

---

## CHRNA1 Polyclonal Antibody, Biotin Conjugated

(Catalog # A53583)

---

### Background

After binding acetylcholine, the AChR responds by an extensive change in conformation that affects all subunits and leads to opening of an ion-conducting channel across the plasma membrane.

### Description

CHRNA1 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

P02708

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Human Acetylcholine receptor subunit alpha protein (21-255AA)

### 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

CHRNA1 ACHRA, CHNRA

### Application

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

---