

PRODUCT INSERT

MONOCLONAL ANTIBODIES TO HUMAN IL-16

| Product Code | Form | Volume | Tests | Excitation (nm) | Peak Emission (nm) | Matching Isotype Controls | |
|--------------|----------|--------|----------|-----------------|--------------------|---------------------------|-------------|
| MHCIL1600 | Purified | 1.0 ml | 200 min. | N/A | N/A | Mouse IgG2a Purified | Code MG2a00 |
| MHCIL1601 | FITC | 1.0 ml | 100 min. | 488 | 525 | Mouse IgG2a FITC | Code MG2a01 |
| MHCIL1601-3 | FITC | 3.0 ml | 300 min. | | | | |
| MHCIL1604 | R-PE | 0.5 ml | 50 min. | 488 | 575 | Mouse IgG2a R-PE | Code MG2a04 |
| MHCIL1604-3 | R-PE | 3.0 ml | 300 min. | | | | |

PRODUCT DESCRIPTION

Mouse monoclonal antibody to human IL-16

Clone: 14.1

Isotype: Mouse IgG2a

Lot No.: See label **Expiration:** See label

Buffer: Phosphate buffered saline (PBS)

Preservative: 0.1% *sodium azide*. Sodium azide is an extremely toxic and dangerous compound particularly when combined with acids or metals. Solutions containing sodium azide should be disposed of properly.

Stabilizer: For conjugated products only, a highly purified grade of BSA has been added as a stabilizing agent.

STORAGE & HANDLING

Store reagents at 2-8°C. For fluorochrome-conjugated antibodies only, light exposure should be avoided. Use dim light during handling, incubation with cells and prior to analysis. It is recommended that investigators dilute only the quantity to be used within one week.

PRODUCT CHARACTERIZATION

Antigen Specificity: According to the literature this antibody recognizes IL-16¹⁻⁹. This antibody recognizes both the 80 kDa precursor and 17 kDa secreted forms of IL-16.

Leukocyte Workshop Status: N/A

PRODUCT QUALITY CONTROL

Each lot is tested by flow cytometry using peripheral blood leukocytes. Based on this testing it is recommended that 5 µl of antibody be used per 1 x 10⁶ cells. Because results may vary, it is suggested that each investigator determine the optimal amount of antibody to be used for each application.

REFERENCES:

- Center, D.M., and W.W. Cruikshank. 1982. *J. Immunol.* 128: 2563.
- Cruikshank, W.W., and D.M. Center. 1982. *J. Immunol.* 128: 2569.
- Biddison, W.E., D.D. Taub, W.W. Cruikshank, D.M. Center, E.W. Connor, and K. Honma. 1997. *J. Immunol.* 158: 3046.
- Theodore, A.C., D.M. Center, J. Nicoll, G. Fine, H. Kornfeld, and W.W. Cruikshank. 1996. *J. Immunol.* 157: 1958.
- Maciaszek, J.W., N.A. Parada, W.W. Cruikshank, D.M. Center, H. Kornfeld, and G. Viglianti. 1997. *J. Immunol.* 158: 5.
- Cruikshank, W.W., K. Lim, A.C. Theodore, J. Cook, G. Fine, P.F. Weller, and D.M. Center. 1996. *J. Immunol.* 157: 5240.
- Center, D.M., H. Kornfeld, and W.W. Cruikshank. 1996. *Immunol. Today* 17: 476.
- Zhang, Y., D.M. Center, D.M. Wu, W.W. Cruikshank, J. Yuan, D.W. Andrews, and H. Kornfeld. 1998. *J. Biol. Chem.* 273: 1144.
- Laberge, S., P. Ernst, O. Ghaffar, W.W. Cruikshank, H. Kornfeld, D.M. Center, and Q. Hamid. 1997. *Am. J. Respir. Cell Mol. Biol.* 17: 193.

FOR RESEARCH USE ONLY. . . NOT FOR THERAPEUTIC OR IN VITRO DIAGNOSTIC USE

www.invitrogen.com

Invitrogen Corporation • 542 Flynn Rd • Camarillo • CA 93012 • Tel: 800.955.6288 • E-mail: techsupport@invitrogen.com

PI: L11271

(Rev 12/08) DCC-08-1818

Important Licensing Information - These products may be covered by one or more Limited Use Label Licenses (see the Invitrogen Catalog or our website, www.invitrogen.com). By use of these products you accept the terms and conditions of all applicable Limited Use Label Licenses. Unless otherwise indicated, these products are for research use only and are not intended for human or animal diagnostic, therapeutic or commercial use.