



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 Granzyme B:Biotin

With HRP-conjugated Secondary Antibody

Catalog No.
DS-MB-01405

Target Species
Human

Isotype
IgG1

Introduction

Synonyms: CGL1, CSPB, CTLA1, GRB

Preparation

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

Immunogen: Purified human Granzyme B

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

Formulation

Product Type: Monoclonal Antibody

Product Form: Ig fraction - liquid

Buffer Solution: Phosphate buffered saline

Preservative Stabilizers: 0.01% Thiomersal

Approx. Protein Concentrations: IgG concentration 0.5 mg/ml

Specificity

DS-MB-01405 recognizes the serine protease Granzyme B, which is expressed within the granules of cytolytic lymphocytes (CTLs).

Granzyme B plays a key role in the induction of apoptosis by CTLs. After delivery to the target cell Granzyme B activates the cascade of caspases that finally results in cell death.

This antibody has been reported to function as a detection reagent in ELISA assays for soluble Granzyme B in conjunction with DS-MB-01406 as capture reagent.

Species Crossreactivity: Chimpanzee, Monkey and Rhesus Monkey. Please note, antibody reactivity and working conditions may vary between species.

**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

Storage

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

Applications

Table Summary of antibody applications and working conditions

Options / Functions	YES	NO	Not determined	Recommended Work dilution or concentration
ELISA (1)	•			1ug/ml
Western Blotting			•	
IHC – Paraffin			•	
IHC - Frozen			•	
Flow Cytometry			•	
Immunoprecipitation			•	
Functional Assays			•	

Note: Other applications are not tested yet. Optimal dilutions should be determined by each laboratory for each application.
(1) This antibody has been reported to function as a detection reagent in ELISA assays for soluble Granzyme B in conjunction with DS-MB-01406 as capture reagent.

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.**



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

References

1. Spaeny-Dekking, E. H. A. *et al.* (1998) Extracellular Granzymes A and B in humans: Detection of native species during CTL responses *in vitro* and *in vivo*. *J. Immunol.* 160: 3610 - 3616.

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