Glowing products for science

CD63 Recombinant Monoclonal Mouse Antibody (2304.63) - Biotium Choice

CD63 (clone 2304.63) is a validated recombinant mouse monoclonal antibody that recognizes the human tetraspanin protein CD63. The antibody belongs to the Biotium Choice list of select antibodies that have been formulated and validated in-house for optimal performance.

afatum 8

Product Description

CD63 (clone 2304.63) is a recombinant mouse monoclonal antibody that recognizes the human tetraspanin protein CD63. This antibody belongs to the Biotium Choice list of select antibodies that have been formulated and validated in-house for optimal performance. This antibody has been validated in flow cytometry and immunofluorescence microscopy. The antibody is available unlabeled or conjugated to CF® Dyes.

The tetraspanins are integral membrane proteins expressed on cell surface and granular membranes of hematopoietic cells and are components of multi-molecular complexes with specific integrins. The tetraspanin CD63 is a lysosomal membrane glycoprotein that translocates to the plasma membrane after platelet activation. CD63 is expressed on activated platelets, monocytes and macrophages, and is weakly expressed on granulocytes, T cell and B cells. It is located on the basophilic granule membranes and on the plasma membranes of lymphocytes and granulocytes. CD63 is a member of the TM4 superfamily of leukocyte glycoproteins that includes CD9, CD37 and CD53, which contain four transmembrane regions. CD63 may play a role in phagocytic and intracellular lysosome-phagosome fusion events. CD63 deficiency is associated with Hermansky-Pudlak syndrome and is strongly expressed during the early stages of melanoma progression.

Tetraspanin proteins CD9, CD63, and CD81, are the most common proteins used as exosome markers. Tetraspanins are plasma membrane proteins with many proposed functions, including activation and sorting of other membrane proteins. They are also thought to play a role in the targeting of proteins to multivesicular bodies (MVBs) and exosomes. These tetraspanins are broadly expressed on many cell types and can therefore be detected on many types of exosomes, but their expression levels vary depending on the cell type of origin.

As an isotype control, we recommend <u>Isotype Control, Monoclonal Mouse IgG1 Kappa (IGG1/1331)</u> for this antibody.

Exosome Stains and Antibodies

Biotium also offers ExoBriteTM CD63 Flow Antibody and ExoBriteTM CD63 Western Antibody that are validated for optimal detection of isolated exosomes by flow cytometry or western blotting. Learn more about our products for exosome research.

Discover Sharper Signals and Unmatched Panel Flexibility with Biotium Choice Antibodies Powered by CF® Dyes and Astral Leap™ Tandem Dyes

Biotium Choice antibodies are carefully curated and validated in-house to offer exceptional signal-to-noise. Labeled with our advanced CF® Dyes and Astral LeapTM tandem dyes, they are our top-recommended antibodies for flow cytometry and other applications.

Biotium Choice Antibody Features

- Robust and validated clones against common targets
- Developed and optimized for flow cytometry and other applications
- Conjugated to bright, photostable <a>CF® <a>Dyes for superior signal and clarity
- Also available with <u>Astral Leap™ tandem dyes</u> for expanding multiplexing
- New antibody clones and dye conjugates continuously in development

View our full catalog of Biotium Choice antibodies

Call us: 800-304-5357 Email: btinfo@biotium.com

Product attributes

Antibody number #P013 SwissProt P08962 Biotium Choice Antibody Primary Antibody type Blotium Choice Flow Antibodies, Primary Clonality Recombinant Monoclonal Host species Mouse Clone 2304.63 Isotype IgG1, kappa Antibody reactivity (target) CD63 Synonyms Gp55; granulophysin; Lysosomal-associated membrane protein 3 (LAMP-3); Mast cell antigen AD1; melanoma 1 antigen; Melanoma-associated antigen MLA1; Melanoma-associated antigen; ME491; MLA1; MeA; Ocular melanoma-associated antigen; OMA81H; PTLGP40; Tetraspanin-30; TSPAN30 Species reactivity Baboon, Cynomolgus monkey, Human, Non-human primates Human gene symbol CD63 Entrez gene ID 967 Unigene 445570 Molecular weight 26 kDa (core protein); 30-60 kDa (glycosylated) Antibody target cellular localization Exosomes/EVs, Lysosomes, Plasma membrane, Membrane/vesicular, Multivesicular bodies Cell/tissue expression Exosomes, Platelats, Granulocytes, Lymphocytes, Monocytes/macrophages Verified antibody applications Positive control U87MG, SK-MEL-28, HL60, THP-1, NIH/3T3, or MCF-7 cells. Human melanoma, spleen or lymphoma tissue. Antibody research areas Exosomes/EVs Antibody research areas Exosomes/EVs Antibody research areas Exosomes/EVs Antibody research areas Exosomes/EVs Shelf life Guaranteed for at least 24 months from date of receipt when stored as recommended Volume per assay 5 uL/test Storage Conditions For research use only (RUO)					
Biotium Choice Antibody Antibody type Biotium Choice Flow Antibodies, Primary Clonality Recombinant Monoclonal Host species Mouse Clone 2304.63 Isotype IgG1, kappa Antibody reactivity (target) Synonyms Qp55; granulophysin; Lysosomal-associated membrane protein 3 (LAMP-3); Mast cell antigen AD1; melanoma 1 antigen; Melanoma-associated antigen MLA1; Melanoma-associated antigen MLA1; Melanoma-associated antigen MLA1; Melanoma-associated antigen; OMA81H; PTLGP40; Tetraspanin-30; TSPAN30 Species reactivity Baboon, Cynomolgus monkey, Human, Non-human primates Human gene symbol CD63 Entrez gene ID 967 Unigene 445570 Molecular weight Antibody target cellular localization Cell/tissue expression Exosomes/EVs, Lysosomes, Plasma membrane, Membrane/vesicular, Multivesicular bodies Cell/tissue expression Exosomes, Platelets, Granulocytes, Lymphocytes, Monocytes/macrophages Verified antibody applications Positive control U87MG, SK-MEL-28, HL60, THP-1, NIH/3T3, or MCF-7 cells. Human melanoma, spleen or lymphoma tissue. Antibody application notes Immunofluorescence: 1 ug/mL, Western Blot: 0.1 ug/mL, Flow cytometry: 0.5-1 ug/million cells Antibody/conjugate formulation Shelf life Guaranteed for at least 24 months from date of receipt when stored as recommended Volume per assay 5 uL/test Storage Conditions Note: store unconjugated formats at -10 °C to -35 °C, Store conjugates at 2 °C to 8 °C	Antibody number	#P013			
Antibody type Biotium Choice Flow Antibodies, Primary Clonality Recombinant Monoclonal Host species Mouse Clone 2304.63 Isotype IgG1, kappa Antibody reactivity (target) CD63 Synonyms gp55; granulophysin; Lysosomal-associated membrane protein 3 (LAMP-3); Mast cell antigen AD1; melanoma 1 antigen; Melanoma-associated antigen ML41; Melanoma-associated antigen; OMA81H; PTLGP40; Tetraspanin-30; TSPAN30 Species reactivity Baboon, Cynomolgus monkey, Human, Non-human primates Human gene symbol CD63 Entrez gene ID 967 Unigene 445570 Molecular weight 26 kDa (core protein); 30-60 kDa (glycosylated) Exosomes/EVs, Lysosomes, Plasma membrane, Membrane/vesicular, Multivesicular bodies Cell/tissue expression Exosomes, Platelets, Granulocytes, Lymphocytes, Monocytes/macrophages Verified antibody application Flow (surface) (verified), IF (verified) Positive control U87MG, SK-MEL-28, HL60, THP-1, NIH/3T3, or MCF-7 cells. Human melanoma, spleen or lymphoma tissue. Antibody application notes Exosomes/EVs Antibody/conjugate Conjugates: PBS/0.1% BSA/0.05% azide, Purified, BSA-free: 5 mg/mL in PBS without azide Shelf life Guaranteed for at least 24 months from date of receipt when stored as recommended Volume per assay 5 uL/test Storage Conditions Note: store unconjugated formats at -10 °C to -35 °C, Store conjugates at 2 °C to 8 °C	SwissProt	P08962			
Clonality Recombinant Monoclonal Host species Mouse Clone 2304.63 Isotype IgG1, kappa Antibody reactivity (target) Synonyms Gp55; granulophysin; Lysosomal-associated membrane protein 3 (LAMP-3); Mast cell antigen AD1; melanoma 1 antigen; Melanoma-associated antigen ML41; Melanoma-associated antigen ML41; Melanoma-associated antigen ML41; Melanoma-associated antigen ME491; MLA1; NGA; Ocular melanoma-associated antigen ME491; MLA1; NGA; Ocular melanoma-associated antigen; OMA81H; PTLGP40; Tetraspanin-30; TSPAN30 Species reactivity Baboon, Cynomolgus monkey, Human, Non-human primates Human gene symbol CD63 Entrez gene ID 967 Unigene 445570 Molecular weight 26 kDa (core protein); 30-60 kDa (glycosylated) Antibody target cellular localization Exosomes/EVs, Lysosomes, Plasma membrane, Membrane/vesicular, Multivesicular bodies Cell/tissue expression Exosomes, Platelets, Granulocytes, Lymphocytes, Monocytes/macrophages Verified antibody applications Positive control U87MG, SK-MEL-28, HL60, THP-1, NIH/3T3, or MCF-7 cells. Human melanoma, spleen or lymphoma tissue. Antibody application notes Immunofluorescence: 1 ug/mL, Western Blot: 0.1 ug/mL, Flow cytometry: 0.5-1 ug/million cells Antibody/conjugate Conjugates: PBS/0.1% BSA/0.05% azide, Purified, BSA-free: 5 mg/mt. in PBS without azide Shelf life Guaranteed for at least 24 months from date of receipt when stored as recommended Volume per assay 5 uL/test Storage Conditions Note: store unconjugated formats at -10 °C to -35 °C, Store conjugates at 2 °C to 8 °C	Biotium Choice Antibody	Primary			
Host species Mouse Clone 2304.63 Isotype IgG1, kappa Antibody reactivity (target) CD63 Synonyms	Antibody type	Biotium Choice Flow Antibodies, Primary			
Clone 2304.63 Isotype IgG1, kappa Antibody reactivity (target) CD63 Synonyms Gp55; granulophysin; Lysosomal-associated membrane protein 3 (LAMP-3); Mast cell antigen AD1; melanoma 1 antigen; Melanoma-associated antigen MLA1; Melanoma-associated antigen MLA1; Melanoma-associated antigen MLA1; Melanoma-associated antigen; OMA81H; PTLGP40; Tetraspanin-30; TSPAN30 Species reactivity Baboon, Cynomolgus monkey, Human, Non-human primates Human gene symbol CD63 Entrez gene ID 967 Unigene 445570 Molecular weight 26 kDa (core protein); 30-60 kDa (glycosylated) Antibody target cellular localization Exosomes/EVs, Lysosomes, Plasma membrane, Membrane/vesicular, Multivesicular bodies Cell/tissue expression Exosomes, Platelets, Granulocytes, Lymphocytes, Monocytes/macrophages Verified antibody applications Flow (surface) (verified), IF (verified) Positive control U87MG, SK-MEL-28, HL60, THP-1, NIH/3T3, or MCF-7 cells. Human melanoma, spleen or lymphoma tissue. Antibody application notes Immunofluorescence: 1 ug/mL, Western Blot: 0.1 ug/mL, Flow cytometry: 0.5-1 ug/million cells Antibody/conjugate Conjugates: PBS/0.1% BSA/0.05% azide, Purified, BSA-free: 5 mg/mt. in PBS without azide Shelf life Guaranteed for at least 24 months from date of receipt when stored as recommended Volume per assay 5 uL/test Storage Conditions Note: store unconjugated formats at -10 °C to -35 °C, Store conjugates at 2 °C to 8 °C	Clonality	Recombinant Monoclonal			
Isotype IgG1, kappa Antibody reactivity (target) CD63 Synonyms gp55; granulophysin; Lysosomal-associated membrane protein 3 (LAMP-3); Mast cell antigen AD1; melanoma 1 antigen; Melanoma-associated antigen MLA1; Melanoma-associated antigen MLA1; Melanoma-associated antigen WE491; MLA1; MGA; Coular melanoma-associated antigen; OMA81H; PTLGP40; Tetraspanin-30; TSPAN30 Species reactivity Baboon, Cynomolgus monkey, Human, Non-human primates Human gene symbol CD63 Entrez gene ID 967 Unigene 445570 Molecular weight 26 kDa (core protein); 30-60 kDa (glycosylated) Antibody target cellular Exosomes/EVs, Lysosomes, Plasma membrane, Membrane/vesicular, Multivesicular bodies Cell/tissue expression Exosomes, Platelets, Granulocytes, Lymphocytes, Monocytes/macrophages Verified antibody applications Positive control U87MG, SK-MEL-28, HL60, THP-1, NIH/3T3, or MCF-7 cells. Human melanoma, spleen or lymphoma tissue. Antibody application notes Exosomes/EVs Antibody/conjugate Conjugates: PBS/0.1% BSA/0.05% azide, Purified, BSA-free: 5 mg/mt. in PBS without azide Shelf life Guaranteed for at least 24 months from date of receipt when stored as recommended Volume per assay 5 uL/test Storage Conditions Note: store unconjugated formats at -10 °C to -35 °C, Store conjugates at 2 °C to 8 °C	Host species	Mouse			
Antibody reactivity (target) Synonyms Qp55; granulophysin; Lysosomal-associated membrane protein 3 (LAMP-3); Mast cell antigen AD1; melanoma 1 antigen; Melanoma-associated antigen MLA1; Melanoma-associated antigen; OMA81H; PTLGP40; Tetraspanin-30; TSPAN30 Species reactivity Baboon, Cynomolgus monkey, Human, Non-human primates Human gene symbol CD63 Entrez gene ID 967 Unigene 445570 Molecular weight Antibody target cellular localization Exosomes/EVs, Lysosomes, Plasma membrane, Membrane/vesicular, Multivesicular bodies Cell/tissue expression Exosomes/EVs, Lysosomes, Plasma membrane, Membrane/vesicular, Multivesicular bodies Verified antibody applications Pilow (surface) (verified), IF (verified) Werified antibody application notes U87MG, SK-MEL-28, HL60, THP-1, NIH/3T3, or MCF-7 cells. Human melanoma, spleen or lymphoma tissue. Antibody application notes Immunofluorescence: 1 ug/mL, Western Blot: 0.1 ug/mL, Flow cytometry: 0.5-1 ug/million cells Antibody/conjugate conjugates: PBS/0.1% BSA/0.05% azide, Purified, BSA-free: 5 mg/mL in PBS without azide Shelf life Guaranteed for at least 24 months from date of receipt when stored as recommended Volume per assay 5 uL/test Storage Conditions Note: store unconjugated formats at -10 °C to -35 °C, Store conjugates at 2 °C to 8 °C	Clone	2304.63			
Synonyms gp55; granulophysin; Lysosomal-associated membrane protein 3 (LAMP-3); Mast cell antigen AD1; melanoma 1 antigen; Melanoma-associated antigen MLA1; Melanoma-associated antigen MLA1; Melanoma-associated antigen WLA1; Melanoma-associated antigen WLA1; Melanoma-associated antigen; OMA81H; PTLGP40; Tetraspanin-30; TSPAN30 Species reactivity	Isotype	IgG1, kappa			
(LAMP-3): Mast cell antigen AD1; melanoma 1 antigen; Melanoma-associated antigen MLA1; Melanoma-associated antigen MLA1; Melanoma-associated antigen MEA91; MLA1; MGA; Ocular melanoma-associated antigen; OMA81H; PTLGP40; Tetraspanin-30; TSPAN30 Species reactivity Baboon, Cynomolgus monkey, Human, Non-human primates Human gene symbol CD63 Entrez gene ID 967 Unigene 445570 Molecular weight Antibody target cellular lexicalization Cell/tissue expression Exosomes/EVs, Lysosomes, Plasma membrane, Membrane/vesicular, Multivesicular bodies Cell/tissue expression Exosomes, Platelets, Granulocytes, Lymphocytes, Monocytes/macrophages Verified antibody applications Positive control U87MG, SK-MEL-28, HL60, THP-1, NIH/3T3, or MCF-7 cells. Human melanoma, spleen or lymphoma tissue. Antibody application notes Immunofluorescence: 1 ug/mL, Western Blot: 0.1 ug/mL, Flow cytometry: 0.5-1 ug/million cells Antibody/conjugate formulation Shelf life Guaranteed for at least 24 months from date of receipt when stored as recommended Volume per assay 5 uL/test Storage Conditions Note: store unconjugated formats at -10 °C to -35 °C, Store conjugates at 2 °C to 8 °C	Antibody reactivity (target)	CD63			
Human gene symbol CD63 Entrez gene ID 967 Unigene 445570 Molecular weight 26 kDa (core protein); 30-60 kDa (glycosylated) Antibody target cellular localization Exosomes, Plasma membrane, Membrane/vesicular, Multivesicular bodies Cell/tissue expression Exosomes, Platelets, Granulocytes, Lymphocytes, Monocytes/macrophages Verified antibody applications Positive control U87MG, SK-MEL-28, HL60, THP-1, NIH/3T3, or MCF-7 cells. Human melanoma, spleen or lymphoma tissue. Antibody application notes Immunofluorescence: 1 ug/mL, Western Blot: 0.1 ug/mL, Flow cytometry: 0.5-1 ug/million cells Antibody/conjugate formulation Exosomes/EVs Antibody/conjugate Gonjugates: PBS/0.1% BSA/0.05% azide, Purified, BSA-free: 5 mg/mL in PBS without azide Volume per assay 5 uL/test Storage Conditions Note: store unconjugated formats at -10 °C to -35 °C, Store conjugates at 2 °C to 8 °C	Synonyms	(LAMP-3); Mast cell antigen AD1; melanoma 1 antigen; Melanoma-associated antigen MLA1; Melanoma-associated antigen ME491; MLA1; NGA; Ocular melanoma-associated			
Entrez gene ID 967 Unigene 445570 Molecular weight 26 kDa (core protein); 30-60 kDa (glycosylated) Antibody target cellular lexosomes/EVs, Lysosomes, Plasma membrane, Membrane/vesicular, Multivesicular bodies Cell/tissue expression Exosomes, Platelets, Granulocytes, Lymphocytes, Monocytes/macrophages Verified antibody applications Positive control U87MG, SK-MEL-28, HL60, THP-1, NIH/3T3, or MCF-7 cells. Human melanoma, spleen or lymphoma tissue. Antibody application notes lmmunofluorescence: 1 ug/mL, Western Blot: 0.1 ug/mL, Flow cytometry: 0.5-1 ug/million cells Antibody/conjugate formulation complete formulation guaranteed for at least 24 months from date of receipt when stored as recommended Volume per assay 5 uL/test Storage Conditions	Species reactivity	Baboon, Cynomolgus monkey, Human, Non-human primates			
Unigene 445570 Molecular weight 26 kDa (core protein); 30-60 kDa (glycosylated) Antibody target cellular lexosomes/EVs, Lysosomes, Plasma membrane, Membrane/vesicular, Multivesicular bodies Cell/tissue expression Exosomes, Platelets, Granulocytes, Lymphocytes, Monocytes/macrophages Verified antibody applications Positive control U87MG, SK-MEL-28, HL60, THP-1, NIH/3T3, or MCF-7 cells. Human melanoma, spleen or lymphoma tissue. Antibody application notes Immunofluorescence: 1 ug/mL, Western Blot: 0.1 ug/mL, Flow cytometry: 0.5-1 ug/million cells Antibody/conjugate formulation Exosomes/EVs Antibody/conjugate Conjugates: PBS/0.1% BSA/0.05% azide, Purified, BSA-free: 5 mg/mL in PBS without azide Shelf life Guaranteed for at least 24 months from date of receipt when stored as recommended Volume per assay 5 uL/test Storage Conditions Note: store unconjugated formats at -10 °C to -35 °C, Store conjugates at 2 °C to 8 °C	Human gene symbol	CD63			
Molecular weight 26 kDa (core protein); 30-60 kDa (glycosylated) Antibody target cellular localization Exosomes/EVs, Lysosomes, Plasma membrane, Membrane/vesicular, Multivesicular bodies Cell/tissue expression Exosomes, Platelets, Granulocytes, Lymphocytes, Monocytes/macrophages Verified antibody applications Flow (surface) (verified), IF (verified) Brostitive control U87MG, SK-MEL-28, HL60, THP-1, NIH/3T3, or MCF-7 cells. Human melanoma, spleen or lymphoma tissue. Antibody application notes Immunofluorescence: 1 ug/mL, Western Blot: 0.1 ug/mL, Flow cytometry: 0.5-1 ug/million cells Antibody/conjugate formulation Conjugates PBS/0.1% BSA/0.05% azide, Purified, BSA-free: 5 mg/mL in PBS without azide Shelf life Guaranteed for at least 24 months from date of receipt when stored as recommended Volume per assay 5 uL/test Storage Conditions Note: store unconjugated formats at -10 °C to -35 °C, Store conjugates at 2 °C to 8 °C	Entrez gene ID	967			
Antibody target cellular localization Exosomes/EVs, Lysosomes, Plasma membrane, Membrane/vesicular, Multivesicular bodies Cell/tissue expression Exosomes, Platelets, Granulocytes, Lymphocytes, Monocytes/macrophages Verified antibody applications Positive control U87MG, SK-MEL-28, HL60, THP-1, NIH/3T3, or MCF-7 cells. Human melanoma, spleen or lymphoma tissue. Antibody application notes Immunofluorescence: 1 ug/mL, Western Blot: 0.1 ug/mL, Flow cytometry: 0.5-1 ug/million cells Antibody/conjugate Conjugates: PBS/0.1% BSA/0.05% azide, Purified, BSA-free: 5 mg/mL in PBS without azide Shelf life Guaranteed for at least 24 months from date of receipt when stored as recommended Volume per assay 5 uL/test Storage Conditions Note: store unconjugated formats at -10 °C to -35 °C, Store conjugates at 2 °C to 8 °C	Unigene	445570			
Nembrane/vesicular, Multivesicular bodies	Molecular weight	26 kDa (core protein); 30-60 kDa (glycosylated)			
Monocytes/macrophages Verified antibody applications Flow (surface) (verified), IF (verified) Positive control U87MG, SK-MEL-28, HL60, THP-1, NIH/3T3, or MCF-7 cells. Human melanoma, spleen or lymphoma tissue. Antibody application notes Immunofluorescence: 1 ug/mL, Western Blot: 0.1 ug/mL, Flow cytometry: 0.5-1 ug/million cells Antibody research areas Excsomes/EVs Antibody/conjugate formulation Conjugates: PBS/0.1% BSA/0.05% azide, Purified, BSA-free: 5 mg/mL in PBS without azide Shelf life Guaranteed for at least 24 months from date of receipt when stored as recommended Volume per assay 5 uL/test Note: store unconjugated formats at -10 °C to -35 °C, Store conjugates at 2 °C to 8 °C					
applications Positive control U87MG, SK-MEL-28, HL60, THP-1, NIH/3T3, or MCF-7 cells. Human melanoma, spleen or lymphoma tissue. Antibody application notes Immunofluorescence: 1 ug/mL, Western Blot: 0.1 ug/mL, Flow cytometry: 0.5-1 ug/mllilion cells Antibody research areas Exosomes/EVs Antibody/conjugate formulation Conjugates: PBS/0.1% BSA/0.05% azide, Purified, BSA-free: 5 mg/mL in PBS without azide Shelf life Guaranteed for at least 24 months from date of receipt when stored as recommended Volume per assay 5 uL/test Storage Conditions Note: store unconjugated formats at -10 °C to -35 °C, Store conjugates at 2 °C to 8 °C	Cell/tissue expression				
Human melanoma, spleen or lymphoma tissue. Antibody application notes Immunofluorescence: 1 ug/mL, Western Blot: 0.1 ug/mL, Flow cytometry: 0.5-1 ug/million cells Antibody research areas Exosomes/EVs Antibody/conjugate formulation Conjugates PBS/0.1% BSA/0.05% azide, Purified, BSA-free: 5 mg/mL in PBS without azide Shelf life Guaranteed for at least 24 months from date of receipt when stored as recommended Volume per assay 5 uL/test Storage Conditions Note: store unconjugated formats at -10 °C to -35 °C, Store conjugates at 2 °C to 8 °C		Flow (surface) (verified), IF (verified)			
cytometry: 0.5-1 ug/million cells Antibody research areas Exosomes/EVs Antibody/conjugate formulation Shelf life Guaranteed for at least 24 months from date of receipt when stored as recommended Volume per assay 5 uL/test Storage Conditions Note: store unconjugated formats at -10 °C to -35 °C, Store conjugates at 2 °C to 8 °C	Positive control				
Antibody/conjugate formulation Conjugates: PBS/0.1% BSA/0.05% azide, Purified, BSA-free: 5 mg/ml. in PBS without azide Shelf life Guaranteed for at least 24 months from date of receipt when stored as recommended Volume per assay 5 uL/test Storage Conditions Note: store unconjugated formats at -10 °C to -35 °C, Store conjugates at 2 °C to 8 °C	Antibody application notes				
formulation mg/mL in PBS without azide Shelf life Guaranteed for at least 24 months from date of receipt when stored as recommended Volume per assay 5 uL/test Storage Conditions Note: store unconjugated formats at -10 °C to -35 °C, Store conjugates at 2 °C to 8 °C	Antibody research areas	Exosomes/EVs			
stored as recommended Volume per assay 5 uL/test Storage Conditions Note: store unconjugated formats at -10 °C to -35 °C, Store conjugates at 2 °C to 8 °C					
Storage Conditions Note: store unconjugated formats at -10 °C to -35 °C, Store conjugates at 2 °C to 8 °C	Shelf life				
conjugates at 2 °C to 8 °C	Volume per assay	5 uL/test			
Regulatory status For research use only (RUO)	Storage Conditions				
	Regulatory status	For research use only (RUO)			
Product origin Product may contain either bovine serum albumin (BSA) from bovine serum (Bos taurus), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.	Product origin	bovine serum (Bos taurus), or recombinant BSA produced in			

Conjugation	Ex/Em	Conc.	Size	Catalog No.	Dye Features
<u>CF®405M</u> 416/452 nm	416/452 nm	100 ug/mL	25 tests (125 uL)	P013-405M-125	CF®405M Features
			100 tests (500 uL)	P013-405M-500	
<u>CF®488A</u> 490/516 n	490/516 nm	100 ug/mL	25 tests (125 uL)	P013-488A-125	CF®488A Features
			100 tests (500 uL)	P013-488A-500	
<u>CF® 568</u>	562/584 nm	100 ug/mL	25 tests (125 uL)	P013-568-125	CF®568 Features
			100 tests (500 uL)	P013-568-500	
CF® 583R	585/609 nm	100 ug/mL	25 tests (125 uL)	P013-583R-125	CF®583R Features
			100 tests (500 uL)	P013-583R-500	
CF® 647Plus	652/668 nm	100 ug/mL	25 tests (125 uL)	P013-647P-125	CF®647Plus Features
			100 tests (500 uL)	P013-647P-500	
<u>CF®700</u>	696/721 nm	100 ug/mL	25 tests (125 uL)	P013-700-125	CF®700 Features
			100 tests (500 uL)	P013-700-500	
R-PE 496, 546, 565/576 nm	496, 546, 565/576 nm	100 ug/mL	25 tests (125 uL)	P013-RPE-125	
			100 tests (500 uL)	P013-RPE-500	
APC 651	651/660 nm	100 ug/mL	25 tests (125 uL)	P013-APC-125	
			100 tests (500 uL)	P013-APC-500	
Purified (BSA-free)	N/A	5 mg/mL	1 mg	P013-1MG	
			5 mg	P013-5MG	

This datasheet was generated on November 4, 2025 at 01:10:18 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/cd63-recombinant-monoclonal-mouse-antibody-230463-biotium-choice/