

# MAB-hMR

Neutralizing IgG monoclonal antibody to human mannose receptor

Catalog # mab-hmr

For research use only, not for diagnostic or therapeutic use

Version # 12J16-MM

## PRODUCT INFORMATION

### Content

100 µg purified MAb-hMR, provided azide-free and lyophilized

**Clone:** 15-2

**Isotype:** Mouse IgG1

**Formulation:** 0.2 µm filtered antibody solution in PBS, containing 0.1% bovine serum albumin

### Antibody resuspension

Add 1 ml of sterile water to obtain a concentration of 0.1 mg/ml.

### Storage

- Product is shipped at room temperature. Store lyophilized antibody at -20°C. Product is stable for 1 year.

- Reconstituted antibody is stable 1 month when stored at 4°C and 6 months when aliquoted and stored at -20°C. Avoid repeated freeze-thaw cycles.

## DESCRIPTION

MAB-hMR is a monoclonal antibody specific for human mannose receptor (CD206). MAB-hMR has been selected for its ability to efficiently neutralize the biological activity of the human mannose receptor.

## BACKGROUND

The mannose receptor (MR), also known as CD206, is a type I transmembrane protein and a member of the C-type lectin family. MR is primarily present on the surface of macrophages and dendritic cells and plays a role in both innate and adaptive immunity. MR binds high-mannose structures on a wide range of gram positive and gram negative bacteria, yeasts, parasites and mycobacteria.

MR was identified as one of the first fungal pattern recognition receptor (PPR) inducing NF-κB activation and the production of cytokines, such as IL-8 and IL-6<sup>1</sup>. In adaptive immunity, it has been shown that MR is important in antigen uptake by dendritic cells which essential for the induction of antigen-specific T-cell responses<sup>2</sup>.

## APPLICATIONS

- MAb-hMR is a neutralizing antibody, it blocks human mannose receptor-induced cellular activation. MAB-MR prevented binding of tissue-type plasminogen activator (t-PA) to the immobilized human MR<sup>3</sup>. The exact concentration of antibody required to neutralize human mannose receptor activity is dependent on the agonist, cell type and growth conditions.

- MAB-hMR can be used for flow cytometry. For flow cytometry, MAb-hMR has been tested at 500 - 2000 ng/10<sup>6</sup> cells with PE goat anti-mouse kappa light chain secondary antibody for indirect immunofluorescence staining.

***Note:** We recommend to test the intracellular and cell surface expression of hMR.*

- MAB-hMR can be used for Western blotting. MAB-hMR recognized the 175 kDa mannose receptor in a crude sample extracts<sup>3</sup>.

## REFERENCES

**1. Willment JA. & Brown GD., 2007.** Dectin-1 mediates the biological effects of beta-glucans. *J Exp Med.* 197: 1119- 24. **2. Autenrieth SE. & Autenrieth IB., 2009.** Variable antigen uptake due to different expression of the macrophage mannose receptor by dendritic cells in various inbred mouse strains. *Immunology.* 127(4):523-9. **3 Barrett-Bergshoeff M. et al., 1997.** Monoclonal antibodies against the human mannose receptor that inhibit the binding of tissue-type plasminogen activator. *Thromb Haemost.* 77(4):718-24.

---

### TECHNICAL SUPPORT

Toll free (US): 888-457-5873

Outside US: (+1) 858-457-5873

Europe: +33 562-71-69-39

E-mail: [info@invivogen.com](mailto:info@invivogen.com)

Website: [www.invivogen.com](http://www.invivogen.com)



3950 Sorrento Valley Blvd. Suite 100  
San Diego, CA 92121 - USA