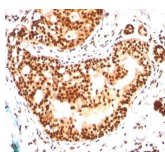


# CD98 Monoclonal Mouse Antibody (IPO-T10)



## 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.

**Catalog number key for antibody number 0337, Anti-CD98 (IPO-T10)**

## Product attributes

<b>Antibody number</b>	#0337
<b>Antibody reactivity (target)</b>	CD98
<b>Antibody type</b>	Primary
<b>Host species</b>	Mouse
<b>Clonality</b>	Monoclonal
<b>Clone</b>	IPO-T10
<b>Isotype</b>	IgM, kappa
<b>Molecular weight of antigen</b>	Heterodimer (125 kDa) made of a disulfide-linked glycosylated heavy chain of approximately 80-90 kDa and a non-glycosylated light chain
<b>Synonyms</b>	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
<b>Human gene symbol</b>	SLC3A2
<b>Entrez gene ID</b>	6520
<b>SwissProt</b>	P08195
<b>Unigene</b>	502769
<b>Immunogen</b>	Molt 13 T cell line
<b>Antibody target cellular localization</b>	Plasma membrane
<b>Expected antibody applications</b>	Immunofluorescence (expected), Flow, surface (published for clone), Functional studies (published for clone)
<b>Species reactivity</b>	Human
<b>Positive control</b>	HL-60, Jurkat, MG63, HUT-78, K562, YT, U937, Hep-G2 cells, Tonsil
<b>Shipping condition</b>	Room temperature
<b>Storage Conditions</b>	Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C
<b>Shelf life</b>	Guaranteed for at least 24 months from date of receipt when stored as recommended
<b>Regulatory status</b>	For research use only (RUO)
<b>Antibody/conjugate formulation</b>	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
<b>Antibody research areas</b>	Transporters
<b>Product origin</b>	Product may contain either bovine serum albumin (BSA) from bovine serum ( <i>Bos taurus</i> ), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel	Dye Features
BNC04	CF@405S	404/431	405	DAPI (microscopy), AF405	<a href="#">CF@405S Features</a>
BNC88	CF@488A	490/515	488	GFP, FITC	<a href="#">CF@488A Features</a>
BNC68	CF@568	562/583	532, 561	RFP, TRITC	<a href="#">CF@568 Features</a>
BNC94	CF@594	593/614	561	Texas Red@	<a href="#">CF@594 Features</a>
BNC40	CF@640R	642/662	633-640	Cy@5	<a href="#">CF@640R Features</a>
BNC47	CF@647	650/665	633-640	Cy@5	<a href="#">CF@647 Features</a>
BNC74	CF@740	742/767	633-685	775/50	<a href="#">CF@740 Features</a>
BNCB	Biotin	N/A	N/A	N/A	
BNUB	Purified	N/A	N/A	N/A	
BNUM	Purified, BSA-free	N/A	N/A	N/A	

Alexa Fluor, Pacific Blue, Pacific Orange, and Texas Red are trademarks or registered trademarks of Thermo Fisher Scientific; Cy is a registered trademark of Cytiva; IRDye, LI-COR, and Odyssey are registered trademarks of LI-COR Bioscience.

## 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)

This datasheet was generated on June 9, 2026 at 09:17:37 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/monoclonal-anti-cd98-ipo-t10/>