

## GBX2 Polyclonal Antibody

(Catalog # A69507)

### Background

May act as a transcription factor for cell pluripotency and differentiation in the embryo.

### Description

GBX2 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

P52951

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Human Homeobox protein GBX-2 protein (65-248AA)

### Storage

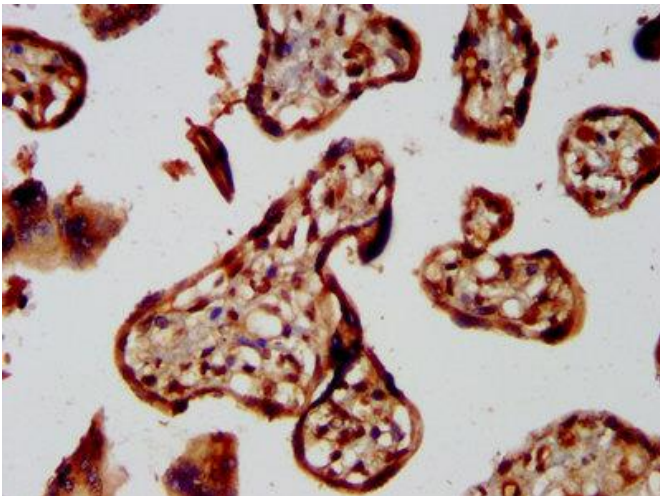
Shipped at 4°C. Upon receipt, store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

### Alternative Names

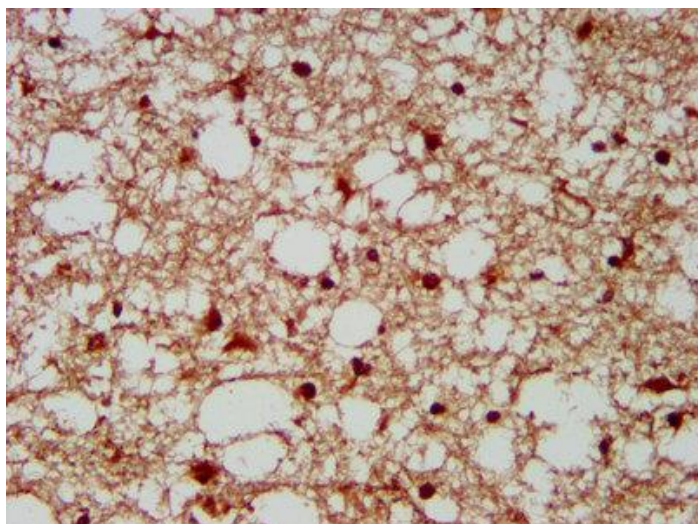
Homeobox protein GBX-2, Gastrulation and brain-specific homeobox protein 2, GBX2

### Application

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



IHC image of GBX2 Polyclonal Antibody diluted at 1:300 and staining in paraffin-embedded human placenta 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 GBX2 Polyclonal Antibody diluted at 1:300 and staining in paraffin-embedded human brain 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.