



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 Goat Anti Mouse Apolipoprotein AI (Apo AI)

Catalog No.
MD-19-0033

Species
Mouse

Description

Goat Antibody to Mouse Apolipoprotein AI (Apo AI)

Preparation

Host Animal: Goat
Immunogen: Mouse Apolipoprotein AI from pooled mouse plasma high density lipoprotein.

Specificity

Recognizes mouse and rat Apo AI. Does not react well with human Apo AI.

Formulation

Format: Neat, Liquid
Concentration: Not determined
Titer: 1:40 (RID)
Buffer: Not applicable
Preservative: 0.1% Sodium azide, 100U/ml penicillin, 10ug/ml streptomycin, 25ng/ml amphotericin B

Applications

Table Summary of antibody applications and working conditions

Options / Functions	YES	NO	Not determined	Recommended Work dilution or concentration
ELISA			•	
Radial Immunodiffusion Assay	•			
Western Blot			•	

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

**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 (up to 1 year) at -70°C .

References

- 1)Furbee, J.W., et al., (2002), "In vivo contribution of LCAT to apolipoprotein B lipoprotein cholesteryl esters in LDL receptor and apolipoprotein E knockout mice", J. Lipid Res., **43**: 428-437
- 2)Furbee, J.W., et al., (2001), "Alteration of plasma HDL cholesteryl ester composition with transgenic expression of a point mutation (E149A) of human LCAT", J. Lipid Res., **42**: 1626-1635

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