



## 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](http://www.raybiotech.com)  
Email: [info@raybiotech.com](mailto:info@raybiotech.com)

# Certificate of Analysis and Data Sheet

## MOUSE ANTI BOVINE IL-10

### With HRP-conjugated Secondary Antibody

**Catalog No.**  
DS-MB-02649

**Target Species**  
Bovine

**Isotype**  
IgG2b

### **Preparation**

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

**Immunogen:** Plasmid cDNA encoding bovine IL-10.

**Fusion Partners:** Spleen cells from immunized Balb/c mice were fused with cells of the mouse SP2/0 myeloma cell line.

### **Specificity**

DS-MB-02649 recognizes bovine IL-10. This antibody has been shown not to inhibit the biological activity of IL-10.

### **Formulation**

**Product Type:** Monoclonal Antibody

**Product Form:** Purified IgG - liquid

**Buffer Solution:** Phosphate buffered saline

**Preservative Stabilizers:** 0.09% Sodium Azide (NaN<sub>3</sub>)

**Approx. Protein Concentrations:** IgG concentration 1.0 mg/ml

### **Storage**

Store at +4°C or at -20°C if preferred.

Storage in frost-free freezers is not recommended.

This product should be stored undiluted. Avoid repeated freezing and thawing as this may denature the antibody. Should this product contain a precipitate we recommend microcentrifugation before use.

**Shelf Life:** 18 months from date of dispatch.

### **Reference**

1. Kwong, L. S. et al. (2002) Development of an ELISA for bovine IL-10. Vet. Immunol. Immunopathol. 85: 213 - 223.

**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](http://www.raybiotech.com)  
Email: [info@raybiotech.com](mailto:info@raybiotech.com)

### Applications

<b>Options</b> <b>Functions</b>	<b>YES</b>	<b>NO</b>	<b>Not determined</b>	<b>Recommended Work dilution or concentration</b>
Elisa (1)	•			5 ug/ml – 10 ug/ml
Western Blotting			•	
Immunoprecipitation			•	
Flow cytometry			•	

Note: Other applications are not tested yet. Optimal dilutions should be determined by each laboratory for each application.  
(1) This reagent may be used as a capture antibody in a sandwich ELISA assay for bovine IL-10 in combination with DS-MB-02650 as detection reagent.

### Secondary Antibody Applications

<b>Options</b> <b>Functions</b>	<b>YES</b>	<b>NO</b>	<b>Not determined</b>	<b>Recommended Work dilution or concentration</b>
Immunoassay (ELISA, Western blot)	•			1:5,000-1:10,000

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