



Mouse (monoclonal) Anti-Human α -chain of the Interleukin-2 Receptor CD25

PRODUCT ANALYSIS SHEET

Catalog Number:	AHS2512
Lot Number:	See product label
Quantity/Volume:	100 μ g/100 μ L
Clone Number:	143-13
Isotype:	IgG ₁
Form of Antibody:	Purified immunoglobulin in phosphate buffered saline, pH 7.2.
Preservation:	None, 0.22 micron sterile filtered.
Purification:	Purified from ascites by Protein G affinity chromatography.
Immunogen	Recognizes an antigen of 55-60 kDa.
Myeloma/Fusion Partners:	Mouse NSI cell line with mouse spleen lymphocytes.
Specificity:	Reacts with Epitope Cluster B of the human Interleukin-2 receptor alpha chain.
Applications:	Flow cytometry and immunofluorescence.
Suggested Working Dilutions:	Use approximately 0.5–1.0 μ g to label 10 ⁶ cells. The optimal antibody concentration should be determined for each specific application.
Storage:	Store in aliquots at –20°C. Avoid repeated freeze/thaw cycles.

This product is for research use only. Not for use in diagnostic procedures.

www.invitrogen.com

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

PI AHS2512

(Rev 11/08) DCC-08-1089

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.

References:

- Knapp, W. et al. (Eds.) Leucocyte Typing IV, Oxford University Press, Oxford, 1989.
- Plana, M., O. Vinas, O. De la Calle-Martin, F. Lozano, J. Ingles-Esteve, M. Romero, J. Alberola-Ila, J. Yague, R. Vilella, and J. Vives (1991) Induction of Interleukin-2 (IL-2) and Interferon-gamma and enhancement of IL-2 receptor expression by a CD26 monoclonal antibody. *Eur. J. Immunol.* 21:1085-1088.
- Oravecz, T., G. Roderiquez, J. Koffi, J. Wang, M. Ditto, D.C. Bou-Habib, P. Lusso, and M.A. Norcross (1995) CD26 expression correlates with entry, replication and cytopathicity of monocyctotropic HIV-1 strains in a T-cell line. *Nat. Med.* 1(9):919-926.
- Bluman, E.M., K.J. Bartynski, B.R. Avalos, and M.A. Caligiuri (1996) Human natural killer cells produce abundant macrophage inflammatory protein-1 alpha in response to monocyte-derived cytokines. *J. Clin. Invest.* 97(12):2722-2727.
- Blauvelt, A., H. Asada, M.W. Saville, V. Klaus-Kovtun, D. J. Altman, R. Yarchoan, and S.I. Katz (1997) Productive infection of dendritic cells by HIV-1 and their ability to capture virus are mediated through separate pathways. *J. Clin. Invest.* 100(8):2043-2053.
- Canfield, S.M. and A.Y. Khakoo (1999) The nonintegrin laminin binding protein (p67 LBP) is expressed on a subset of activated human T lymphocytes and, together with the integrin very late activation antigen-6, mediates avid cellular adherence to laminin. *J. Immunol.* 163(6):3430-3440.
- Hladik, F., G. Lentz, E. Delpit, A. McElroy, and M.J. McElrath (1999) Coexpression of CCR5 and IL-2 in human genital but not blood T cells: Implications for the ontogeny of the CCR5(+) Th1 phenotype. *J. Immunol.* 163(4):2306-2313.

This product is for research use only. Not for use in diagnostic procedures.

www.invitrogen.com

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

PI AHS2512

(Rev 11/08) DCC-08-1089

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.