

---

## ANKRD49 Polyclonal Antibody

(Catalog # A66554)

---

### Background

Induces HBG1 expression (PubMed:16131492, PubMed:11162141). May have a role in spermatogenesis where it promotes autophagy in response to serum starvation, via the NF-kappaB pathway.

### Description

ANKRD49 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

Q8WVL7

### Purification

Protein G purified

### Immunogen

Recombinant Human Ankyrin repeat domain-containing protein 49 protein (4-239AA)

### Storage

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

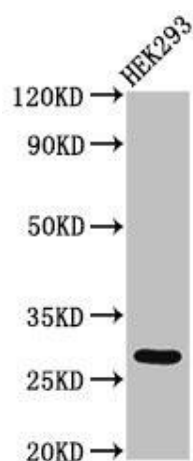
### Alternative Names

Ankyrin repeat domain-containing protein 49, Fetal globin-inducing factor, ANKRD49, FGIF

### Application

ELISA, WB, IHC; Recommended dilution: WB 1:500 - 1:5000, IHC 1:20 - 1:200

---



#### Western Blot

Positive WB detected in: HEK293 whole cell lysate

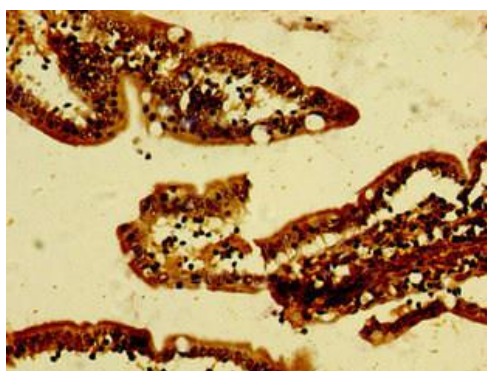
All lanes: ANKRD49 Polyclonal Antibody at 3ug/ml

Secondary

Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 28 KDa

Observed band size: 28 KDa



Immunohistochemistry of paraffin-embedded human small intestine tissue using ANKRD49 Polyclonal Antibody at dilution of 1:100