
Allergen Ara h 1 Polyclonal Antibody

(Catalog # A64330)

Description

Allergen Ara h 1 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Peanut

Isotype

IgG

Uniprot ID

P43237

Purification

>95%, Protein G purified

Immunogen

Recombinant Peanut Allergen Ara h 1, clone P17 protein (26-216AA)

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

Allergen Ara h 1, clone P17

Application

ELISA
