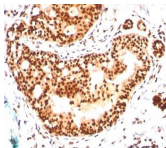


CD98 Monoclonal Mouse Antibody (IPO-T10)



Product Description

CD98 exists as a heterodimer containing a disulphide-linked glycosylated heavy chain and a non-glycosylated light chain. It is a member of the solute carrier family and encodes a cell surface, transmembrane protein. The protein exists as the heavy chain of a heterodimer, covalently bound through disulfide bonds to one of several possible light chains. The encoded transporter plays a role in regulation of intracellular calcium levels and transports L-type amino acids. Alternatively spliced transcript variants, encoding different isoforms, have been characterized. This antibody is available purified with BSA/azide at 200 ug/mL, or BSA/azide-free at 1 mg/mL.

References

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

- Int Immunol (2000) 12(7), 1051-1061. (Flow, functional studies)
- Immunology (1999) 96(1): 35-47. (functional studies)

Product attributes	
Antibody number	#0337
Antibody reactivity (target)	CD98
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	IPO-T10
Isotype	IgM, kappa
Molecular weight	Heterodimer (125 kDa) made of a disulfide-linked glycosylated heavy chain of approximately 80-90 kDa and a non-glycosylated light chain
Synonyms	4F2, 4F2HC, 4T2HC, CD98HC; MDU1; NACAE; Slc3a2; Solute Carrier Family 3 (activators of dibasic and neutral amino acid transport) Member 2, E16, LAT1, MPE16, hLAT1
Human gene symbol	SLC3A2
Entrez gene ID	6520
SwissProt	P08195
Unigene	502769
Immunogen	Molt 13 T cell line
Antibody target cellular localization	Plasma membrane
Expected antibody applications	Immunofluorescence (expected), Flow, surface (published for clone), Functional studies (published for clone)
Species reactivity	Human
Positive control	HL-60, Jurkat, MG63, HUT-78, K562, YT, U937, Hep-G2 cells, Tonsil
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
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	Transporters