



## 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 Atrial Natriuretic Peptide (ANP) a.a. 1-28

**Catalog No.**  
MD-17-0005

**Species**  
Human

### **Description**

Rabbit Antibody to Rat Atrial Natriuretic Peptide (ANP) (amino acid 1-28)

### **Preparation**

**Host Animal:** Rabbit  
**Immunogen:** Synthetic human ANP peptide (1-28) poly Lysine conjugated (SLRRSSCFGGRIDRIGAQSGLGCNSFRY).  
**Purification:** Not applicable

### **Specificity**

Synthetic rat ANP peptide (1-28)

### **Formulation**

**Format:** Neat, Lyophilized.  
Reconstitute in 20ul double distilled water.  
**Concentration:** Not determined  
Titer: 1:90,000 (ELISA)  
**Buffer:** Not applicable  
**Preservative:** None  
**Concentration:** Not Determined

### **Storage**

Store lyophilized product at 2-8°C. After reconstitution, store at -20°C. Avoid multiple freeze/thaw cycles.

### **Applications**

Suitable for use in RIA (1:3,000). 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.**