

---

## MUC16 Polyclonal Antibody, Biotin Conjugated

(Catalog # A54128)

---

### Description

MUC16 Polyclonal Antibody, Biotin Conjugated. Biotin. Raised in: Rabbit.

### Formulation

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

### Specificity

Human

### Isotype

IgG

### Uniprot ID

Q8WXI7

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Human Mucin-16 protein (21712-21973AA)

### 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

MUC-16, Ovarian cancer-related tumor marker CA125, MUC16, CA125

### Application

ELISA; Recommended dilution: ELISA 1:500-1:1000

---