

Qdot[®] Mouse IgG1 Isotype Control Conjugate

Product	Form	Volume	Qdot [®]	Tests	Peak		Recommended
			Nanocrystal Conc.		Excitation (nm)	Emission (nm)	
Q10073	Qdot [®] 605	0.1 ml	1 μM	100 min.	405 (488)*	605	605/20
Q10297	Qdot [®] 655	0.1 mL	1 μM	100 min	405 (488)*	655	655/20
Q10298	Qdot [®] 800	0.1 mL	1 μM	100 min	405 (488)*	800	780/60

*Qdot[®] nanocrystals excite optimally in the UV to 405 nm range, but can also be excited with wavelengths shorter than their emission maximum, such as with a 488 nm laser.

Product Description

Mouse IgG1 isotype control

Lot No: See label **Expiration:** See label

Buffer: 50 mM borate, 1 M betaine, pH 8.3

Preservative: 0.05% sodium azide. Sodium azide is an extremely toxic and dangerous compound particularly when combined with acids or metals. Properly dispose of solutions containing sodium azide.

Storage and Handling

Store reagents at 2–8°C. **Do not freeze.** Because Qdot[®] nanocrystals are conjugated to biological materials, some loss of activity may be observed with prolonged storage.

Qdot[®] nanocrystals are photostable, and do not need to be protected from light. However, if using Qdot[®] conjugates in combination with conventional fluorochrome conjugated antibodies, minimize light exposure during handling, incubation with cells, and prior to analysis. We recommend analysis of cells within 18 hours of staining. If dilute reagent is used, dilute only the quantity of reagent to be used within one day.

The Qdot[®] conjugate contains cadmium and selenium in an inorganic crystalline form. Dispose of the material in compliance with all applicable local, state, and federal regulations for disposal of these classes of material. For more information on the composition of these materials, consult the Material Safety Data Sheet.

Qdot[®] Nanocrystals

Qdot[®] nanocrystals are nanometer-scale atom clusters of semiconductor material which exhibit narrow and symmetrical emission bandwidths with very long Stokes shifts. The nanocrystals can be excited with any wavelength below their emission maximum, but are best excited by UV or violet light. Qdot[®] nanocrystals demonstrate an intrinsic brightness and photostability that can be many times greater than observed with other classes of fluorophores. These advantages make Qdot[®] nanocrystals powerful tools for antibody staining.^{1,2}

Product Use

Staining: Stain cells in any standard staining buffer, such as phosphate-buffered saline (PBS) with 1% bovine serum albumin (BSA). Use the amount of isotype control necessary to match the amount of Qdot[®] primary antibody conjugate used: 1 μL of Qdot[®] antibody conjugate per 1 × 10⁶ cells in a 100 μL staining volume represents a 10 nM final concentration of Qdot[®] nanocrystals. Qdot[®] nanocrystal conjugates may be mixed with other antibodies, but use diluted conjugates on the day of dilution. Because conditions may vary, you may need to determine the optimal amount of antibody used for each application. Qdot[®] nanocrystal conjugates can be used for surface staining applications with most conventional sample preparation reagents, such as Caltag[®] Cal-Lyse[™] and Caltag[®] FIX & PERM[®] reagents, with minimal affect on fluorescence. We have observed that some batches of BD FACS[™] Lysing Solution have been found to interfere with Qdot[®] nanocrystal fluorescence.

Instrument setup: Qdot[®] nanocrystals are excited optimally with UV or 405 nm light, although excitation can be obtained with any wavelength below the emission wavelength of a given nanocrystal. Make sure the cytometer has an appropriate emission filter for the Qdot[®] nanocrystal being used. The table above has filter recommendations; alternate filters can be used as long as they capture the emission maximum, but filter width impacts spectral overlap corrections. **Note:** Qdot[®] nanocrystals can be used on cytometers that do not have UV or violet excitation sources as long as they have appropriate emission filters. Be sure to check for Qdot[®] nanocrystal emission in any channel that can capture nanocrystal emission off of other lasers on the cytometer.

Product Quality Control

Each lot has been tested by flow cytometry using human peripheral blood leukocytes (PBLs). This testing was performed using 2 μL of antibody per 1 × 10⁶ cells in a 100 μL staining volume (20 nM final concentration of Qdot[®] nanocrystals). Qdot[®] nanocrystal concentration is assigned based on optical density. See reverse for representative flow cytometry data.

References

1. Telford, W. G. 2004. *Cytometry Part A* 61A:9.
2. Perfetto, S. P., P. K. Chattopadhyah, and M. Roederer. 2004. *Nature Reviews - Immunology* 4: 648.

Related Products

Catalog no.	Product Name	Unit Size
GAS-010	Cal-Lyse [™] Whole Blood Lysing Solution	25 mL
GAS-010S-100	Cal-Lyse [™] Whole Blood Lysing Solution	100 mL
HYL-250	High-Yield Lyse Fixative	500 mL
GAS001S-5	FIX & PERM [®] Reagent A (Individual)	5 mL
GAS001S-100	FIX & PERM [®] Reagent A (Bulk)	100 mL
GAS002S-5	FIX & PERM [®] Reagent B (Individual)	5 mL
GAS002S-100	FIX & PERM [®] Reagent B (Bulk)	100 mL
GAS-003	FIX & PERM [®] Reagents	50 tests
GAS-004	FIX & PERM [®] Reagents	200 tests

Limited Use Label License No. 223: Labeling and Detection Technology

The manufacture, use, sale or import of this product may be subject to one or more patents or pending applications owned or licensed by Invitrogen Corporation. The purchase of this product conveys to the buyer the non-transferable right to use the purchased amount of the product and components of the product in research conducted by the buyer (whether the buyer is an academic or for-profit entity) in a manner consistent with the accompanying product literature. The buyer cannot sell or otherwise transfer (a) this product (b) its components or (c) materials made using this product or its components to a third party or otherwise use this product or its components or materials made using this product or its components for Commercial Purposes. The buyer may transfer information or materials made through the use of this product to a scientific collaborator, provided that such transfer is not for any Commercial Purpose, and that such collaborator agrees in writing (a) to not transfer such materials to any third party, and (b) to use such transferred materials and/or information solely for research and not for Commercial Purposes. Commercial Purposes means any activity by a party for consideration and may include, but is not limited to: (1) use of the product or its components in manufacturing; (2) use of the product or its components to provide a service, information, or data; (3) use of the product or its components for therapeutic, diagnostic or prophylactic purposes; or (4) resale of the product or its components, whether or not such product or its components are resold for use in research. For products that are subject to multiple limited use label licenses, the most restrictive terms apply. Invitrogen Corporation will not assert a claim against the buyer of infringement of patents that are owned or controlled by Invitrogen Corporation and/or Molecular Probes, Inc. which cover this product based upon the manufacture, use or sale of a therapeutic, clinical diagnostic, vaccine or prophylactic product developed in research by the buyer in which this product or its components was employed, provided that neither this product nor any of its components was used in the manufacture of such product. If the purchaser is not willing to accept the limitations of this limited use statement, Invitrogen is willing to accept return of the product with a full refund. For information on purchasing a license to this product for purposes other than research, contact Molecular Probes, Inc., Business Development, 29851 Willow Creek Road, Eugene, OR 97402, Tel: (541) 465-8300. Fax: (541) 335-0354.

Qdot, Caltag, FIX & PERM, and Cal-Lyse are trademarks of Invitrogen Corporation. FACS is a trademark of BD Biosciences.

Copyright 2007–2009, Molecular Probes, Inc. All rights reserved. This information is subject to change without notice.

For research use only.

Invitrogen Corporation • 5791 Van Allen Way • Carlsbad • CA 92008 • Tel: 800.955.6288 • www.invitrogen.com • E-mail: tech_support@invitrogen.com