



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

Rabbit Anti Disintegrin and Metalloproteinase-12 (ADAM-12), N-terminal

Catalog No. MD-15-0106	Species Human	Isotype Rabbit IgG
----------------------------------	-------------------------	------------------------------

Preparation

Host Animal: Rabbit
Immunogen: Synthetic peptide based on the N-terminal end of human ADAM-12.
Purification: Immunoaffinity purified using synthetic peptide

Specificity

Metalloproteinase-12. Reacts with human, porcine and rat. Does not react with other ADAMs.

Formulation

Format: Affinity Purified, Liquid
Concentration: 1 mg/ml
Affinity Constant: Not determined.
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 product. 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
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	.			1:1,000 (colormetric substrates) 1:5,000 (chemiluminescent substrates)
Immunohistochemistry			.	
IHC			.	
Flow Cytometry			.	
Immunofluorescence Microscopy			.	
Competitive RIA			.	

Note: Other applications are not tested yet. Optimal dilutions should be determined by each laboratory for each application.

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