

---

## AGA2 Polyclonal Antibody, Biotin Conjugated

(Catalog # A50795)

---

### Background

Receptor binding subunit of the a-agglutinin heterodimer. *S. cerevisiae* a and alpha cells express the complementary cell surface glycoproteins a-agglutinin and alpha-agglutinin, respectively, which interact with one another to promote cellular aggregation during mating.

### Description

AGA2 Polyclonal Antibody, Biotin Conjugated. Biotin. Raised in: Rabbit.

### Formulation

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

### Specificity

*Saccharomyces cerevisiae*

### Isotype

IgG

### Uniprot ID

P32781

### Purification

>95%, Protein G purified

### Immunogen

Recombinant *Saccharomyces cerevisiae* A-agglutinin-binding subunit protein (19-87AA)

### 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

YGL032C, A-agglutinin-binding subunit, AGA2

### Application

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

---