



RayBiotech, Inc.

3607 Parkway Lane suite 200
Norcross, GA 30092
Tel: 770-729-2992, 1-888-494-8555
Fax: 770-206-2393
Website: www.raybiotech.com
Email: info@raybiotech.com

Certificate of Analysis and Data Sheet

Mouse Anti Bovine Interleukin-4:Biotin

With HRP-conjugated Secondary Antibody

Catalog No.
DS-MB-01657

Species
Bovine

Isotype
IgG2b

Preparation

Preparation: Purified IgG prepared by affinity chromatography on Protein G from tissue culture supernatant

Immunogen: Recombinant bovine IL-4

Fusion Partners: Spleen cells from immunized BALB/c mice were fused with cells of the SP2/O myeloma cell line.

Specificity

DS-MB-01657 is specific for bovine interleukin 4 (IL-4).

IL-4 is expressed by a range of cells including T and B cells, macrophages and monocytes. It plays an important role in the regulation of bovine T and B cell responses.

Formulation

Product Type: Monoclonal Antibody

Product Form: Purified IgG conjugated to Biotin - liquid

Preservatives & Stabilizers: 0.09% Sodium Azide, 1% Bovine Serum Albumin

Buffer Solution: Phosphate buffered saline pH7.4

Approx. Protein Concentrations: IgG concentration 1.0 mg/ml

Storage

Store at -20°C only. This product should be stored undiluted. Storage in frost-free freezers is not recommended. Avoid repeated freezing and thawing as this may denature the antibody. Should this product contain a precipitate we recommend microcentrifugation before use.

**The products are furnished for LABORATORY RESEARCH USE ONLY.
Not for diagnostic or therapeutic use.**



RayBiotech, Inc.

3607 Parkway Lane suite 200
Norcross, GA 30092
Tel: 770-729-2992, 1-888-494-8555
Fax: 770-206-2393
Website: www.raybiotech.com
Email: info@raybiotech.com

Applications

Table Summary of antibody applications and working conditions

Options / Functions	YES	NO	Not determined	Recommended Work dilution or concentration
ELISA	•			1ug/ml – 5ug/ml
Western Blotting			•	
Functional Assays			•	

Note: Other applications are not tested yet. Optimal dilutions should be determined by each laboratory for each application.

Secondary Antibody Applications

Options / Functions	YES	NO	Not determined	Recommended Work dilution or concentration
Immunoassay (ELISA, Western blot)	•			1:5000-1:10000

**The products are furnished for LABORATORY RESEARCH USE ONLY.
Not for diagnostic or therapeutic use.**