

PRODUCT INFORMATION



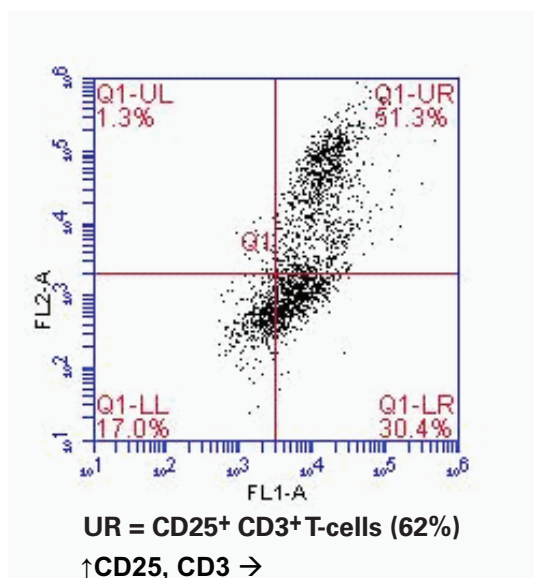
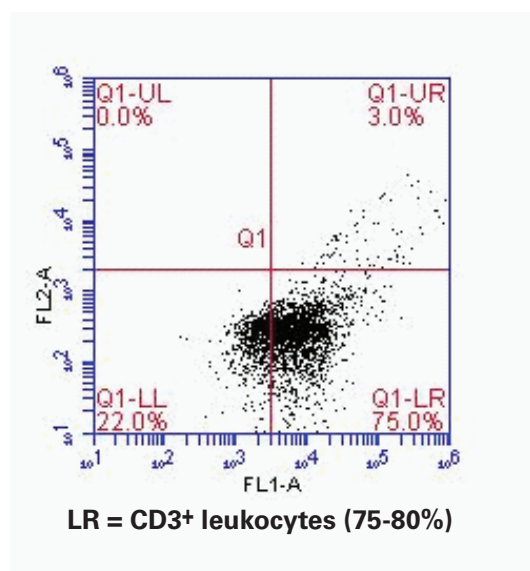
CD25 Monoclonal PE Antibody (Clone 143-13)

Item No. 10339

Overview and Properties

Contents: 100 µg of affinity purified, PE labeled monoclonal antibody
Synonyms: Interleukin-2 Receptor α Subunit, IL-2Ra, p55, Tac
Species Reactivity: (+) Human CD25
Uniprot No.: P01589
Form: Liquid
Storage: 4°C (as supplied)
Stability: ≥1 year
Storage Buffer: TBS, pH 7.4 with 50% glycerol, 5 mg/ml BSA, and 0.02% sodium azide
Clone: 143-13
Host: Mouse
Isotype: IgG₁
Applications: Flow cytometry (5-10 µl per test), frozen tissue sections, and immunofluorescence; assay dependent dilution. Does not work for western blotting.

Images



WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 09/26/2018

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

Interleukin-2 (IL-2) elicits its effects by binding to a multi-subunit cell-surface receptor that is inducible by various mitogens.¹⁻³ The highest affinity for IL-2 is achieved when all three subunits (IL-2R α /CD25, IL-2R β /CD122, and IL-2R γ /CD132) are expressed together forming a heterotrimeric complex.⁴ CD25 is inducible in T-cells by antigen, mitogen, or antibodies bound to the T-cell receptor CD3.^{2,5}

Binding of IL-2 to IL-2R on NK cells results in their conversion to LAK cells, which release several other cytokines (TNF, INF- γ , and GM-CSF).^{4,6} Activated T-cells will respond to IL-2 by proliferating and secreting cytokines and cytolytic molecules. Activated β -cells will proliferate and secrete antibodies.

References

1. Ruscetti, F.W. and Gallo, R.C. Human T-lymphocyte growth factor: Regulation of growth and function of T lymphocytes. *Blood* **57(3)**, 379-394 (1981).
2. Robb, R.J., Munck, A., and Smith, K.A. T cell growth factor receptors. Quantification, specificity, and biological relevance. *J. Exp. Med.* **154**, 1455-1474 (1981).
3. Leonard, W.J., Krönke, M., Pfeffer, N.J., *et al.* Interleukin 2 receptor gene expression in normal human T lymphocytes. *Proc. Natl. Acad. Sci. USA* **82**, 6281-6285 (1985).
4. Waldmann, T.A. The IL-2/IL-2 receptor system: A target for rational immune intervention. *Immunol. Today* **14**, 264-269 (1993).
5. Venuta, S., Mertelsmann, R., Welte, K., *et al.* Production and regulation of interleukin-2 in human lymphoblastic leukemias studied with T-cell monoclonal antibodies. *Blood* **61(4)**, 781-789 (1983).
6. Smith, K.A. The interleukin 2 receptor. *Annu. Rev. Cell Biol.* **5**, 397-425 (1989).
7. Kaplan, G., Cohn, Z.A., and Smith, K.A. Rational immunotherapy with interleukin 2. *Biotechnology* **10**, 157-162 (1992).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM