

Biosciences Human CD Marker Chart



Reviewed by H1DA

bdbiosciences.com/go/humancdmarkers

Table 1: Human CD Marker Chart (CDs 1-100). Columns include Alternative Name, Ligands & Associated Molecules, and expression in various cell types (T Cell, B Cell, Dendritic Cell, etc.).

Table 2: Human CD Marker Chart (CDs 101-200). Columns include Alternative Name, Ligands & Associated Molecules, and expression in various cell types (T Cell, B Cell, Dendritic Cell, etc.).

Table 3: Human CD Marker Chart (CDs 201-360). Columns include Alternative Name, Ligands & Associated Molecules, and expression in various cell types (T Cell, B Cell, Dendritic Cell, etc.).

Visual representation of cell types and their key markers. Includes T Cell (CD3, CD4, CD8), B Cell (CD20), Dendritic Cell (CD123), NK Cell (CD56), Stem Cell/Precursor (CD34), Macrophage/Monocyte (CD33), Granulocyte (CD66), Platelet (CD61, CD62), Erythrocyte (CD235), Endothelial Cell (CD146), and Epithelial Cell (CD236).

*Positive refers to any expression of this CD marker either in all cells, a subset, or upon activation. For Research Use Only. Not for use in diagnostic or therapeutic procedures. BD, BD Logo and all other trademarks are property of Becton, Dickinson and Company. © 2010 BD

The new CDs were established at the H1DA9 meeting held in Barcelona 2010. For additional information regarding CD molecules, please visit www.hcdm.org.