



Glowing products for science

## CD16 Mouse Monoclonal Antibody (3G8) - Biotium Choice

CD16 (clone 3G8) is a validated mouse monoclonal antibody that recognizes CD16 (also known as FC gamma receptor III), the low-affnity receptor for IgG. The antibody belongs to the Biotium Choice list of select antibodies that have been validated and optimized in-house for optimal performance.



### Product Description

CD16 (clone 3G8) is a mouse monoclonal antibody that recognizes IgG receptor CD16, also known as FC gamma receptor III. This antibody belongs to the Biotium Choice list of select antibodies that have been validated and optimized in-house for optimal performance. The antibody is available conjugated to CF® Dyes, Astral™ tandem dyes, R-PE and PerCP. They are supplied in PBS, 0.1% BSA, 0.05% azide.

This antibody recognizes CD16 (Fc gamma receptor III), the low-affinity receptor for IgG with an apparent molecular weight of 50-80 kDa. CD16 is represented by two similar genes, CD16A (Fc gamma receptor IIIA), which exists as a hetero-oligomeric polypeptide-anchored form in macrophages and NK cells and CD16B (Fc gamma receptor IIIB), which exist as a monomeric GPI-anchored form in neutrophils.

Furthermore, there are two known polymorphisms of CD16B, NA-1 and NA-2. Individuals homozygous for NA-2 show a lower phagocytic capacity compared with NA-1. CD16 binds IgG in the form of immune complexes and shows preferential binding of IgG1 and IgG3 isotypes and minimal binding of IgG2 and IgG4. Upon IgG binding, both CD16 isoforms initiate signal transduction cascades that lead to a variety of responses including antibody-dependent cell-mediated cytotoxicity (ADCC), phagocytosis, degranulation and proliferation.

As an isotype control, we recommend [Isotype Control, Monoclonal Mouse IgG1 Kappa \(IGG1/1331\)](#) for this antibody.

### 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 Leap™ 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 [CF® Dyes](#) for superior signal and clarity
- Also available with [Astral Leap™ tandem dyes](#) 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](tel:800-304-5357)

#### Product attributes

Antibody number	P011
Antibody reactivity (target)	CD16
Biotium Choice Antibody	Primary
Antibody type	Biotium Choice Flow Antibodies, Primary
Host species	Mouse
Clonality	Monoclonal
Clone	3G8
Isotype	IgG1, kappa
Molecular weight	29 kDa (predicted); 50-65 kDa (observed)
Synonyms	FCG3A, Fc gamma receptor 3
Human gene symbol	CD16
Entrez gene ID	2214
SwissProt	P08637
Unigene	372679
Immunogen	Human
Antibody target cellular localization	Plasma membrane
Verified antibody applications	Flow (surface) (verified)
Species reactivity	Human
Antibody application notes	Recommended concentration for flow cytometry: 5 uL per million cells/0.1 mL
Positive control	Human peripheral blood mononuclear cells (PBMCs)
Shipping condition	Room temperature
Storage Conditions	Protect fluorescent conjugates from light
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Regulatory status	For research use only (RUO)
Volume per assay	5 uL/test
Antibody/conjugate formulation	Conjugates: PBS/0.1% BSA/0.05% azide
Antibody research areas	Immunology
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.
Cell/tissue expression	Lymphocytes

## CD16 Monoclonal Mouse Antibody (3G8) - Biotium Choice

Conjugation	Ex/Em	Conc.	Size	Catalog No.	Dye Features
<a href="#">CF@405M</a> 100 tests (500 uL)	416/452 nm <a href="#">P011-405M-500</a>	100 ug/mL	25 tests (125 uL)	<a href="#">P011-405M-125</a>	<a href="#">CF@405M Features</a>
<a href="#">CF@488A</a> 100 tests (500 uL)	490/516 nm <a href="#">P011-488A-500</a>	50 ug/mL	25 tests (125 uL)	<a href="#">P011-488A-125</a>	<a href="#">CF@488A Features</a>
<a href="#">CF@568</a> 100 tests (500 uL)	562/584 nm <a href="#">P011-568-500</a>	50 ug/mL	25 tests (125 uL)	<a href="#">P011-568-125</a>	<a href="#">CF@568 Features</a>
<a href="#">RPE-Astral™616</a> 100 tests (500 uL)	496, 546, 566/617 nm <a href="#">P011-R616-500</a>	200 ug/mL	25 tests (125 uL)	<a href="#">P011-R616-125</a>	<a href="#">RPE-Astral™616 Features</a>
<a href="#">CF@647Plus</a> 100 tests (500 uL)	652/668 nm <a href="#">P011-647P-500</a>	25 ug/mL	25 tests (125 uL)	<a href="#">P011-647P-125</a>	
<a href="#">CF@700</a> 100 tests (500 uL)	696/721 nm <a href="#">P011-700-500</a>	100 ug/mL	25 tests (125 uL)	<a href="#">P011-700-125</a>	<a href="#">CF@700 Features</a>
<a href="#">APC-Astral™813</a> 100 tests (500 uL)	633, 638/813 nm <a href="#">P011-A813-500</a>	100 ug/mL	25 tests (125 uL)	<a href="#">P011-A813-125</a>	<a href="#">APC-Astral™813 Features</a>
<a href="#">R-PE</a> 100 tests (500 uL)	496, 546, 565/576 nm <a href="#">P011-RPE-500</a>	100 ug/mL	25 tests (125 uL)	<a href="#">P011-RPE-125</a>	
<a href="#">PerCP</a> 100 tests (500 uL)	482/675 nm <a href="#">P011-PCP-500</a>	100 ug/mL	25 tests (125 uL)	<a href="#">P011-PCP-125</a>	

This datasheet was generated on December 7, 2025 at 03:25:06 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/cd16-monoclonal-mouse-antibody-3g8/>