

## GLCE Polyclonal Antibody

(Catalog # A55279)

---

### Background

Converts D-glucuronic acid residues adjacent to N-sulfate sugar residues to L-iduronic acid residues, both in maturing heparan sulfate (HS) and heparin chains. This is important for further modifications that determine the specificity of interactions between these glycosaminoglycans and proteins.

### Description

GLCE Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

### Formulation

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

### Specificity

Human, Mouse, Rat

### Isotype

IgG

### Uniprot ID

O94923

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Human D-glucuronyl C5-epimerase protein (29-617AA)

### Storage

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

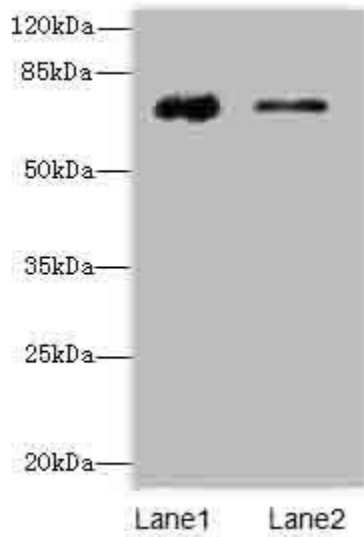
### Alternative Names

Heparan sulfate C5-epimerase, Hsepi, Heparin/heparan sulfate:glucuronic acid C5-epimerase, Heparosan-N-sulfate-glucuronate 5-epimerase, GLCE, KIAA0836

### Application

ELISA, WB, IHC; Recommended dilution: WB:1:1000-1:5000, IHC:1:20-1:200

---



Western blot

All lanes: GLCE Polyclonal Antibody at 12ug/ml

Lane 1: Rat liver tissue

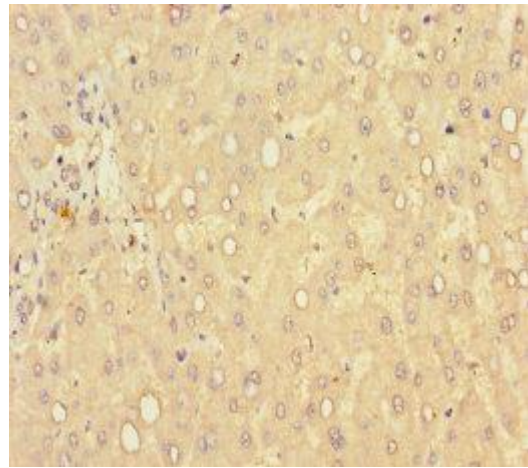
Lane 2: Mouse kidney tissue

Secondary

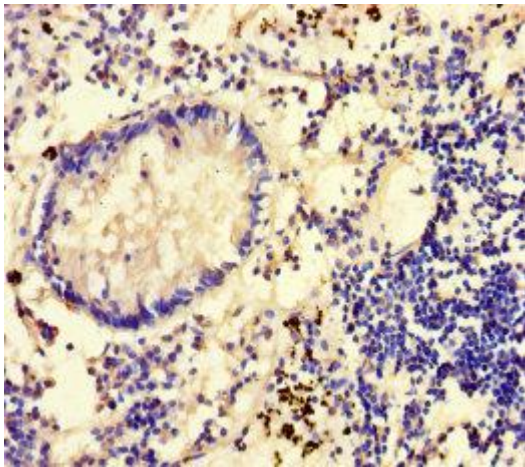
Goat polyclonal to Rabbit IgG at 1/10000 dilution

Predicted band size: 70 kDa

Observed band size: 70 kDa



Immunohistochemistry of paraffin-embedded human liver tissue at dilution of 1:100



Immunohistochemistry of paraffin-embedded human appendix tissue at dilution of 1:100