



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 Mouse Apolipoprotein E (Apo E)

Catalog No.
MD-19-0078

Target Species
Mouse

Preparation

Host Animal: Rabbit
Immunogen: Purified Mouse Apolipoprotein E from pooled mouse plasma high and very low density lipoproteins
Purification: Not applicable
Heat inactivated
Product is 0.22 μ m filtered.

Specificity

Species specific: extremely low recognition of human Apo E. No data available regarding rat activity.

Formulation

Format: Neat, Liquid
Concentration: Not determined
Titer: 1:40 (RID)
Buffer: Not applicable
Preservative: 0.05% Sodium azide

Applications

Suitable for use in Radial immunodiffusion, immuno double-diffusion, rocket IEP and Western blot. Starting dilution for Western blot: 1:500 - 1:1,000. 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.

Storage

Upon receipt, aliquot and store (up to 1 year) 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
Email: info@raybiotech.com

References:

1. Hayhurst, G.P., et al., (2001), "Hepatocyte Nuclear Factor 4 alpha (Nuclear Receptor 2A1) is essential for maintenance of Hepatic Gene Expression and Lipid Homeostasis", Molecular and Cellular Biology, **21**(4): 1393-1403
2. Rees, D., et al., (1999), "Apolipoprotein A-I stimulates secretion of Apolipoprotein E by Foam Cell Macrophages", J. Biol. Chem., **274**(39): 27925-27933
3. Vaisman, B.L., et al., (2001), "ABCA1 overexpression leads to hyperalphalipoproteinemia and increased biliary cholesterol excretion in transgenic mice", J. Clin. Invest., **108**: 303-309

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