



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

Sheep anti Mouse Albumin

Catalog No.
MD-14-0016

Species
Human

Preparation

Host Animal: Sheep
Immunogen: Mouse albumin purified from serum. Specific by IEP
Purification: Ion exchange chromatography. Zone electrophoresis is performed to ensure that only a gamma-mobility component is present. Product is 0.2 μ m filtered.

Specificity

Gives a single arc when tested by IEP against mouse serum. Identity has been confirmed by double diffusion (Ouchterlony) against mouse serum and a known anti-mouse albumin. Typical cross-reactivity against the sera of the species listed below has been assessed by double diffusion using 10 μ l of antisera against 1 μ l of species sera. The results are intended as a guideline only. It is possible that cross-reactivity against species listed as negative may be observed when using different test conditions or techniques.

Cat	+	Chicken	+	Cow	-	Dog	+	Goat	-	Guinea Pig	+
Horse	+	Human	+	Mouse	+	Pig	+	Rabbit	+	Rat	+

Formulation

Host Animal: Sheep
Format: Purified, Liquid
Concentration: 27.8mg/ml (OD280nm E^{1%} = 14.5)
Buffer: Glycine buffered saline, pH 7.4
Preservative: 0.099% Sodium azide, 0.1% E-Amino-n-Caproic Acid, 0.01% Benzamidine, 1mM EDTA

**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

Storage

Store at 2-8°C. Slight precipitation can occur on storage which may be removed by centrifugation, and should not affect performance characteristics.

Avoid multiple freeze-thaw cycles.

Centrifuge before opening to ensure complete recovery of vial contents.

Applications

Table Summary of antibody applications and working conditions

Options Functions	YES	NO	Not determined	Recommended Work dilution or concentration
Radial Immunodiffusion (RID)	•			10ul antiserum/cm ² in gel vs 5ul mouse serum, 1:50 - 1:200
Double Diffusion	•			10ul antiserum vs 3ul mouse serum (1:50 dilution)
Immunoelectrophoresis (IEP)	•			100ul antiserum vs 5ul mouse Serum (1:50 dilution)

Note: (1) Other applications are not tested yet. Optimal dilutions should be determined by each laboratory for each application.

(2) The use of 3% PEG 6000 with 1.2% agarose in a suitable buffer (such as TBE or Tris-barbital pH >8.2) is recommended.

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