 1:200 and staining in paraffin-embedded human small intestine 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.



IHC image of MEIS3 Polyclonal Antibody diluted at 1:200 and staining in paraffin-embedded human gastric cancer 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.