

# CD14 (human) monoclonal antibody (biG 2/RoMo-1)

The CD14 glycoprotein, gp 55, is present on most monocytic and macrophage like cell types: monocytes, macrophages, Kupffer cells, pleural phagocytic cells and dendritic reticular cells. CD14 is also observed on granulocytes and activated or transformed B-cells. Furthermore CD14 is present in a soluble form in human serum, urine and other body fluids. The CD14 molecule has been reported to be a receptor for endotoxin.

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 1

[View Online »](#)

## Ordering Information

[Order Online »](#)

ALX-804-498-C100	100µg
------------------	-------

## Manuals, SDS & CofA

[View Online »](#)

# Handling & Storage

Handling                                      Avoid freeze/thaw cycles.

Long Term Storage                      -20°C

Shipping                                      Blue Ice

# Regulatory Status

RUO - Research Use Only

# Product Details

Application                                      ELISA, Flow Cytometry, IP

Application Notes                              Does **not** inhibit binding of LPS to CD14.

Clone    biG 2/RoMo-1

Formulation                                      Lyophilized from PBS. Contains no preservatives.

Host    Mouse

Immunogen                                      Human monocytes.

Isotype    IgG2a

Purity Detail                                      Protein G-affinity purified.

Reconstitution                                      Reconstitute with 100µl distilled water.

Species Reactivity                                      Human

Specificity    Recognized epitope: aa 147-152 of human CD14.

Technical Info / Product Notes                                      Binding titer at human CD14 transfected CHO-cells (FACS) is 1:5000.

UniProt ID    P08571

Worry-free Guarantee                                      This antibody is covered by our [Worry-Free Guarantee](#) .



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)