

Human IgA

IgA standard

Catalog # ctrl-iga

For research use only, not for diagnostic or therapeutic use

Version #10A13-MM

PRODUCT INFORMATION

Content

0.5 mg of human IgA, provided as an essentially salt-free, lyophilized powder. No preservative or amine-containing buffer salts added.

Storage

Product is shipped at room temperature. Store lyophilized powder at 2-8°C. Product should be reconstituted to a minimum of 1 mg/ml with saline (150mM NaCl) solution. After reconstitution, aliquot and store at -20°C. Avoid repeated freeze-thaw cycles.

DESCRIPTION

Source

Pooled normal human colostrum

Isotype

IgA

Purification

Ion exchange chromatography

Purity

>95% as judged by SDS-polyacrylamide gel electrophoresis

CHARACTERIZATION

To ensure lot-to-lot consistency, each batch of product is tested by ELISA and immunoelectrophoresis to conform with characteristics of a standard reference reagent. In accordance with current Good Manufacturing and Good Laboratory Practices (cGMP/cGLP), any protein of human origin should be handled pursuant to your organization's documented safety procedures and as if it is capable of transmitting infection. This product has NOT been tested for viral, bacterial, or other infectious agents such as, but not limited to, HIV, HbsAg, HCV.

APPLICATIONS

- Enzyme-Linked-Immunosorbent-Assay (ELISA)
- SDS-PAGE and immunoblotting

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