



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 HUMAN MMP-2

With HRP-conjugated Secondary Antibody

Catalog No.	Species	Isotype
DS-MB-01925	Human	Mouse IgG ₁

Preparation

Preparation: Purified IgG prepared by affinity chromatography
Immunogen: Synthetic peptide APSPIIKFPGDVAPKTDK corresponding to gelatinase
Fusion Partners: Spleen cells from immunized Balb/c mice were fused with cells of the P3X63Ag8.653 myeloma cell line.

Specificity

This product recognizes human matrix metalloproteinase-2 (MMP-2), also known as gelatinase A.

Formulation

Product Type: Monoclonal Antibody
Product Form: Purified IgG - liquid
Buffer Solution: 0.1M glycine HCl, 1M TRIS HCl, pH 7-8
Preservative & Stabilizers: 0.09% Sodium Azide
Approx. Protein Concentrations: IgG concentration 1.0mg/ml

Storage

18 months from date of dispatch.
Store at 4°C or at -20°C if preferred.
This product should be stored undiluted. Storage in frost free freezers is not recommended. Should this product contain a precipitate we recommend microcentrifugation before use.
Avoid repeated freezing and thawing as this may denature the antibody.

**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			•	
Western Blotting	•			
Immunohistology – frozen(*)	•			1:100 – 1:1000
Immunohistology - paraffin			•	
Immunohistology - resin			•	
Immunoprecipitation			•	
Flow Cytometry			•	
Immunodiffusion			•	
Neutralization			•	

Note: Other applications are not tested yet. Optimal dilutions should be determined by each laboratory for each application.
 (*) Histology Positive Control Tissue Trophoblast

Secondary Antibody Applications

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

Reference

1. Soini, Y. et al. (1994) 72kD and 92kD type IV collagenase, type IV collagen, and laminin mRNAs in breast cancer: a study by in situ hybridization. J. Histochem. Cytochem. 42: 945-951.
2. Autio Harmainen, H. et al. (1993) Expression of 72 kilodalton type IV collagenase (gelatinase A) in benign and malignant ovarian tumors. Lab. Invest. 69: 312-321.

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