

Goat F(ab)'₂ Anti-Human Kappa

(κ chain specific)

Catalog # fab-igak

For research use only, not for diagnostic or therapeutic use

Version # 08D25-MT

PRODUCT INFORMATION

Content

0.5 mg of goat F(ab)'₂ anti-human Kappa 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

Antisera of goats hyperimmunized with human myeloma proteins containing κ light chains

Cross Absorption

Pooled human myeloma proteins with λ light chains

Purification

Affinity chromatography on pooled human Igs with κ light chains covalently linked to agarose

Specificity

Reacts with human κ light chains 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/or flow cytometry for conformance to characteristics of a standard reference reagent.

APPLICATIONS

- Flow cytometry
- Direct immunofluorescent staining of IgA⁺ human B lymphocytes
- Enzyme-Linked-Immunesorbent-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

Website: www.invivogen.com



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