



## 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

# Rabbit Anti Matrix Metalloproteinases 2 (MMP-2) (Gelatinase A), Whole Molecule

**Catalog No.**  
MD-15-0048

**Species**  
Human

**Isotype**  
IgG

## ***Preparation***

**Immunogen:** Native human MMP-2.  
**Purification:** Protein G affinity purified

## ***Specificity***

Rabbit antibody to the Matrix Metalloproteinase Gelatinase-A (72kD Gelatinase, MMP-2). Binds to Gelatinase-A but does not cross react with the other MMP family members (MMP-1, MMP-3, MMP-9, etc).

## ***Formulation***

**Format:** Affinity Purified, Liquid  
**Concentration:** 1mg/ml  
**Buffer:** PBS containing 50% glycerol  
**Preservative:** 0.05% Sodium Azide

## ***Storage***

Short-term (up to 6 months) store at 2-8°C as an undiluted liquid. Long term store at -20°C.

**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

**Table Summary of antibody applications and working conditions**

Options Functions	YES	NO	Not determined	Recommended Work dilution or concentration
ELISA			.	
Western Blotting	.			1:1,000 for colorimetric substrates such as BCIP/NBT. 1:5,000 for chemiluminescent substrates
Immunohistology - frozen			.	
Immunohistology - paraffin			.	
Immunohistology - resin			.	
Dot Blot			.	
Flow Cytometry			.	
Immunofluorescence staining			.	
IFA			.	

**Note:**

- Other applications are not tested yet. Optimal dilutions should be determined by each laboratory for each application.
- When used against the reduced protein, identifies bands at 72kDa and 68kDa (the pro-form and active form).
- Binds to both native and reduced forms of MMP-2. Gelatinase-A is constitutively produced in quiescent cells and tissues, and the enzyme has a high specific activity against denatured collagen.
- The low protein levels produced (pg/ml) often require concentration of cell culture media to visualize the bands by Western blotting. Gelatinase-A and Gelatinase-B may be enriched from conditioned culture media by binding to gelatin-agarose, and eluted by 10% DMSO.

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