



Mouse (monoclonal) Anti-Human T-Cell Subset CD28 Unconjugated, azide free

PRODUCT ANALYSIS SHEET

| | |
|-------------------------------------|----------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|
| Catalog Number: | AHS2812 |
| Lot Number: | See product label |
| Quantity/Volume: | 0.1 mg/0.1 mL |
| Clone Number: | 152-2E10 |
| 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: | Allogenic stimulated human peripheral blood mononuclear cells. |
| Myeloma/Fusion Partners: | Mouse spleen cells fused with the NS1 cell line. |
| Specificity: | <p>This monoclonal antibody recognizes CD28, a cell surface glycoprotein expressed by mature CD3⁺ thymocytes, most peripheral T lymphocytes, and plasma cells. CD28 serves as a receptor for the ligands CD80 and CD86. The binding of these ligands to CD28 stimulates T-cell effector functions, including T-cell dependent antibody production. This monoclonal antibody immunoprecipitates an antigen with M_r=41 kDa, as determined by SDS PAGE performed under reducing conditions.</p> <p>The specificity of this antibody was studied at the Fifth Leukocyte Typing Workshop.</p> |
| Applications: | This monoclonal antibody is suitable for use in immunoprecipitation, flow cytometry, and immunofluorescence. |
| Suggested Working Dilutions: | For flow cytometry, use approximately 0.25-1.0 µg to label up to 10 ⁶ cells. The optimal concentration should be determined for each specific application. |
| Storage: | Store at -20°C. Upon initial thawing, apportion the antibody into working aliquots and store at -20°C. Avoid repeated freeze/thaw cycles. The addition of a preservative such as 0.1% sodium azide is also recommended for long term storage. |
| References: | <p>Fox, D.A., J. Saunders, and L.A. Diaz (1995) Classification and functional properties of normal T-cell subsets. In Leukocyte Typing V, edited by S.F. Schlossman, L. Boumsell, W. Gilks, J.M. Harlan, T. Kishimoto, C. Morimoto, J. Ritz, S. Shaw, R. Silverstein, T. Springer, T.F. Tedder, and R.F. Todd. Oxford University Press. pp. 439, 442.</p> <p>Lara-Marquez, M.L., M.S. O'Dorisio, T.M. O'Dorisio, M.H. Shah, and B. Karacay (2001) Selective gene expression and activation-dependent regulation of vasoactive intestinal peptide receptor type 1 and type 2 in human T cells. J. Immunol. 166 (4):2522-2530.</p> |

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 AHS2812

(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.