Revised: 15–September–2005

Long-Wavelength Calcium Indicators

Calcium Green™ • Oregon Green® 488 BAPTA • Calcium Yellow™ • Calcium Orange™ • Calcium Crimson™

Introduction

The long-wavelength calcium indicators Calcium Green[™], Oregon Green[®] 488 BAPTA, Calcium Yellow[™], Calcium Orange[™], and Calcium Crimson[™] are visible light-excitable probes derived from fluorescein, Oregon Green® 488 dye, lucifer yellow, tetramethylrhodamine, and Texas Red®, respectively. Molecular Probes prepares these indicators in the water-soluble potassium salt form, cell-permeant acetoxymethyl (AM) ester form, and in some cases as compartmentalization-resistant dextran conjugates (Table 1). Upon binding to calcium, these indicators exhibit an increase in fluorescence emission intensity with little shift in wavelength (Figure 1). The spectral characteristics of the long-wavelength indicators have three major advantages: 1) Their emissions are in regions of the spectrum where cellular autofluorescence and scattering backgrounds are often less of a problem; 2) The energy of the excitation light is low, reducing the potential for cellular photodamage; and 3) The wavelengths required for optimal excitation are compatible with those produced by laser-based instrumentation, such as confocal laser scanning microscopes. Long-wavelength Ca2+ indicators also provide increased options for simultaneous measurements of other physiological parameters, such as pH and membrane potential,^{1,2} and for experiments involving photoactivatable "caged" probes.³

Although Calcium GreenTM-1 is structurally similar to fluo-3, it is more fluorescent at low calcium concentrations, facilitating the determination of baseline Ca²⁺ levels and increasing the visibility of resting cells (Figure 1A). Calcium GreenTM-1 and Oregon Green[®] 488 BAPTA-1 are currently the indicators of choice for multiphoton excitation imaging of calcium dynamics in brain slice preparations⁴ and intact live brains.⁵

In contrast to Calcium GreenTM-1, Calcium GreenTM-2 has two fluorescent reporter groups, which are believed to quench one another in the absence of calcium. Calcium GreenTM-2 undergoes a much larger increase in fluorescence emission upon calcium binding than does Calcium GreenTM-1 and its lower affinity for calcium makes it particularly suited to measuring relatively high spikes of calcium, up to 25 μ M (Table 1 and Figure 1B).

Oregon Green[®] 488 BAPTA-1 and -2 have generally similar properties to Calcium Green[™]-1 and -2, respectively. The major difference is that both the fluorescence excitation and emission maxima of the Oregon Green[®] 488 BAPTA indicators are shorter by ~10 nm (Table 1), resulting in more efficient 488 nm excitation by argon-ion lasers. The extinction coefficient for Oregon Green[®] 488 BAPTA-1 absorption at 488 nm is ~93% of its maximum value, whereas for Calcium Green[™]-1 it is only about 45% of maximum.

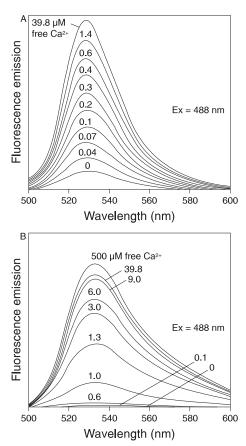


Figure 1. Emission spectra of A) Calcium Green[™]-1 and B) Calcium Green[™]-2.

Cell-impermeant Calcium YellowTM has an exceptionally large Stokes shift, with excitation/emission maxima of ~435/550 nm. Like Calcium GreenTM-1, Calcium YellowTM is fluorescent at low Ca²⁺ concentrations. This indicator has a high affinity for Ca²⁺ (K_d = 185 nM) and exhibits an approximately 19-fold increase in fluorescence intensity upon binding calcium in buffer solutions. Calcium YellowTM is compatible with standard filter sets suitable for lucifer yellow.

Calcium OrangeTM and Calcium CrimsonTM are spectrally similar to tetramethylrhodamine and Texas Red[®] (Table 1). The long-wavelength spectral characteristics of these indicators allow them to be used in combination with fluorescein and ultravioletexcitable dyes.⁶ The wavelength separation between the excitation spectrum of Calcium OrangeTM and the action spectrum of genetically-altered rhodopsin pigments has enabled researchers to detect and image light-induced Ca²⁺ transients in *Drosophila* photoreceptors.⁷

Table 1. Spectral characteristics of long-wavelength calcium indicators.

-		0 0				F _{Ca} / F _{free} ** K _d ††		
Indicator	Salt *	AM †	Dextran ‡	Ex/Em § (nm)	F _{Ca} /F _{free} **	K _d ††		
Calcium Green™-1	C3010	C3011, C3012	C3713, C3714, C6765, C6766	506/531	~14	190 nM		
Calcium Green™-2	C3730	C3732		503/536	~100	550 nM		
Calcium Green™-5N	C3737	C3739		506/532	~38	14 µM		
Oregon Green [®] 488 BAPTA-1	06806	06807	06798	494/523	~14	170 nM		
Oregon Green [®] 488 BAPTA-2	06808	06809		494/523	~100	580 nM		
Oregon Green [®] 488 BAPTA-6F	023990	023991		494/524	~14	3 µM		
Oregon Green [®] 488 BAPTA-5N	06812	06813		494/521	~44	20 µM		
Calcium Yellow™	C36202			435/550	~19	185 nM		
Calcium Orange™	C3013	C3015		549/576	~3	185 nM		
Calcium Crimson™		C3018		590/615	~2.5	185 nM		

* Catalog numbers for water-soluble potassium salt forms. † Catalog numbers for cell-permeant acetoxymethyl (AM) esters. ‡ Catalog numbers for dextran conjugates. § Excitation and Emission maxima for Ca²⁺-bound indicator. ** Fluorescence intensity of the Ca²⁺-bound indicator relative to its Ca²⁺-free form. All these indicators exhibit fluorescence enhancement on binding Ca²⁺ with essentially no change in excitation or emission wavelengths. †† Dissociation constant for Ca²⁺ determined *in vitro* at 22°C in 100 mM KCl, 10 mM MOPS, pH 7.2. K_d values depend on temperature, ionic strength, pH, presence of other ions (e.g. Mg²⁺) and other factors. Values reported are determined using the acid form of the indicator. K_d values for dextran conjugates may vary between production lots; lot-specific values are printed on the vial in most cases.

Calcium GreenTM-5N, Oregon Green[®] 488 BAPTA-5N, and Oregon Green[®] 488 BAPTA-6F have the lowest Ca²⁺ binding affinities in this series of indicators (Table 1) and are not expected to respond to normal changes in intracellular magnesium. These low-affinity indicators can be used to detect high-amplitude calcium transients, such as those evoked by neurotransmitters ⁸ and IP₃.⁹ The high Ca²⁺ dissociation rates of these low-affinity indicators are advantageous for rapid kinetic tracking of Ca²⁺ release.¹⁰⁻¹²

High molecular weight dextran conjugates can be loaded into cells using our Influx[™] pinocytic cell-loading reagent (I14402), by direct injection via a recording electrode, by electroporation, or by simply delivering the conjugate just outside cell bodies by pressure or electrophoretic application, which leads to spontaneous uptake.^{13, 14} The dextran conjugates show a dramatic reduction in both leakage and compartmentalization—a particular problem for the acidic forms of these probes in plant cells.¹⁵ We offer Calcium Green[™]-1 conjugated to 3000, 10,000, 70,000, and 500,000 MW dextrans. Our 500,000 MW dextran conjugate of Calcium Green[™]-1 should be effectively excluded from cellular organelles, including the nucleus.

Materials

These products are provided as lyophilized solids and should be stored at $\leq -20^{\circ}$ C, desiccated, and protected from light until use. Allow products to warm to room temperature before opening.

The salts may be reconstituted in aqueous buffers or distilled water; store aqueous stock solutions at $\leq -20^{\circ}$ C and protected from light. The salts are stable for at least six months in most physiological buffers.

AM esters are susceptible to hydrolysis, particularly in solution. They should be reconstituted just before use in highquality, anhydrous dimethylsulfoxide (DMSO). DMSO stock solutions may be stored desiccated at \leq -20°C and protected from light. Under these conditions, AM esters should be stable for several months. To prevent repeated exposure to moisture, AM esters are supplied in 500 µg units prealiquoted into ten vials of 50 µg each so that the entire contents do not have to be dissolved at the same time. Dilute aqueous solutions of the AM esters for cell loading should be used on the same day that they are prepared.

Application

The following AM ester loading protocol is provided as an introductory guide only. AM ester loading protocols for a variety of cell types and vesicle preparations can be obtained from the literature.¹⁶

Suggested AM Ester Loading Protocol

1. Prepare a 2 to 5 mM stock solution of the AM ester in high-quality, anhydrous DMSO, as described in *Materials*; the molecular weight (MW) is indicated on the product label. The nonionic detergent Pluronic[®] F-127 is sometimes used to increase the aqueous solubility of AM esters. For the convenience of our customers, Molecular Probes offers Pluronic F-127 in three forms: 1 mL of a 20% (w/v) solution in DMSO (P3000MP), 30 mL of a sterile 10% (w/v) solution in water (P6866), and 2 g of the solid (P6867). A 20% Pluronic F-127 solution can be used in place of DMSO to prepare solutions of these calcium indicators; however, long-term storage of AM esters in the presence of Pluronic F-127 is not recommended.

2. On the day of the experiment, either dissolve the lyophilized solid in DMSO or thaw an aliquot of the indicator stock solution to room temperature. Prepare a working solution of 1 to 10 μ M in the buffer of choice; the exact concentration of indicator required for cell loading must be determined empirically. To avoid calcium buffering, toxicity and other artifacts of overloading, one should generally use the lowest probe concentration that yields sufficient signal.

3. Incubate cells with the AM ester for 20 minutes to one hour at or below room temperature. Adherent cultures do not need to be lifted for loading. Some investigators report that decreasing the loading temperature reduces indicator compartmentalization. The zwitterionic Calcium Crimson[™] has been found to be particularly susceptible to this problem.

4. Wash cells to remove excess probe that either has not been loaded or may be noncovalently associated with the membrane. This step is important when labeling with Calcium Orange[™] and Calcium Crimson[™], which are fluorescent prior to ester cleavage.

Response Calibration

The excitation (λ_{EX}) and emission (λ_{EM}) wavelengths and calcium responses of these long-wavelength calcium indicators are shown in Table 1. The following bandpass filter sets are recommended for fluorescence microscopy applications:

- Oregon Green[®] 488 BAPTA indicators: Omega sets XF100 or XF23; Chroma sets 41001 or 31001
- Calcium Green[™] indicators: Omega sets XF104 or XF23; Chroma sets 41028 or 31001
- Calcium Yellow[™] indicators: Omega sets XF14-2, XF15, XF77-2; Chroma sets 31010 or 31038
- Calcium Orange[™] indicators: Omega sets XF108 or XF32; Chroma sets 41002 or 31002
- Calcium Crimson[™] indicators: Omega sets XF102 or XF43; Chroma sets 41004 or 31004

Omega[®] filters are supplied by Omega Optical Inc. (www.omegafilters.com). Chroma filters are supplied by Chroma Technology Corp. (www.chroma.com). Calibration is typically accomplished using the acid form of the indicator in solutions of precisely known free calcium concentrations, such as those generated using Molecular Probes' Calcium Calibration Buffer Kits (C3008MP, C3009, C3723) and Calcium Calibration Standard Kit (C6775). The Calcium Calibration Buffer Kits, which contain CaEGTA solutions that have been precisely prepared using Roger Tsien's "pH-metric" method,¹⁷ can be used to obtain calibration curves and K_ds for high- to moderate-affinity calcium indicators. In contrast, our Calcium Calibration Standard Kit provides 11 solutions, containing from zero to 1 mM calcium, for calibrating our low-affinity calcium indicators such as Calcium GreenTM-5N and Oregon Green[®] 488 BAPTA-5N.

To determine either the free calcium concentration of a solution or the K_d of a single-wavelength calcium indicator, the following equation is used:

$$\left[Ca^{2+}\right]_{\text{free}} = K_{d} \left[\frac{F - F_{\min}}{F_{\max} - F}\right]$$

where F is the fluorescence of the indicator at experimental calcium levels, F_{min} is the fluorescence in the absence of calcium and F_{max} is the fluorescence of the calcium-saturated probe. The dissociation constant (K_d) is a measure of the affinity of the probe for calcium.

The Ca²⁺ -binding and spectroscopic properties of fluorescent indicators vary quite significantly in cellular environments compared to calibration solutions. *In situ* response calibrations of intracellular indicators typically yield K_d values significantly higher than *in vitro* determinations.¹⁸ *In situ* calibrations are performed by exposing loaded cells to controlled Ca²⁺ buffers in the presence of ionophores such as A-23187 (A1493), 4-bromo A-23187 (B1494) and ionomycin (I24222).¹⁹ Alternatively, cell permeabilization agents such as digitonin or Triton[®] X-100 can be used to expose the indicator to the controlled Ca²⁺ levels of the extracellular medium.¹⁶ Calcium GreenTM-1, Oregon Green[®] 488 BAPTA-1, Calcium OrangeTM, and Calcium CrimsonTM are quite fluorescent at low free calcium concentrations, making visualization of resting cells and determination of F_{min} values easier than with fluo-3.

References

Biophys J 68, 739 (1995);
Cell Calcium 19, 337 (1996);
Neuron 15, 755 (1995);
Methods 18, 231 (1999);
Nature 385, 161 (1997);
Methods 18, 204 (1999);
Neuron 13, 837 (1994);
Neurosci Lett 163, 149 (1993);
J Physiol 476, 17 (1994);
Biophys J 70, 222 (1996);
Biophys J 70, 896 (1996);
Biophys J 70, 1006 (1996);
Real-Time Imaging 8, 175 (2002);
Neuron 27, 25 (2000);
J Microscopy 166, 57 (1992);
Cell Biology: A Laboratory Handbook, Second Edition, Volume 3, Celis J.E., Ed., Academic Press (1998) pp. 363-374;
Methods Enzymol 172, 230 (1989);
Cell Calcium 28, 213 (2000);
Cell Calcium 21, 233 (1997).

Product List Current prices may be obtained from our Web site or from our Customer Service Department.

Cat #	Product Name	Unit Size
C3018	Calcium Crimson™, AM *cell permeant* *special packaging*	
C3011MP	Calcium Green™-1, AM *cell permeant*	10
C3012	Calcium Green™-1, AM *cell permeant* *special packaging*	
C6765	Calcium Green™-1 dextran, potassium salt, 3000 MW, anionic	
C3713	Calcium Green™-1 dextran, potassium salt, 10,000 MW, anionic	
C3714	Calcium Green™-1 dextran, potassium salt, 70,000 MW, anionic	5 mg
C6766	Calcium Green™-1 dextran, potassium salt, 500,000 MW, anionic	5 mg
C3010MP	Calcium Green™-1, hexapotassium salt *cell impermeant*	500 µg
C3732	Calcium Green™-2, AM *cell permeant* *special packaging*	10 x 50 µg
C3730	Calcium Green™-2, octapotassium salt *cell impermeant*	500 µg
C36202	Calcium Yellow™, tetrapotassium salt *cell impermeant*	500 µg
C3739	Calcium Green™-5N, AM *cell permeant* *special packaging*	10 x 50 µg
C3737	Calcium Green™-5N, hexapotassium salt *cell impermeant*	500 µg
C3015	Calcium Orange™, AM *cell permeant* *special packaging*	10 x 50 µg
C3013	Calcium Orange™, tetrapotassium salt *cell impermeant*	500 µg
06807	Oregon Green [®] 488 BAPTA-1, AM *cell permeant* *special packaging*	
06798	Oregon Green® 488 BAPTA-1 dextran, potassium salt, 10,000 MW, anionic	
06806	Oregon Green® 488 BAPTA-1, hexapotassium salt *cell impermeant*	
06809	Oregon Green [®] 488 BAPTA-2, AM *cell permeant*	
06808	Oregon Green® 488 BAPTA-2, octapotassium salt *cell impermeant*	
06813	Oregon Green® 488 BAPTA-5N, AM *cell permeant* *special packaging*	10
06812	Oregon Green® 488 BAPTA-5N, hexapotassium salt *cell impermeant*	
023990	Oregon Green® 488 BAPTA-6F, hexapotassium salt *cell impermeant*	
023991	Oregon Green® 488 BAPTA-6F, AM *cell impermeant **special packaging *	
020001		10 X 00 µg

Contact Information

Further information on Molecular Probes products, including product bibliographies, is available from your local distributor or directly from Molecular Probes. Customers in Europe, Africa and the Middle East should contact our office in Paisley, United Kingdom. All others should contact our Technical Service Department in Eugene, Oregon.

Please visit our website-probes.invitrogen.com-for the most up-to-date information.

Molecular Probes, Inc. 29851 Willow Creek Road, Eugene, OR 97402 Phone: (541) 465-8300 • Fax: (541) 335-0504

Customer Service: 6:00 am to 4:30 pm (Pacific Time) Phone: (541) 335-0338 • Fax: (541) 335-0305 • probesorder@invitrogen.com

Toll-Free Ordering for USA:

Order Phone: (800) 438-2209 • Order Fax: (800) 438-0228

Technical Service: 8:00 am to 4:00 pm (Pacific Time) Phone: (541) 335-0353 • Toll-Free (800) 438-2209 Fax: (541) 335-0238 • probestech@invitrogen.com

Invitrogen European Headquarters

Invitrogen, Ltd. 3 Fountain Drive Inchinnan Business Park Paisley PA4 9RF, UK Phone: +44 (0) 141 814 6100 • Fax: +44 (0) 141 814 6260 Email: euroinfo@invitrogen.com Technical Services: eurotech@invitrogen.com

Molecular Probes products are high-quality reagents and materials intended for research purposes only. These products must be used by, or directly under the supervision of, a technically qualified individual experienced in handling potentially hazardous chemicals. Please read the Material Safety Data Sheet provided for each product; other regulatory considerations may apply.

Limited Use Label License

For research use only. Not intended for any animal or human therapeutic or diagnostic use. 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). 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 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. Invitrogen Corporation will not assert a claim against the buyer of infringement of the above patents 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) 335-0504.

Several Molecular Probes products and product applications are covered by U.S. and foreign patents and patents pending. All names containing the designation [®] are registered with the U.S. Patent and Trademark Office.

Copyright 2005, Molecular Probes, Inc. All rights reserved. This information is subject to change without notice.