



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 Chicken Interleukin-2:Azide Free

With HRP-conjugated Secondary Antibody

Catalog No. DS-MB-01639	Target Species Chicken	Isotype IgG2b
-----------------------------------	----------------------------------	-------------------------

Preparation

- **Immunogen:** Recombinant chicken IL-2
- **Purification:** Purified IgG prepared by affinity chromatography on Protein G from tissue culture supernatant
- **Fusion Partners:** Spleen cells from immunized mice were fused with cells of the mouse Sp2/0-Ag myeloma cell line.

Formulation

- **Product Type:** Monoclonal Antibody
- **Product Form:** Purified IgG - liquid
- **Buffer:** Phosphate buffered saline pH 7.4
- **Preservative Stabilizers:** None present
- **Approx. Protein Concentration:** IgG concentration 1.0 mg/ml

Specificity

- DS-MB-01639 recognizes natural chicken IL-2 and recombinant chicken IL-2 derived from CHO cells.
- DS-MB-01639 neutralizes the proliferative activity of rchIL-2 and recognizes a different epitope to the Raybiotech, Inc product: DS-MB-01640
- **Species Cross Reactivity:** Reacts with: Mammals
- Antibody reactivity and working conditions may vary between species.

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.
- **Shelf Life:** 18 months from date of shipment

**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	•			5 µg/ml – 10 µg/ml
Western Blotting	•			Non-reducing conditions
Immunohistochemistry			•	
Functional Assays	•			2 µg/ml – 20 µg/ml
Flow Cytometry			•	

Note: Other applications are not tested yet. Optimal dilutions should be determined by the user. Membrane permeabilisation is required for this application.

Secondary Antibody Applications

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

References

1. Rothwell, L., Hamblin, A. and Kaiser, P. (2001) Production and characterisation of monoclonal antibodies specific for chicken interleukin-2. *Vet. Immunol. Immunopathol.* 83: 149-160.

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