

Goat Anti-Human IgA - HRP

α chain specific

Catalog # hrp-iga

For research use only, not for diagnostic or therapeutic use

Version # 08D25-MT

PRODUCT INFORMATION

Content

1 ml of goat anti-human IgA - HRP (horseradish peroxidase conjugate) in 50% glycerol/50% PBS, pH 7.4. No preservative added.

Storage

Product is shipped at room temperature. Store at 2-8°C or long-term at -20°C.

DESCRIPTION

Source

Isolated from antisera of goats hyperimmunized with human IgA paraproteins

Cross Absorption

Human IgM, IgD and IgG

Purification

Affinity chromatography on pooled human IgA covalently linked to agarose

Specificity

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

CHARACTERIZATION

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

APPLICATIONS

- Enzyme-Linked-Immunosorbent-Assay (ELISA)
- Western blotting
- Dot- and slot-immunoblotting

WORKING DILUTIONS

ELISA: 1:4,000 - 1:8,000

Other applications: Since applications vary, you should determine the optimum working dilution of the product that is appropriate for your specific need.

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
Website: www.invivogen.com

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