

# Anti-mDectin-1-IgG

Neutralizing monoclonal antibody against mouse Dectin-1

Catalog # mabg-mdect

For research use only, not for diagnostic or therapeutic use

Version # 16C04-MM

## PRODUCT INFORMATION

### Content

100 µg purified anti-mDectin-1-IgG, provided azide-free and lyophilized

**Clone:** R1-8g7

**Isotype:** Rat IgG2a

**Formulation:** 0.2 µm filtered solution in 68 mM phosphate buffer with 91 mM glycine, 5% w/v saccharose and stabilizing agents

### 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 6 months.

- 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.

### Quality control

- The absence of bacterial contamination (e.g. endotoxins and peptidoglycans) is controlled using HEK-Blue™ TLR2 and HEK-Blue™ TLR4 cells.

- The antibody is validated for neutralization and flow cytometry using HEK-Blue™ Dectin cells.

## BACKGROUND

Dectin-1 is a member of the C-type lectin receptor (CLR) family and plays an important role in antifungal innate immunity<sup>1</sup>. In humans and mice, Dectin-1 is alternatively spliced into two major isoforms, a full-length A isoform and a 'stalkless' B isoform. Dectin-1 is expressed on phagocytic cells, including macrophages and neutrophils. Dectin-1 is a specific receptor for β-glucans<sup>2</sup>. β-Glucans are glucose polymers found in the cell walls of fungi, including the yeasts *Saccharomyces cerevisiae* and *Candida albicans*. Upon binding to its ligand, Dectin-1 triggers phagocytosis and signaling through the kinase Syk and the adaptors CARD9-Bcl10-Malt1 leading to the production of reactive oxygen species (ROS), the activation of NF-κB and the subsequent production of pro-inflammatory cytokines<sup>3,4</sup>. Dectin-1 signaling has been shown to collaborate with Toll-like receptor 2 (TLR2) signaling to enhance the responses triggered by each receptor<sup>4,5</sup>.

**1. Herre J. et al., 2004.** The role of dectin-1 in antifungal immunity. *Crit Rev Immunol.* 24(3):193-204. **2. Brown GD. et al., 2003.** Dectin-1 mediates the biological effects of beta-glucans. *J Exp Med.* 197: 1119- 24. **3. Gross O. et al., 2006.** Card9 controls a non-TLR signaling pathway for innate anti-fungal immunity. *Nature.* 442:651- 6. **4. Dennehy KM, & Brown GD., 2007.** The role of the beta-glucan receptor Dectin-1 in control of fungal infection. *J Leukoc Biol.*82(2):253-8. **5. Gantner BN. et al., 2003.** Collaborative induction of inflammatory responses by dectin-1 and Toll- like receptor 2. *J Exp Med.* 197:1107- 17. **6. Seifert L. et al., 2015.** Dectin-1 regulates hepatic fibrosis and hepatocarcinogenesis by suppressing TLR4 signaling pathways. *Cell Rep.* 13(9):1909-21.

## TECHNICAL SUPPORT

InvivoGen USA (Toll-Free): 888-457-5873

InvivoGen USA (International): +1 (858) 457-5873

InvivoGen Europe: +33 (0) 5-62-71-69-39

InvivoGen Hong Kong: +852 3-622-34-80

E-mail: info@invivogen.com

## DESCRIPTION

Anti-mDectin-1-IgG is a monoclonal rat IgG2a antibody against mouse Dectin-1 (mDectin-1). It recognizes both mDectin-1 isoforms A and B. This antibody was screened for flow cytometry and neutralization activity. Anti-mDectin-1-IgG is produced in hybridomas and purified by affinity chromatography with protein G/agarose.

## APPLICATION

Anti-mDectin-1-IgG is a neutralizing antibody, it blocks mDectin-1-induced cellular activation. This antibody is also intended for use in flow cytometry. It can also be used for immunohistochemistry<sup>6</sup>.

### Applications tested

1. Anti-mDectin-1-IgG has been tested in neutralizing experiments performed in HEK293 cells that were transfected with the mDectin-1 gene and other genes from the Dectin-1 signaling pathway. These cells also stably express a secreted embryonic alkaline phosphatase (SEAP) reporter gene.

Transfected cells were incubated with 100 ng-10 µg/ml anti-mDectin-1-IgG for 1 hour prior to the addition of the Dectin-1 ligand HKCA (10<sup>8</sup> cells/ml).

Following an overnight incubation, neutralization of Dectin-1 signaling by anti-mDectin-1-IgG was determined by assessing SEAP production using QUANTI-Blue™, a SEAP detection medium. QUANTI-Blue™ turns blue following Dectin-1 stimulation but remains pink if neutralization occurs.

2. For flow cytometry, anti-mDectin-IgG has been tested at 500-2000 ng/10<sup>6</sup> cells with PE goat anti-rat IgG2a secondary antibody for indirect immunofluorescence staining.

## RELATED PRODUCTS

Product	Catalog Code
QUANTI-Blue™	rep-qb1
pUNO1-mDectin1a	puno-mdectin1a
pUNO-mDectin1b	puno-mdectin1b
pNiFty2-SEAP	pnipty2-seap
<b>Dectin-1 Agonists</b>	
Curdlan AL	trl-cura
HKCA (Heat killed <i>C. albicans</i> )	trl-hkca
HKSC (Heat killed <i>S. cerevisiae</i> )	trl-hksc
WGP Dispersable (Whole Glucan particles)	trl-wgp
Zymosan Depleted	trl-dzn