



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 DataSheet Mouse Anti CD14 Monocytes

Catalog No.
MD-11-0022

Source
Cell Culture

Isotype
IgG_{2a}

Description:	Monoclonal Antibody to CD14 Monocytes
Host Animal:	Mouse
Immunogen:	Thymocytes and peripheral blood leucocytes.
Specificity:	Recognizes a (Mr 55 kDa) cell surface glycoprotein expressed on most peripheral blood monocytes. CD14 antigen is expressed on 70-95% of peripheral blood monocytes. It is also found on other cells of myeloid lineage including granulocytes, dendritic reticulum cells and pleural phagocytic cells.
Format:	Purified Liquid
Purification:	Protein G chromatography
Concentration:	100ug/ml (OD280nm)
Buffer:	0.01M PBS, pH7.2 containing 1.0% BSA
Preservative:	0.09% Sodium azide
Storage:	Store at 2–8° C. Avoid multiple freeze/thaw cycles.
Applications:	May be used for enumerating mature monocytes in peripheral blood. We recommend using 1ug to stain 1.0 x 10 ⁶ cells in flow cytometric applications. 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.
References:	<ol style="list-style-type: none">1. Hogg, N.H., et al., (1984), <i>Immunology</i>, 53:7532. Linch, D.C., et al., (1984), <i>Blood</i>, 63:5663. Khim, Y., et al., (1985), <i>Int. J. Cancer</i>, 36:4334. Leucocyte Typing: 3rd Workshop: code No. 203 4th Workshop: code No.M24.

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