

# Goat F(ab)'<sub>2</sub> Anti-Human IgA

α chain specific- Capture antibody

Catalog # fab-iga

For research use only, not for diagnostic or therapeutic use

Version # 08D25-MT

## PRODUCT INFORMATION

### Content

0.5 mg of goat F(ab)'<sub>2</sub> anti-human IgA in 1.0 ml of 100 mM borate buffered saline, pH 8.0. No preservative or amine-containing buffer salts added.

### Storage

Product is shipped at room temperature. Store at 2-8°C.

## DESCRIPTION

### Source

Pepsin digest of goat anti-human IgA

### Cross Absorption

Human IgM, IgG and IgD

### Purification

Gel filtration chromatography of pepsin digested antibody

### Specificity

Reacts with the heavy chain of human IgA as demonstrated by ELISA and flow cytometry. May also react with IgA from other species.

## CHARACTERIZATION

To insure lot-to-lot consistency, each batch of product is tested by ELISA and flow cytometry for conformance to characteristics of a standard reference reagent.

## APPLICATIONS

- Flow cytometry
- Direct immunofluorescent staining of IgA<sup>+</sup> human B lymphocytes
- Enzyme-Linked-Immunosorbent-Assay (ELISA)
- Western blotting
- Dot- and slot-immunoblotting
- Immunohistochemistry

## WORKING DILUTIONS

Optimal dilutions for use of this product vary according to specific applications and should be determined by the individual investigator.

---

### TECHNICAL SUPPORT

Toll free (US): 888-457-5873

Outside US: (+1) 858-457-5873

Europe: +33 562-71-69-39

E-mail: [info@invivogen.com](mailto:info@invivogen.com)

Website: [www.invivogen.com](http://www.invivogen.com)



3950 Sorrento Valley Blvd. Suite 100  
San Diego, CA 92121 - USA