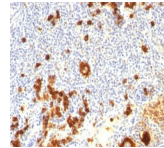


CD15 / FUT4 / Lewis-X Monoclonal Mouse Antibody (FUT4/815 + BRA-4F1)



Product Description

CD15 plays a role in mediating phagocytosis, bactericidal activity, and chemotaxis. It is present on >95% of granulocytes including neutrophils and eosinophils and to a lesser degree on monocytes. In addition, CD15 is expressed in Reed-Sternberg cells and some epithelial cells. CD15 antibody is very useful in the identification of Hodgkin's disease. CD15 is occasionally expressed in large cell lymphomas of both B and T phenotypes which otherwise have a quite distinct histological appearance. This antibody is available purified with BSA/azide at 200 ug/mL, or BSA/azide-free at 1 mg/mL.

Product attributes

Antibody number	#0562
Antibody reactivity (target)	CD15, FUT4, Lewis-x
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	FUT4/815 + BRA-4F1
Isotype	IgM
Molecular weight	~220 kDa (Glycoprotein)
Synonyms	3 Fucosyl N Acetyl Lactosamine; Alpha (1,3) Fucosyltransferase; Alpha 13 fucosyltransferase FucT; ELAM Ligand Fucosyltransferase; ELFT; FCT3A; Fuc-TIV; Fucosyltransferase 4 Alpha 1 3 Fucosyltransferase Myeloid Specific; Fucosyltransferase 4; Galactoside 3 L Fucosyltransferase; Lewis X; LeX; SSEA1; Stage Specific Embryonic Antigen 1
Human gene symbol	FUT4
Entrez gene ID	2526
SwissProt	P22083
Unigene	654379
Immunogen	Purified human neutrophils (FUT4/815); Myelomonocytic leukemia cells (BRA-4F1)
Verified antibody	IHC (FFPE) (verified)
Antibody target cellular localization	Plasma membrane, Perinuclear
Species reactivity	Human
Antibody application notes	Immunohistology (formalin), Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM Tris with 1 mM EDTA, pH 9.0, for 10-20 min followed by cooling at RT for 20 minutes, Immunofluorescence 0.5-1 ug/mL, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user
Positive control	U937 cells, Reed-Sternberg's cells in Hodgkin's lymphoma
Shipping condition	Room temperature
Storage Conditions	Store at 2 to 8 °C, Note: store BSA-free antibodies at -10 to -35 °C
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
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
Antibody research areas	Immunology
Cell/tissue expression	Granulocytes
Tumor expression	Leukemia/lymphoma