



6042 Cornerstone Court West, Suite B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Data Sheet

Anti-Human IgG, Unconjugated Antibody

Mouse, Anti-human, Unconjugated,
Monoclonal
Catalog #: 100736
Lot #: 200427

Host Species | Isotype:

Mouse | IgG

Formulated in: 10 mM Tris-HCl, pH 8.0, 50 mM NaCl, 0.05% (w/v) sodium azide.

Stability: At least 1 year at 4°C. *Avoid freeze/thaw cycles.*

References:

1. Woo, P.C.Y., *et al.*, *Clinical and Vaccine Immunology*. 2004; **11(4)**: 665-668.
2. Hsueh, P.-R. *et al.*, *Clinical Microbiology and Infection* 2004; **10(12)**: 1062-1066

Description: Anti-human, unconjugated antibody recognizing human IgG generated in mouse.

Immunogen: This antibody was produced against human recombinant heavy chain IgG.

Purification: Affinity chromatography.

Applications: Useful in ELISA and lateral flow assays.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com