

CDw78 Monoclonal Mouse Antibody (DF1588)

Product Description

This antibody reacts with a carbohydrate determinant, identified as CDw78 (Workshop V; Code CDw78.2). Epitope recognized by this MAb and Leu21 is the same. It is present on some immature and some peripheral blood B cells, and some B cell lines. Its expression is increased after B-cell activation in vitro. CDw78 is also found on tissue macrophages and on epithelial cells, but not on T cells, NK cells, monocytes, granulocytes and myeloid tissues. It appears that CDw78 is a determinant on a major histocompatibility complex (MHC) class II subpopulation.

Product attributes

Antibody number	#0353
Antibody reactivity (target)	CDw78
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	DF1588
Isotype	IgG1, kappa
Molecular weight	Not Known
Synonyms	Leu21; Ba antigen; LO panBa
Human gene symbol	Not Known
Entrez gene ID	Not Known
SwissProt	Not Known
Unigene	Not Known
Immunogen	Human leucocytes
Antibody target cellular localization	Plasma membrane
Species reactivity	Human
Antibody application notes	For coating for ELISA, order Ab without BSA. Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody. Optimal dilution and staining procedure for a specific application should be determined by user. Recommended starting concentrations for titration are 1-2 ug/mL for most applications, or 1 ug/million cells/100 uL for flow cytometry
Positive control	Raji and human peripheral blood lymphocytes. Human tonsil and lymph node.
Shipping condition	Room temperature
Storage Conditions	Store at 2 to 8 °C. Protect fluorescent conjugates from light. Note: store BSA-free antibodies at -10 to -35 °C
Regulatory status	For research use only (RUO)
Antibody/conjugate formulation	Conjugates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide. HRP conjugates: 0.1 mg/mL in PBS/0.05% BSA. Purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% azide. Purified, BSA-free: 1 mg/mL in PBS without azide
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Cell/tissue expression	Macrophage/monocyte, B-cells, Epithelial cells
Antibody research areas	Immunology