

Erythrocyte Specific Monoclonal Mouse Antibody (SFL23.6)

Product Description

Mouse Anti-Human Erythrocytes Clone SFL23.6 MAb is directed against an erythroid cell surface antigen. It shows a well-defined reactivity with cells of the erythroid lineage at all stages of maturation in the peripheral blood, bone marrow, and fetal liver. Non-erythroid lineages are negative by flow cytometry. Although the exact identity of the antigen has yet to be determined, it has been shown to be distinct from Glycophorin A. This MAb is useful in the diagnosis of erythroleukemia, identification of bone marrow erythroid precursors, gating erythroid nucleated precursor cells from malignant cells in bone marrow specimens.

References

Note: References for this clone sold by other suppliers may be listed for expected applications.

Blood (1985) 66(3), 522-526. (Flow, IF)

Product attributes

Antibody number	#0105
Antibody reactivity (target)	Erythrocytes
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	SFL23.6
Isotype	IgG2b, kappa
Molecular weight	Not Known
Synonyms	RBC, Red blood cells, Erythroid Marker, Erythrocyte specific protein
Human gene symbol	Not Known
Entrez gene ID	Not Known
SwissProt	Not Known
Unigene	Not Known
Immunogen	Hepatocytes from a 20-22 week-old human fetus
Antibody target cellular localization	Plasma membrane
Species reactivity	Human
Expected antibody applications	Flow, surface (published for clone), IF (published for clone)
Antibody application notes	Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Flow cytometry: 0.5-1 ug/million cells, Immunofluorescence: 0.5-1 ug/mL, Optimal dilution for a specific application should be determined by user
Positive control	RBC, Placenta.
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	Red blood cells
Antibody research areas	Hematology