



## 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 ADAM-12-L Cytoplasmic Domain

**Catalog No.**  
MD-15-0107

**Species**

**Isotype**  
IgG

### **Description**

Rabbit Antibody to A Disintegrin and Metalloproteinase-12 (ADAM-12), Meltrin-alpha, Cytoplasmic domain, Long Form

### **Preparation**

**Immunogen:** Synthetic peptide based on the cytoplasmic domain of human ADAM-12.  
**Purification:** Immunoaffinity purified using synthetic peptide.

### **Specificity**

This product recognizes full-length (ADAM-12L) but not soluble form (ADAM-12S).

### **Formulation**

**Format:** Affinity Purified, Liquid  
**Concentration:** 1mg/ml  
**Affinity Constant:** Not determined.  
**Buffer:** PBS containing 50% glycerol  
**Preservative:** 0.05% Sodium Azide

### **Applications**

Recognizes bands of 116kDa (minor band), 92kDa (major band) and cleaved products at 68kDa, 40kDa in reduced Western blots of cell lysates. We recommend a starting concentration of 1:1,000 when using colorimetric substrates and 1:5,000 for chemiluminescent substrates for Western blot. Each laboratory should determine an optimum working titer for use in its particular application. Other applications have not been tested but use in such assays should not necessarily be excluded.

**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)

---

### **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.**