

---

## RSPO3 Polyclonal Antibody

(Catalog #A60722)

---

### Background

Activator of the canonical Wnt signaling pathway by acting as a ligand for LGR4-6 receptors. Upon binding to LGR4-6 (LGR4, LGR5 or LGR6), LGR4-6 associate with phosphorylated LRP6 and frizzled receptors that are activated by extracellular Wnt receptors, triggering the canonical Wnt signaling pathway to increase expression of target genes. Also regulates the canonical Wnt/beta-catenin-dependent pathway and non-canonical Wnt signaling by acting as an inhibitor of ZNRF3, an important regulator of the Wnt signaling pathway. Acts as a ligand for frizzled FZD8 and LRP6. May negatively regulate the TGF-beta pathway.

### Description

RSPO3 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

### Formulation

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

### Specificity

Human

### Isotype

IgG

### Uniprot ID

Q9BXY4

### Purification

Protein G purified

### Immunogen

Recombinant Human R-spondin-3 protein (22-272AA)

### Storage

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

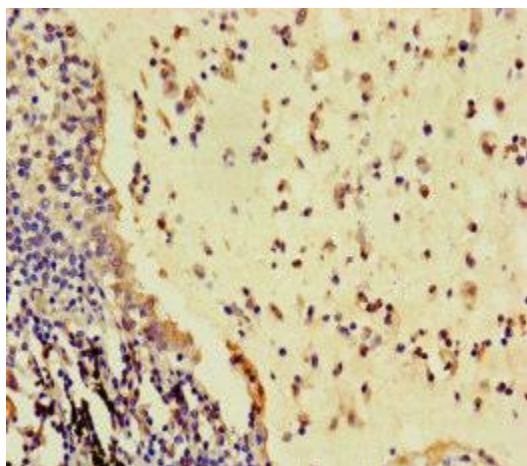
### Alternative Names

R-spondin-3, RSPO3, PWTSR, THSD2, Protein with TSP type-1 repeat, hPWTSR, Roof plate-specific spondin-3, hRspo3, Thrombospondin type-1 domain-containing protein 2

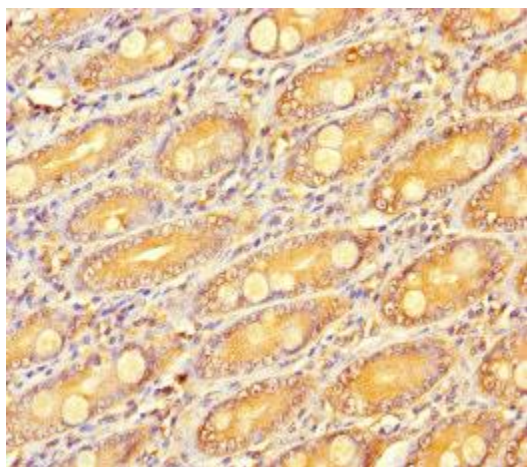
### Application

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

---



Immunohistochemistry of paraffin-embedded human lung cancer using RSPO3 Antibody at dilution of 1:100



Immunohistochemistry of paraffin-embedded human small intestine tissue using RSPO3 Antibody at dilution of 1:100