



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

SHEEP ANTI HUMAN MMP-2

With HRP-Conjugated Second Antibody

Catalog No.	Species	Isotype
DS-PB-01357	Human	Sheep IgG

Preparation

Immunogen: Purified proMMP-2 from human cultured rheumatoid synovial fibroblasts

Preparation: Ion Exchange chromatography

Specificity

This product is specific for human matrix metalloproteinase-2 (MMP-2).

MMP-2 belongs to a family of endopeptidases which are involved in the degradation and remodelling of the ECM. No cross reactivity is seen with MMP 1, 3 or 9 or fibronectin in blots.

Formulation

Product Type: Polyclonal Antibody

Product Form: Purified IgG - liquid

Buffer Solution: Phosphate buffered saline, pH 7.2

Preservative Stabilizers: 0.09% Sodium Azide (NaN₃)

Approx. Protein Concentrations: IgG concentration 1.0 mg/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	•			1/100 – 1/1000
Immunohistology – frozen			•	
Immunohistology – paraffin	•			1/10 – 1/100
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.

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.**