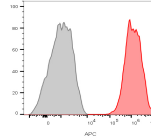


## CD81 / TAPA-1 Recombinant Monoclonal Rabbit Antibody (C81/2885R)

This antibody recognizes CD81, a member of a family of tetraspan transmembrane proteins.



### Product Description

This antibody recognizes a protein of 26 kDa, identified as CD81 (Workshop VI; Code CD81.1). CD81 has a very broad cellular distribution, being expressed on T- and B-lymphocytes, NK cells, thymocytes, eosinophils, fibroblasts, epithelial and endothelial cells. Neutrophils, erythrocytes and platelets are negative, while monocytes are variably positive. CD81 is a member of a family of tetraspanin transmembrane proteins, including CD9, CD37, CD53, CD63, and CD82. It associates with CD19, CD21, Leu 13, and integrins on cell membrane and is involved in signal transduction in B lymphocyte development and cell adhesion. CD81 also acts as a receptor for the envelope protein E2 of chronic hepatitis C virus. Antibodies to CD81 have anti-proliferative effects on different lymphoid cell lines, particularly those derived from large cell lymphomas. Primary antibodies are available purified, or with a selection of fluorescent CF® dyes and other labels. CF® dyes offer exceptional brightness and photostability. See the [CF® Dye Brochure](#) for more information. Note: Conjugates of blue fluorescent dyes like CF®405S and CF®405M are not recommended for detecting low abundance targets, because blue dyes have lower fluorescence and can give higher non-specific background than other dye colors. **Stock status:** Because Biotium offers a large number of antibody and conjugation options, primary antibody conjugates may be made to order. Typical lead times are up to one week for CF® dye and biotin conjugates, and up to 2-3 weeks for fluorescent protein and enzyme conjugates. Please email [order@biotium.com](mailto:order@biotium.com) to inquire about stock status and lead times before placing your order.

### Catalog number key for antibody number 2885, Anti-CD81 / TAPA-1 (C81/2885R)

Antibody #	prefix	Conjugation	Ex/Em	Concentration	Storage Buffer
BNC04	CF®405S	404/431 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide	
BNC05	CF®405M	408/452 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide	
BNC06	CF®405L	395/545 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide	
BNC88	CF®488A	490/515 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide	
BNC14	CF®514	516/548 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide	
BNC43	CF®543	541/560 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide	
BNC55	CF®555	555/565 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide	
BNC68	CF®568	562/583 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide	
BNC94	CF®594	593/614 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide	
BNC40	CF®640R	642/662 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide	
BNC47	CF®647	650/665 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide	
BNC60	CF®660C	667/685 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide	
BNC61	CF®660R	663/682 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide	
BNC80	CF®680	681/698 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide	
BNC81	CF®680R	680/701 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide	
BNC00	CF®700	695/720 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide	
BNC70	CF®770	770/797 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide	
BNCR	R-PE (PE)	496, 546, 565/578 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide	
BNCA	APC	650/660 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide	
BNCP	PerCP	482/677 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide	
BNCB	Biotin	N/A	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide	
BNCAP	Alkaline Phosphatase	N/A	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide	
BNCH	Horseradish Peroxidase	N/A	0.1 mg/mL	PBS, 0.05% BSA, no azide	
BNUB	Purified, with BSA	N/A	0.2 mg/mL	PBS, 0.05% BSA, 0.05% azide	
BNUM	Purified, BSA-free	N/A	1 mg/mL	PBS, no BSA, no azide	

### References

Kishimoto T et al. eds. Leukocyte Typing VI, p187-189, Garland Publishing, New York, 1997. |Takahashi S et al. J Immunol 1990, 145(7):2207-2213.

### Product attributes

Supplied As	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, BSA-free: 1 mg/mL in PBS without azide, Purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% azide
Antibody number	2885
Reactivity (target)	CD81, TAPA-1
Antibody type	Primary
Host species	Rabbit
Clonality	Recombinant Monoclonal
Clone	C81/2885R
Isotype	IgG
Molecular weight	26 kDa
Synonyms	CD81; Target of the antiproliferative antibody 1 (TAPA-1); Tetraspanin-28; Tspan-28
Human gene symbol	CD81
Entrez gene ID	975
SwissProt	P60033
Unigene	54457
Immunogen	B-Cell line derived from a Burkitt lymphoma
Cellular localization	Membrane/cell surface
Species reactivity	Human, Mouse, Rat
Applications	Immunohistology (formalin)
Application notes	Immunohistology (formalin): 2-4 ug/mL for 30 minutes at RT, Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM Tris with 1 mM EDTA pH 9.0 for 10-20 minutes followed by cooling at RT for 20 minutes, Optimal dilution for a specific application should be determined by user
Positive control	Ramos, Daudi, Raji, U266, HUT-78, HUT-102, Jurkat, HL-60, KG1a, A293, A431, MG63, and human lymphocytes. Tonsil or Liver.
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)
Research areas	Immunology