



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 AQUAPORIN 5

With HRP-Conjugated Second Antibody

Catalog No. DS-PB-00186	Species Rat	Isotype Rabbit IgG
-----------------------------------	-----------------------	------------------------------

Preparation

Preparation: Antisera to rat Aquaporin 5 were raised by repeated immunizations of rabbits with highly purified antigen. Purified IgG was prepared from whole serum by affinity chromatography.

Immunogen: This product was generated using a 17 amino acid synthetic peptide corresponding to the C terminus of rat Aquaporin 5 (Raina et al.1995) with the sequence "WEDHREERKKTIELTAH". See Entrez Protein accession number: NP_036911.

Specificity

This product recognizes rat Aquaporin 5 (AQP5) a member of a family of membrane water channels which exist throughout nature and include members CHIP, MIP and other aquaporins. AQP5 is expressed at locations implicating its involvement in generation of tears, saliva and pulmonary secretions.

Formulation

Product Type: Polyclonal Antibody

Product Form: Purified IgG - liquid

Buffer Solution: 100mM Tris. pH7.5

Preservative Stabilizers: 0.05 Sodium Azide (NaN₃)

Approx. Protein Concentrations: Ig concentration 1.0 mg/ml

Protein Molecular Weight: Rat AQP5 has a predicted molecular mass of 28.4kDa.

Storage

18 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.**



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	.			0.5 – 1.0 ug/ml
Western Blotting(*)	.			1 – 10 ug/ml
Immunohistology - frozen			.	
Immunohistology - paraffin			.	
Immunohistology - resin			.	
Immunoprecipitation			.	
Flow Cytometry			.	
Immunofluorescence staining			.	
Neutralization			.	

Note: Other applications are not tested yet. Optimal dilutions should be determined by each laboratory for each application.
 * detects a band/s of approximately 27-28kDa in salivary gland cell lysates under reducing conditions.

Secondary Antibody Applications

Options / Functions	YES	NO	Not determined	Recommended Work dilution or concentration
Immunoassay (ELISA, Western blot)	.			1:5000-1:10000

Reference:

Raina, S. et al. (1995) Molecular cloning and characterization of an aquaporin cDNA from salivary, lacrimal, and respiratory tissues. J Biol Chem. 270: 1908-1912.

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