



## 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 Mouse Apo B48/100

**Catalog No.**  
MD-19-0056

**Target Species**  
Mouse

### ***Description***

Rabbit Antibody to Mouse Apolipoprotein B48/100 (Apo B48/100)

### ***Preparation***

**Immunogen:** Mouse very low density lipoprotein

**Host Animal:** New Zealand Rabbit

**Format:** Neat, liquid

**Purification:** Not applicable, Heat inactivated, Product is 0.22µm filtered.

### ***Specificity***

Recognizes mouse as well as human Apo B48 and B100 as these proteins are highly conserved.

### ***Formulation***

**Buffer:** Not applicable

**Preservative:** 0.05% Sodium azide

**Concentration:** Not determined. Titer: 1:30 (RID)

### ***Storage***

Upon receipt, aliquot and store at -70°C. Avoid multiple freeze/thaw cycles.

**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

<b>Options</b> <b>Functions</b>	<b>YES</b>	<b>NO</b>	<b>Not determined</b>	<b>Recommended Work dilution or concentration</b>
ELISA			•	
Western Blotting			•	
Immunodiffusion	•			
Immunohistochemistry			•	
Immunoblotting	•			

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.**