

NEW PRODUCT SINGLE COLOR ANTIBODIES

CD105-PE

PN B76299

Antibody

Specificity: Human
 Clone: TEA3/17.1.1
 Isotype: IgG1 kappa mouse

Reagent

Line: IOTest
 Format: Liquid
 Size: 1 mL
 Status: ASR

Example of results

Antibody characteristics

The TEA3/17.1.1 monoclonal antibody reacts with endothelial cells from various tissues, monocytes, pre B lymphocytes and proerythroblasts from bone marrow.

The TEA3/17.1.1 monoclonal antibody has been assigned to the CD105 cluster of differentiation at the sixth International Workshop on Human Leucocyte Differentiation Antigens held in Kobe, Japan, in 1996 (WS code : E105).

Antigen characteristics

Other name: Endoglin
 Gene Name: ENG
 Gene Location: 9q33-q34.1
 Molecular Weight: 170-180 kDa

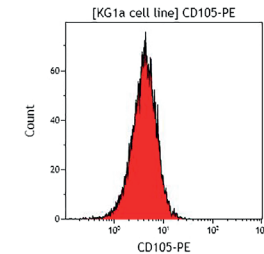
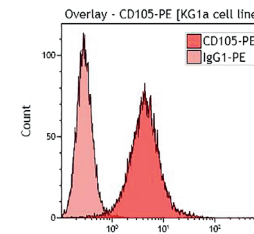
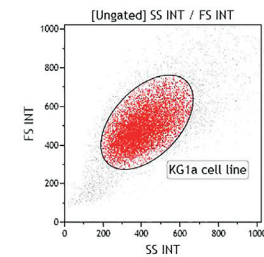
The CD105 antigen is a RGD-containing molecule which is present on the cell surface as a disulphide-linked homodimer. The molecule is a co-receptor for transforming growth factor (TGF) β 1 and β 3 (1). It is expressed mainly on endothelial cells, but also strongly in pre-B lymphocytes and in a subpopulation of monocytes.

CD105 is a regulatory protein of TGF-beta receptor complex. It is a modulator of TGF-beta 1 cell response.

Formats available

	Size	Line	Status	Part #		Size	Line	Status	Part #
PE	1 mL	IOTest	ASR	B76299	PC7	0.5 mL	IOTest	ASR	B43293

ASR: Analyte Specific Reagent. Analytical and performance characteristics are not established.



KG1a cell line

References

- ◆ Lopez-Casillas et al, "Structure and expression of the membrane proteoglycan betaglycan, a component of the TGF-beta receptor system", 1991, Cell, 67, 785-795.
- ◆ Letarte M, Bourdeau A, Vera S, Pece N, Greaves A, Dignat-George F, Mutin M, Kraling B, Linask K, O'Brien ER, et al. EC2 CD105 Workshop Panel Report. In: Kishimoto T, Kikutani H, von dem Borne AEGK, Goyert SM, Mason DY, Miyasaka M, Moretta L, Okumura K, Shaw S, Springer TA, et al, eds. Leukocyte Typing VI. New York, NY: Garland Publishing Inc; 1997:703-708.