

Fluorescent Calcium Indicators

I. Introduction

Calcium flux assays are preferred methods in drug discovery for screening G protein coupled receptors (GPCR). Our Fluo-8® and Rhod-4™ serial calcium detection reagents are the brightest green and red calcium indicators while our Cal-520® and Cal-590™ give the highest signal/background ratio for intracellular calcium detection due to their excellent retention in live cells. AAT Bioquest offers other calcium indicators such as Fluo-4, Fluo-3, Fura-2, Indo-1, Rhod-5N, and Rhod-2 AM in the highest possible quality.

Table 1. Spectral and Ca²⁺-Binding Properties of Calcium Detection Reagents

Ca ²⁺ Indicator	Catalog Numbers		Excitation	Emission	K _d of Ca ²⁺ -Binding
	Salt	AM Ester			
Calbryte™ 520	20656, 20658	20650, 20651, 20653	492 nm	514 nm	1.2 uM
Calbryte™ 520L	20642	20640	492 nm	524 nm	91 uM
Calbryte™ 520XL	20645	N/A	492 nm	524 nm	300 uM
Calbryte™ 590	20706	20700, 20701, 20702	573 nm	588 nm	1.4 uM
Calbryte™ 630	20727	20720, 20721, 20722	608 nm	626 nm	1.2 uM
Cal-520®	21135, 21136, 21140, 21141	21130, 21131	492 nm	514 nm	320 nM
Cal-520FF™	21144	21142, 21143	492 nm	514 nm	9.2 μM
Cal-520N™	21147	21146	492 nm	514 nm	90 μM
Cal-590™	20515, 20518	20510, 20511, 20512	573 nm	588 nm	561 nM
Cal-630™	20535, 20538	20530, 20531, 20532	608 nm	626 nm	792 nM
Cal-670™	20455	N/A	650 nm	675 nm	853 nM
Cal-770™	20460	N/A	750 nm	775 nm	850 nM
Cal 500™	20410	20412	390 nm	500 nm	303 nM
Cal Green™ 1*	20500	20501, 20502	506 nm	531 nm	190 nM
Cal Red™ R525/650	20588	20590, 20591	492 nm	525/650 nm	330 nM
Fluo-8®	21086, 21087, 21088, 21089	21080, 21081, 21082, 21083	490 nm	514 nm	389 nM
Fluo-8FF™	21102, 21103	21104, 21105	490 nm	514 nm	10 μM
Fluo-8H™	21095	21090, 21091	490 nm	514 nm	232 nM
Fluo-8L™	21098, 21099, 21100, 21101	21096, 21097	490 nm	514 nm	1.86 μM
Fluo-4	20555, 20556	20550, 20551, 20552	494 nm	516 nm	345 nM
Fluo-3	21016, 21017, 21018	21010, 21011, 21012, 21013	506 nm	526 nm	325 nM
Fluo-3FF	21019	21014	506 nm	526 nm	10 μM
Fluo-5F	20562	20560	494 nm	516 nm	2.3 μM
Fluo-5N	20567	20566	494 nm	516 nm	90 μM

Fura-2	21025, 21026	21020, 21021 21022, 21023	340/380 nm	510 nm	140 nM
Fura FF	21028	21027	340/380 nm	510 nm	5.5 μ M
Fura-8	21057, 21058	21055, 21056	354/415 nm	524 nm	260 nM
Fura-8 FF	20621	20620	354/415 nm	524 nm	6 μ M
Fura Red	21045, 21047	21046, 21048	436/471 nm	630/652	400 nM
Indo-1	21040, 21044	21030, 21032 21033, 21036	355 nm	400/475 nm	230 nM
Mag-Fluo-4	20400	20401	494 nm	516 nm	90 μ M
OG488 BAPTA-1**	20506	20507	494 nm	523 nm	170 nM
Rhod-4™	21128, 21119 21128, 21129	21120, 21121 21122, 21123	530 nm	555 nm	525 nM
Rhod-2	21067, 21068	21060, 21062 21063, 21064	549 nm	578 nm	570 nM
Rhod-FF	21075, 21076	21077, 21078	549 nm	549 nm	19 μ M
Rhod-5N	21072	21070	551 nm	577 nm	320 μ M

* Cal Green™ 1 is the same molecule to Calcium Green-1. ** OG488 BAPTA-1 is equivalent to Oregon Green 488 BAPTA-1

Table 2. Dextran, Biotin or Biocytin conjugated Fluorescent Calcium Indicators

Cat. #	Product Name	Unit	MW	Ex (nm) ²	Em (nm) ²
20605	Cal-520® -Biotin Conjugate	5x50 μ g	1112.40	492	514
20606	Cal-520® -Biocytin Conjugate	5x50 μ g	1341.55	492	514
20600	Cal-520®-Dextran Conjugate *MW 3,000*	1 mg	~4,000	492	514
20601	Cal-520®-Dextran Conjugate *MW 10,000*	5 mg	~11,000	492	514
20508	Cal-590™-Dextran Conjugate *MW 3,000*	1 mg	~4,000	573	588
20509	Cal-590™-Dextran Conjugate *MW 10,000*	1 mg	~11,000	573	588
20545	Cal-630™-Dextran Conjugate *MW 3,000*	1 mg	~4,000	608	626
20546	Cal-630™-Dextran Conjugate *MW 10,000*	1 mg	~11,000	608	626
20456	Cal-670™-Dextran Conjugate *MW 3,000*	1 mg	~4,000	650	675
20457	Cal-670™-Dextran Conjugate *MW 10,000*	1 mg	~11,000	650	675
20461	Cal-770™-Dextran Conjugate *MW 3,000*	1 mg	~4,000	750	775
20462	Cal-770™-Dextran Conjugate *MW 10,000*	1 mg	~11,000	750	775

II. Storage Conditions

Store at -20°C , protected from light. Expiration date is 12 months from the date of receipt.

III. Use of Calcium indicator AM Esters

1. Load Cells with Calcium Indicator AM Esters:

AM esters are the non-polar esters that readily cross live cell membranes, and rapidly hydrolyzed by cellular esterases inside live cells. AM esters are widely used for loading a variety of polar fluorescent probes into live cell non-invasively. However, cautions must be excised when AM esters are used since they are susceptible to hydrolysis, particularly in solution. They should be reconstituted in high-quality, anhydrous dimethylsulfoxide (DMSO). DMSO stock solutions should be stored desiccated at -20°C and protected from light. Under these conditions, AM esters should be stable for several months.

Following is our recommended protocol for loading AM esters into live cells. This protocol only provides a guideline, and should be modified according to your specific needs.

- a) Prepare a 2 to 5 mM AM esters stock solution in high-quality, anhydrous DMSO.
- b) On the day of the experiment, either dissolve calcium indicators solid in DMSO or thaw an aliquot of the indicator stock solutions to room temperature. Prepare a working solution of 2 to 20 μ M in the buffer of your choice (such as Hanks and Hepes buffer) with 0.04% *Pluronic*® F-127. For most cell lines we recommend the final concentration of calcium indicators be 4-5 μ M. The exact concentration of indicators required for cell loading must be determined empirically. To avoid any artifacts caused by overloading and potential dye toxicity, it is recommended to use the minimal probe concentration that can yield sufficient signal strength.
Note: The nonionic detergent Pluronic® F-127 is sometimes used to increase the aqueous solubility of calcium indicator AM esters. A variety of *Pluronic*® F-127 solutions can be purchased from AAT Bioquest.
- c) If your cells (such as CHO cells) containing the organic anion-transporters, probenecid (2–5 mM) or sulfinpyrazone (0.2–0.5 mM) may be added to the dye working solution (final in well concentration will be 1-2.5 mM for probenecid, or 0.1 -0.25 mM for sulfinpyrazone) to reduce the leakage of the de-esterified indicators.
Note: A variety of ReadiUse™ probenecid including water soluble sodium salt and stabilized solution can be purchased from AAT Bioquest
- d) Add equal volume of the dye working solution (from Step b or c) into your cell plate.
- e) Incubate the dye-loading plate room at temperature or 37 °C for 20 minutes (especially Fluo-8 AM) to 2 hours, and then incubate the plate at room temperature for another 30 minutes.
Note1: Decreasing the loading temperature might reduce the compartmentalization of the indicator.
Note2: Incubate the Cal-520 AM longer than 2 hours gives better signal intensity for some cell lines.
- f) Replace the dye working solution with HHBS or buffer of your choice (containing an anion transporter inhibitor, such as 1 mM probenecid, if applicable) to remove excess probes.
- g) Run the experiments at desired Ex/Em wavelengths (see Table 1).

2. Measure Intracellular Calcium Responses:

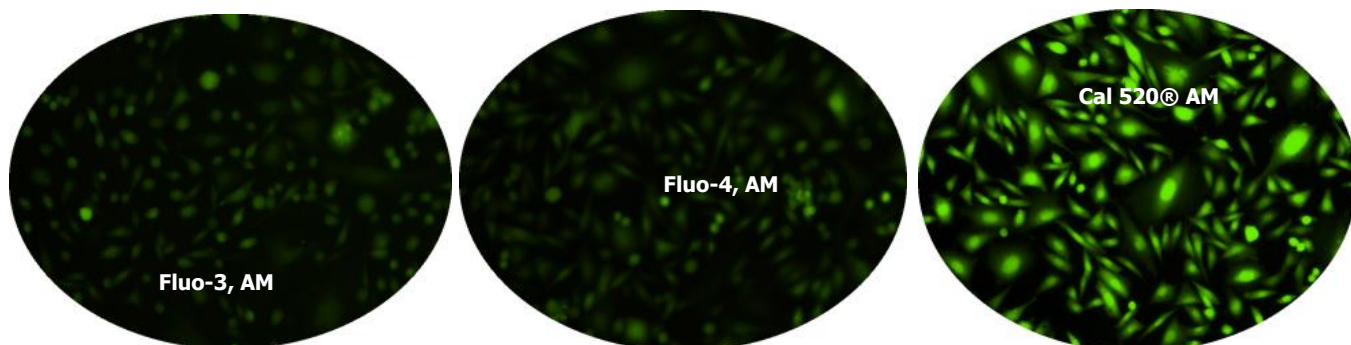


Figure 1. **Response of endogenous P2Y receptor to ATP in CHO-M1 cells without probenecid.** CHO-M1 cells were seeded overnight at 40,000 cells per 100 μ L per well in a 96-well black wall/clear bottom costar plate. 100 μ l of 4 μ M Fluo-3 AM, Fluo-4 AM or Cal 520® AM in HHBS were added into the wells, and the cells were incubated at 37 °C for 2 hour. The dye loading medium were replaced with 100 μ l HHBS, 50 μ l of 300 μ M ATP were added, and then imaged with a fluorescence microscope (Olympus IX71) using FITC channel.

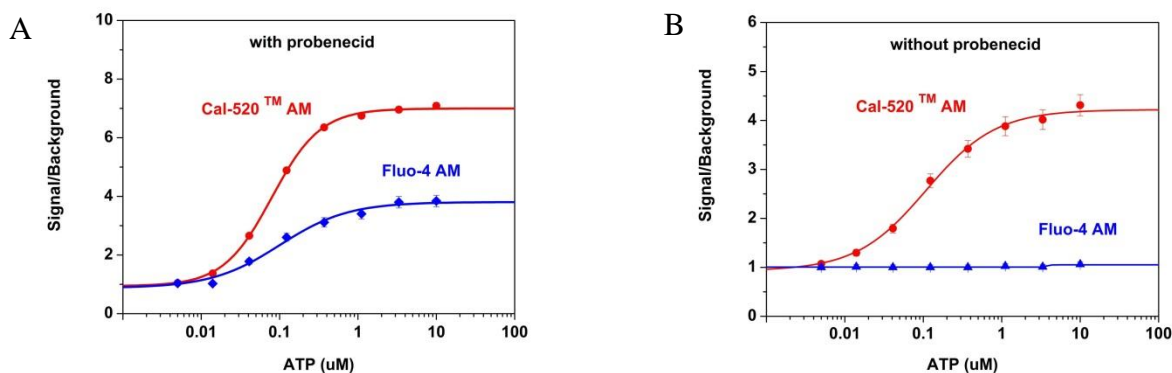


Figure 2. ATP-stimulated calcium response of endogenous P2Y receptor in CHO-K1 cells measured with Cal-520® or Fluo-4 AM. CHO-K1 cells were seeded overnight in 50,000 cells per 100 μ L per well in a 96-well black wall/clear bottom costar plate. 100 μ L of 5 μ M Fluo-4 AM or the Cal-520® AM with (A) or without (B) 2.5 mM probenecid was added into the cells, and the cells were incubated at 37°C for 2 hours. ATP (50 μ L/well) was added by FlexStation (Molecular Devices) to achieve the final indicated concentrations.

IV. Use of Calcium indicator Salts

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:

$$[\text{Ca}]_{\text{free}} = K_d[F - F_{\text{min}}]/F_{\text{max}} - F]$$

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 Ca^{2+} -binding and spectroscopic properties of fluorescent indicators vary quite significantly in cellular environments compared to calibration solutions. *In situ* calibrations of intracellular indicators typically yield K_d values significantly higher than in *in vitro* determinations. *In situ* calibrations are performed by exposing loaded cells to controlled Ca^{2+} buffers in the presence of ionophores such as A-23187, 4-bromo A-23187 and ionomycin. Alternatively, cell permeabilization agents such as digitonin or Triton® X-100 can be used to expose the indicator to the controlled Ca^{2+} levels of the extracellular medium. The K_d values of some calcium reagents are listed in Table 1 for your reference.

V. Use of Calcium indicator Conjugates

Compared to the free ion indicator, dextran conjugates of these same indicators exhibit both reduced compartmentalization and much lower rates of dye leakage. Since the molecular weight of the dextran, net charge, degree of labeling, and nature of the dye may affect the experiment, researchers are advised to consult the primary literature for information specific to the application of interest.

VII. References

1. J.T. Lock, I. Parker, I.F. Smith, A comparison of fluorescent Ca^{2+} indicators for imaging local Ca^{2+} signals in cultured cells, *Cell Calcium* (2015) October, <http://dx.doi.org/10.1016/j.ceca.2015.10.003>
2. Carsten Tischbirek, Antje Birkner, Hongbo Jia, Bert Sakmann, and Arthur Konnerth. Deep two-photon brain imaging with a red-shifted fluorometric Ca^{2+} indicator. *PNAS*. 2015; 112:11377-11382. doi: 10.1073/pnas.1514209112
3. Søren Grubb, Gary L. Aistrup, Jussi T. Koivumäki, Tobias Speerschneider, Lisa A. Gottlieb, Nancy A. M. Mutsaers, Søren-Peter Olesen, Kirstine Calloe, Morten B. Thomsen. Preservation of cardiac function by prolonged action potentials in mice deficient of KCHIP2 *American Journal of Physiology - Heart and Circulatory Physiology* Published 1 August 2015 Vol. 309 no. 3, H481-H489 DOI: 10.1152/ajpheart.00166.2015
4. Emery Smith, Peter Chase, Colleen M. Niswender, Thomas J. Utley, Douglas J. Sheffler, Meredith J. Noetzel, Atin Lamsal, Michael R. Wood, P. Jeffrey Conn, Craig W. Lindsley, Franck Madoux, Mary Acosta, Louis Scampavia, Timothy Spicer, and Peter Hodder. Application of Parallel Multiparametric Cell-Based FLIPR Detection Assays for the Identification of Modulators of the Muscarinic Acetylcholine Receptor 4 (M_4). *J Biomol Screen*. 2015; 20:858-868. doi:10.1177/1087057115581770.
5. Wenxiang Hu, Binlong Qiu, Wuqiang Guan, Qinying Wang, Min Wang, Wei Li, Longfei Gao, Lu Shen, Yin Huang, Gangcai Xie, Hanzhi Zhao, Ying Jin, Beisha Tang, Yongchun Yu, Jian Zhao, and Gang Pei Direct

- Conversion of Normal and Alzheimer's Disease Human Fibroblasts into Neuronal Cells by Small Molecules. *Cell Stem Cell* 17, 204–212, August 6, 2015. <http://dx.doi.org/10.1016/j.stem.2015.07.006>
6. Carsten Tischbirek, Antje Birkner, Hongbo Jia, Bert Sakmann, and Arthur Konnerth. Deep two-photon brain imaging with a red-shifted fluorometric Ca²⁺ indicator. *PNAS*. 2015; 112:11377-11382. doi: 10.1073/pnas.1514209112
 7. Songqing Tang, Taoyong Chen, Mingjin Yang, Lei Wang, Zhou Yu, Bin Xie, Cheng Qian, Sheng Xu, Nan Li, Xuetao Cao and Jianli Wang. Extracellular calcium elicits feedforward regulation of the Toll-like receptor-triggered innate immune response. *Cellular & Molecular Immunology*, (17 August 2015) | doi:10.1038/cmi.2015.59.
 8. Mayumi Tada, Atsuya Takeuchi, Miki Hashizume, Kazuo Kitamura, Masanobu Kano Article. A highly sensitive fluorescent indicator dye for calcium imaging of neural activity in vitro and in vivo. *European Journal of Neuroscience* 9 JAN 2014. DOI: 10.1111/ejn.12476.
 9. Daisuke Kodama, Akifumi Togari. Store-operated calcium entry induced by activation of Gq-coupled alpha1B adrenergic receptor in human osteoblast *Biochemical and Biophysical Research Communications* June (2013) doi: 10.1016/j.bbrc.2013.06.047.
 10. Rie Yamamoto, Shigeharu Ueki, Yuki Moritoki, Yoshiki Kobayashi, Hajime Oyamada, Yasunori Konno, Mami Tamaki, Masamichi Itoga, Masahide Takeda, Wataru Ito, and Junichi Chihara. Adiponectin attenuates human eosinophil adhesion and chemotaxis: implications in allergic inflammation. *Journal of Asthma* 2013. Posted online on July 17, 2013. (doi:10.3109/02770903.2013.816725).
 11. Alkhaldi, Jan Martinek, Brian Panicucci, Christophe Dardonville, Alena Zikova, Harry P. de Koning. Trypanocidal action of bisphosphonium salts through a mitochondrial target in bloodstream form *Trypanosoma brucei*. *International Journal for Parasitology: Drugs and Drug Resistance* 6 (2016) 23e34.
 12. Takahiro Shibata, Katsuhiko Takahashi, Yui Matsubara, Emi Inuzuka, Fumie Nakashima, Nobuaki Takahashi, Daisuke Kozai, Yasuo Mori & Koji Uchida. Identification of a prostaglandin D₂ metabolite as a neurogenesis enhancer targeting the TRPV1 ion channel. *Sci Rep*. 2016; 6: 21261. Published online 2016 Feb 16. doi: 10.1038/srep21261.
 13. Boris Gourévitchh, Jun Cai, Nicholas Mellen. Cellular and network-level adaptations to in utero methadone exposure along the ventral respiratory column in the neonate rat. *Experimental Neurology*. Available online 20 March 2016.
 14. Aditya J. Desai, Maoqing Dong, Laurence J. Miller. Beneficial effects of β-sitosterol on type 1 cholecystokinin receptor dysfunction induced by elevated membrane cholesterol. *Clinical Nutrition*. Available online 15 March 2016.
 15. Wiktor S. Phillips, Mikkel Herly, Christopher A. Del Negro, and Jens C. Rekling. **Organotypic slice cultures containing the preBötzinger complex generate respiratory-like rhythms**. *J Neurophysiol*. 2016; 115:1063-1070.
 16. Jin-Feng Zhao, Song-Kun Shyue, and Tzong-Shyuan Lee. Excess Nitric Oxide Activates TRPV1-Ca²⁺-Calpain Signaling and Promotes PEST-dependent Degradation of Liver X Receptor α. *Int J Biol Sci*. 2016; 12(1): 18–29. doi: 10.7150/ijbs.13549.

Warning: The products shall be only sold to our authorized distributors and end users. Calbryte™ series are covered by US 9,810,700 Cal-520® AM is covered by US 9,097,730, and Cal-590™ AM is covered by US 9,097,730. Fluo-8® AM is covered by US 8,779,165 and US 8,927,224. Neither resale nor transfer to a third party is allowed without written permission from AAT Bioquest. Chemical analysis of the products is strictly prohibited. Please call us at 408-733-1055 or e-mail us at info@aatbio.com if you have any questions.