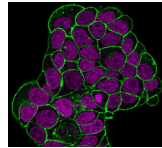


CD47 Monoclonal Mouse Antibody (IAP/964)

This antibody reacts with Ig domain of CD47 protein.



Product Description

This antibody reacts with Ig domain of CD47 protein. CD47, originally named integrin-associated protein (IAP), is a 50 kDa protein containing five membrane-spanning sequences and a short cytoplasmic tail. CD47 plays a role in both cell adhesion by acting as an adhesion receptor for THBS1 on platelets, and in the modulation of integrins. It is important in memory formation and synaptic plasticity in the hippocampus. CD47 may play a role in membrane transport and/or integrin dependent signal transduction. 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 to inquire about stock status and lead times before placing your order. **Catalog number key for antibody number 0964, Anti-CD47 (IAP/964)**

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
BNC88	CF®488A		490/515 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
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
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

Edris B et. al. Proc Natl Acad Sci USA 109:6656-61 (2012). | Willingham SB et. al. Proc Natl Acad Sci USA 109:6662-7 (2012)

Product attributes

Antibody number	0964
Reactivity (target)	CD47
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	IAP/964
Isotype	IgG1, kappa
Molecular weight	~50 kDa
Synonyms	Antigenic Surface Determinant Protein OA3; IAP; Integrin Associated Protein; Integrin Associated Signal Transducer; MER6; OA3; Protein MER6; Rh Related Antigen
Human gene symbol	CD47
Entrez gene ID	961
SwissProt	Q08722
Unigene	446414
Immunogen	Recombinant human IAP protein
Cellular localization	Cytoplasmic, Membrane/cell surface
Species reactivity	Human, Mouse
Applications	Immunofluorescence, Immunohistology (formalin), Flow cytometry
Application notes	Immunohistology formalin-fixed 0.5-1 ug/mL, Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 min followed by cooling at RT for 20 minutes, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Immunofluorescence 1-2 ug/mL, Optimal dilution for a specific application should be determined by user
Positive control	HUVEC, OVCA3 cells, Placenta, Brain and Ovarian Tumors
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	Cell adhesion, Hematology, Neuroscience, Signal transduction