



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 RAT VASOPRESSIN RECEPTOR (ARG)

With HRP-Conjugated Second Antibody

Catalog No.	Species	Isotype
DS-PB-02073	Rat	IgG

Preparation

Preparation: Purified IgG prepared by affinity chromatography.

Immunogen: 21 amino acid peptide from the 3rd cytoplasmic domain of rat AVP-V2.

Specificity

This product is specific for Arginine Vasopressin Receptor (AVPR-V2). The immunizing peptide has no significant similarity with AVP V1a or V1b receptors.

Species Cross Reactivity: Based on sequence similarity, is expected to react with Pig, Mouse, Human, Bovine. Antibody reactivity and working conditions may vary between species.

Formulation

Product Type: Polyclonal Antibody

Product Form: Purified IgG - liquid

Buffer Solution: TRIS buffered saline

Approx. Protein Concentrations: IgG concentration 1.0 mg/ml

Preservative Stabilizers: 0.05% Sodium Azide (NaN₃), 0.2% Bovine Serum Albumin

Storage

12 months from date of dispatch.

Store at 4°C or at -20°C if preferred.

Storage in frost-free freezers is not recommended. This product should be stored undiluted.

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.**



Applications

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

Table Summary of antibody applications and working conditions

Options / Functions	YES	NO	Not determined	Recommended Work dilution or concentration
ELISA	•			
Western Blotting (*)	•			
Immunohistology - frozen	•			
Immunohistology - paraffin			•	
Immunohistology - resin			•	
Immunofluorescence			•	
Flow Cytometry			•	
Immunodiffusion			•	
Neutralization			•	

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

(*) A glycosylated band at approximately 62 kD is observed using membranes from rat inner medulla. N-Glycanase reduces the size of this band to 36 kD (deglycosylated form).

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.**