

Premo™ FUCCI Cell Cycle Sensor *BacMam 2.0*

Catalog nos. P36237, P36238

Table 1 Contents and storage

Material	Amount	Concentration	Storage*	Stability
Premo™ geminin-GFP (G2/M reagent) (Component A)	1 mL (P36237) or 200 µL (P36238)	~1 × 10 ⁸ viral particles/mL	• 2–8°C • Desiccate • Protect from light • DO NOT FREEZE	When stored as directed, this kit is stable for at least 6 months.
Premo™ Cdt1-RFP (G1/S reagent) (Component B)	1 mL (P36237) or 200 µL (P36238)	~1 × 10 ⁸ viral particles/mL		
*These storage conditions are appropriate when storing the entire kit upon receipt. After preparing stock solutions, optimal storage conditions may change. For storing prepared stock solutions, follow recommendations included in this product information sheet.				
Number of assays: Sufficient material is supplied for 50–100 coverslips, based on the protocol below.				
Approximate fluorescence excitation/emission maxima: Premo™ geminin-GFP: 485/520 nm; Premo™ Cdt1-RFP: 555/584 nm.				

Introduction

In 2008, Miyawaki and colleagues developed the Fluorescence Ubiquitination Cell Cycle Indicator (FUCCI), a fluorescent protein (FP)-based sensor that employs a red (RFP) and a green (GFP) fluorescent protein fused to different regulators of the cell cycle: Cdt1 and geminin.¹ These two constructs, Cdt1 and geminin, are ubiquitinated by specific ubiquitin E3 ligases targeting them to the proteasome for degradation. The temporal regulation of the activity of these E3 ligases results in the biphasic cycling of geminin and Cdt1 through the cell cycle. In the G1 phase of the cell cycle, geminin is broken down and only Cdt1 tagged with RFP may be visualized, thus identifying cells in the G1 phase with red fluorescent nuclei. In the S, G2, and M phases, however, Cdt1 is degraded and only geminin tagged with GFP remains, thus identifying cells in these phases with green fluorescent nuclei. During the G1/S transition, as Cdt1 levels decrease and geminin levels increase, both proteins are present in the cells, allowing GFP and RFP fluorescence to be observed—when green and red images are overlaid, the cells appear with yellow fluorescent nuclei. This dynamic color change from red-to-yellow-to-green represents the progression through cell cycle and division (Figures 1 and 2, next page).

The Premo™ FUCCI Cell Cycle Sensor combines the Cdt1 and geminin FP constructs with the powerful BacMam gene delivery system. The genetically encoded and pre-packaged reagents enable immediate usage and eliminate the need to purify plasmid or to use lipids, dye-loading chemicals, or other potentially harmful treatments to transduce cells. Additionally, BacMam technology permits defined optimization as expression levels can easily be titrated by adding more or less virus to cells in culture. Cellular transduction is efficient and reproducible in most cell types, including primary and stem cells, without apparent cytotoxic effects. BacMam 2.0 greatly expands the efficiency and utility of this

popular gene delivery platform.²⁻⁴ Cell types previously not compatible with the technology (primary neurons), or cells that were poorly transduced with version 1.0 (some stem cells, CHO) can now be transduced quantitatively in a simple, one-step process. The improved performance is due to inclusion of elements that greatly enhance transduction efficiency and expression levels: a pseudotyped capsid protein for more efficient cell entry and genetic elements (enhanced CMV promoter and Woodchuck Post-Transcriptional Regulatory Element) that boost expression levels. To date, over 90 cell types have been shown to be effectively transduced using BacMam delivery technology. For the most up-to-date list of cells and transduction efficiencies, visit www.lifetechnologies.com/bacmamcompatible.

Premo™ FUCCI Cell Cycle Sensor is designed for live-cell imaging of cell cycle progression and can be used to assess the effect of drugs, siRNA, or other factors on the transition of cells through the cell cycle. The fluorescence from geminin-GFP and Cdt1-RFP have been demonstrated to be resistant to fixation with 4% formaldehyde and permeabilization with 0.1% Triton® X-100, thereby enabling processing of labeled cells with antibodies to other cellular targets. Each kit contains all of the components needed to label cells with the Premo™ FUCCI Cell Cycle Sensor using a transduction volume of 1 mL; however, the protocol can easily be adjusted for larger or smaller volumes. The workflow is straightforward: just add the reagent directly to your cells and incubate overnight to allow for the expression of fluorescent proteins. The next day, visualize cell cycle progression in populations of cells using traditional fluorescence microscopy (Figure 3, next page).

Figure 1 Dynamic color change of the Premo™ FUCCI Cell Cycle Sensor

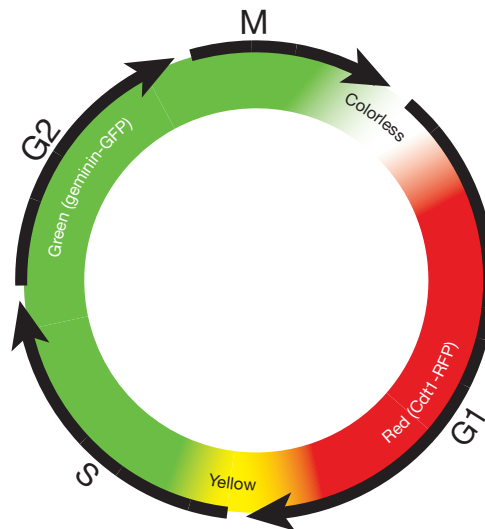


Figure 2 Fluorescence excitation and emission spectra for Premo™ geminin-GFP (panel A) and Premo™ Cdt1-RFP (panel B)

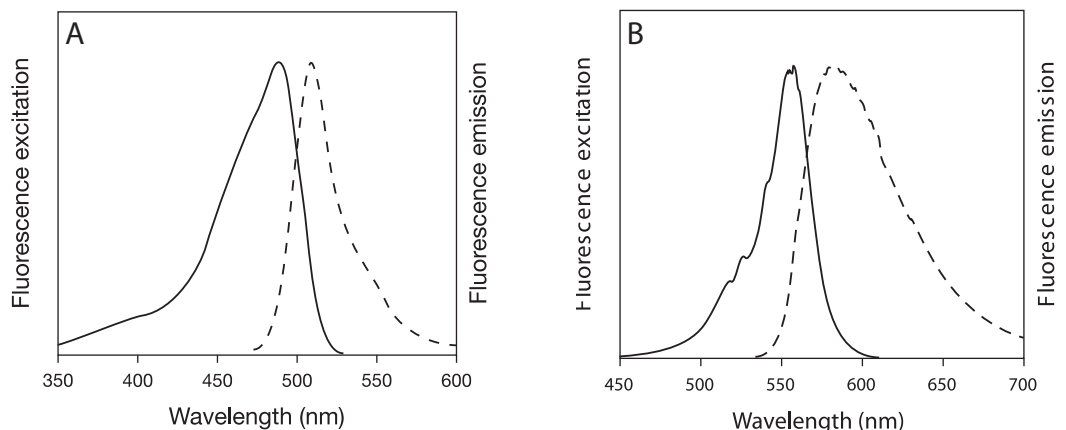
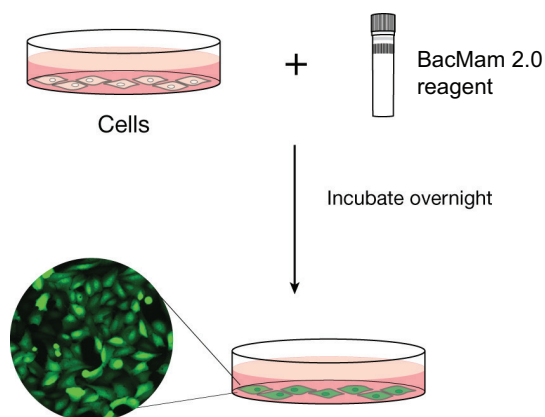


Figure 3 Workflow for Premo™ FUCCI Cell Cycle Sensor *BacMam 2.0*



Before Starting

Working with Premo™ FUCCI Cell Cycle Sensor Reagents

- Premo™ FUCCI Cell Cycle Sensor reagents work with most cell types between 20 and 100 particles per cell (PPC).
- For best results, transduce cells at a confluence of no more than 70%.
- The BacMam Enhancer (Cat. no. B10107) is generally not required for BacMam 2.0 reagents. However, its use has been shown to boost expression in some challenging cell types such as Jurkat.
- For optimal results you may need to alter the PPC, volume, cell density, temperature, or incubation time. Following the PPC, adjusting the volume is the next best parameter to change to optimize protein expression.

Experimental Protocols

Labeling Cells with Premo™ FUCCI Cell Cycle Sensor

The following protocol is based on a 1 mL labeling volume and ~50,000 cells plated in a 35-mm dish or for 1 well of a 6-well culture plate and a PPC (particles per cell) of 40. For applications that require a larger number of cells such as flow cytometry and high-content screening (HCS), we recommend plating cells in a 10-cm dish or T-75 flask and increasing the labeling volume to 5 mL with a proportionate increase in the volume of the virus.

- 1.1** Plate cells at a desired density and allow sufficient time for cells to adhere. BacMam reagents work best when used on cells of a low passage number plated at a low density. Plating density should take into account the desired confluence at the time of cell labeling and accommodate additional growth during the overnight incubation required for expression of geminin-GFP and Cdt1-RFP.

- 1.2 Calculate the volume of Premo™ geminin-GFP (Component A) and Premo™ Cdt1-RFP (Component B) using the equation below.

$$\text{Volume of Premo™ geminin-GFP or Premo™ Cdt1-RFP reagent (mL)} = \frac{(\text{number of cells})(\text{PPC})}{(1 \times 10^8)}$$

where the number of cells is the estimated total number of cells at the time of cell labeling, PPC (particles per cell) is the number of viral particles per cell, and 1×10^8 is the number of viral particles per mL of the reagent.

For example, to label 50,000 cells with a PPC of 40:

$$\text{mL of Premo™ geminin-GFP or Premo™ Cdt1-RFP reagent} = \frac{(50,000)(40)}{(1 \times 10^8)} = 0.02 \text{ mL (20 } \mu\text{L)}$$

Note: Sakaue-Sawano, *et. al.* report no deleterious effects of these modified proteins on the cell cycle;¹ however, we encourage optimizing the experimental conditions to avoid overexpressing geminin-GFP and Cdt1-RFP. Factors that affect labeling efficiency include PPC, labeling volume, and the incubation time. We have found an PPC of 40 to be a good starting point for expression studies. Table 2 provides initial guidelines for these factors; however, we recommend that you optimize for your specific cell type and application.

Table 2 Initial guidelines and factors for optimization of Premo™ FUCCI Cell Cycle Sensor labeling efficiency

Cell type	PPC	Labeling volume
Standard cell lines (HeLa, A549, CHO)	40	1 mL
Cell lines that are difficult to transfect	80	0.5 mL
	40	0.5 mL

- 1.3 Mix each Premo™ reagent (Components A and B) by inversion to ensure a homogenous solution.
- 1.4 Add the volume of Premo™ reagent calculated in step 1.2 directly to the cells in complete cell medium and mix gently.
- 1.5 Return the cells to the culture incubator overnight (≥ 16 hours).

Imaging and Analysis

Note: At this stage, you may trypsinize the cells and store frozen at -80°C or in liquid nitrogen for future use.

- 2.1 Image and analyze using appropriate instrument filter sets. Refer to Figure 2 (page 2) for spectral characteristics of geminin-GFP and Cdt1-RFP.

Premo™ FUCCI Cell Cycle Sensor is primarily designed for live-cell imaging of cell cycle progression, but the fluorescence from geminin-GFP and Cdt1-RFP have been demonstrated to be resistant to fixation with 4% formaldehyde and permeabilization with 0.1% Triton® X-100, enabling processing of labeled cells with antibodies to other cellular targets. You can accurately correlate Premo™ FUCCI Cell Cycle Sensor expression with cell cycle phase by measuring DNA content. For live cell imaging, the cell-permeant nucleic acid stains Hoechst 33342 and HCS NuclearMask™ Blue stains are spectrally compatible with the Premo™ FUCCI Cell Cycle Sensor fluorescence.

Frequently Asked Questions

Q: Will BacMam 2.0 transduce my cells?

A: The first generation BacMam reagents were shown to efficiently transduce over 90 cell types, including stable cell lines and primary cells. For the most up-to-date list of cells and transduction efficiencies, refer to www.lifetechnologies.com/bacmamcompatible. With BacMam 2.0, it is now possible to efficiently transduce primary neurons and stem cells.

Q: How long does expression last?

A: The duration of transgene expression depends on many factors, including transduction levels, cell division rates, mRNA, and protein stability. In most transformed cell lines, such as HeLa and CHO, expression lasts about 5 days. In cells that divide more slowly or show contact inhibition, such as some stem cells, primary cells, and neurons, we have observed bright staining and transgene expression for more than 2 weeks. For non-dividing, terminally differentiated cells we have observed expression for 2 to 4 weeks.

Q: Can I transduce with more than one BacMam reagent at a time?

A: Yes, this is one of the advantages of the system. For instance the Premo™ FUCCI Cell Cycle Sensor and the BacMam Kv7.2/7.3 Potassium Ion Channel reagent are based on optimized ratios of 2 BacMam constructs that give rise to a two-color cell cycle sensor and a functional heterotetrameric K channel, respectively.

Q: Will BacMam transduction hurt my cells?

A: BacMam transduction is well-tolerated, even at very high number of viral particles to cell ratios (>1,000). However, we have occasionally observed apparent cytotoxic effects by some BacMam reagents at very high virus levels; this may be due to the nature of the transgene. For this reason, we recommend using no more virus than is needed.

Q: If I freeze my cells after transduction, how long can I store them without reducing expression levels?

A: Our data show that transduced cells can be stored in liquid nitrogen for several months without reducing the level of transgene expression.

Q: Can transduction be optimized if my cells are difficult to transduce?

A: Yes. Try varying particle-to-cell ratio (PPC), incubation volume, temperature or duration, and cell density (if adherent cells are transduced). For adherent cells, we recommend a confluence of about 70%. Following the PPC, adjusting the volume is the next best parameter to change to optimize protein expression.

Q: Can a cell be transduced more than once?

A: Yes. Because transduction is well-tolerated, you can add more BacMam reagent after a few days if expression levels need to be boosted or if a different BacMam-based assay is needed.

Q: It's a virus—is it safe to use?

A: Yes. Baculoviruses are insect viruses that do not replicate in mammalian cells and are generally used under the safety precautions common for standard cell-based reagents.

References

1. Cell 132, 487 (2008); 2. Biochem Biophys Res Comm 349, 1220 (2006); 3. J Biotechnol 131, 1 (2007); 4. Mol Ther 17, 1585 (2009).

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

Cat. no.	Product Name	Unit Size
P36237	Premo™ FUCCI Cell Cycle Sensor *BacMam 2.0* *set of 2 vials*	1 mL set
P36238	Premo™ FUCCI Cell Cycle Sensor *BacMam 2.0* *set of 2 vials*	200 µL set
<i>Related Products</i>		
B10107	BacMam Enhancer Kit	1 kit
H3570	Hoechst 33342, trihydrochloride, trihydrate *10 mg/mL solution in water*	10 mL
H10325	HCS NuclearMask™ Blue stain *for 10 × 96-well plates* *2000X concentrate*	65 µL

Purchaser Notification

Corporate Headquarters

5791 Van Allen Way
Carlsbad, CA 92008
USA
Phone: +1 760 603 7200
Fax: +1 760 602 6500
Email: techsupport@lifetech.com

European Headquarters

Inchinnan Business Park
3 Fountain Drive
Paisley PA4 9RF
UK
Phone: +44 141 814 6100
Toll-Free Phone: 0800 269 210
Toll-Free Tech: 0800 838 380
Fax: +44 141 814 6260
Tech Fax: +44 141 814 6117
Email: euroinfo@invitrogen.com
Email Tech: eurotech@invitrogen.com

Japanese Headquarters

LOOP-X Bldg. 6F
3-9-15, Kaigan
Minato-ku, Tokyo 108-0022
Japan
Phone: +81 3 5730 6509
Fax: +81 3 5730 6519
Email: jpinfo@invitrogen.com

Additional international offices are listed at
www.lifetechnologies.com

These high-quality reagents and materials must be used by, or directly under the supervision of, a technically qualified individual experienced in handling potentially hazardous chemicals. Read the Safety Data Sheet provided for each product; other regulatory considerations may apply.

Obtaining Support

For the latest services and support information for all locations, go to www.lifetechnologies.com.

At the website, you can:

- Access worldwide telephone and fax numbers to contact Technical Support and Sales facilities
- Search through frequently asked questions (FAQs)
- Submit a question directly to Technical Support (techsupport@lifetech.com)
- Search for user documents, SDSs, vector maps and sequences, application notes, formulations, handbooks, certificates of analysis, citations, and other product support documents
- Obtain information about customer training
- Download software updates and patches

SDS

Safety Data Sheets (SDSs) are available at www.lifetechnologies.com/sds.

Certificate of Analysis

The Certificate of Analysis provides detailed quality control and product qualification information for each product. Certificates of Analysis are available on our website. Go to www.lifetechnologies.com/support and search for the Certificate of Analysis by product lot number, which is printed on the product packaging (tube, pouch, or box).

For Research Use Only. Not for human or animal therapeutic or diagnostic use.

LIFE TECHNOLOGIES CORPORATION AND/OR ITS AFFILIATE(S) DISCLAIM ALL WARRANTIES WITH RESPECT TO THIS DOCUMENT, EXPRESSED OR IMPLIED, INCLUDING BUT NOT LIMITED TO THOSE OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, OR NON-INFRINGEMENT. TO THE EXTENT ALLOWED BY LAW, IN NO EVENT SHALL LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) BE LIABLE, WHETHER IN CONTRACT, TORT, WARRANTY, OR UNDER ANY STATUTE OR ON ANY OTHER BASIS FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING BUT NOT LIMITED TO THE USE THEREOF.

Limited Warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at www.lifetechnologies.com/termsandconditions. If you have any questions, please contact Life Technologies at www.lifetechnologies.com/support.

Limited Use Label License: Research Use Only

The purchase of this product conveys to the purchaser the limited, non-transferable right to use the purchased amount of the product only to perform internal research for the sole benefit of the purchaser. No right to resell this product or any of its components is conveyed expressly, by implication, or by estoppel. This product is for internal research purposes only and is not for use in commercial services of any kind, including, without limitation, reporting the results of purchaser's activities for a fee or other form of consideration. For information on obtaining additional rights, please contact outlicensing@lifetech.com or Out Licensing, Life Technologies Corporation, 5791 Van Allen Way, Carlsbad, California 92008.

Limited Use Label License No. 21: Bac-to-Bac® and Bac-to-Bac® HT

This product is sold under patent license from Monsanto for research purposes only and no license for commercial use is included. Requests for licenses for commercial manufacture or use should be directed to Director, Monsanto Corporate Research, 800 N. Lindbergh, St. Louis, Missouri 63167.

Limited Use Label License No. 127: GFP with Heterologous Promoter

This product is sold under license from Columbia University. Rights to use this product are limited to research use only, and expressly exclude the right to manufacture, use, sell or lease this product for use for measuring the level of toxicity for chemical agents and environmental samples in cells and transgenic animals. No other rights are conveyed. Not for human use or use in diagnostic or therapeutic procedures. Inquiry into the availability of a license to broader rights or the use of this product for commercial purposes should be directed to Columbia Innovation Enterprise, Columbia University, Engineering Terrace-Suite 363, New York, New York 10027.

Limited Use Label License No. 198: Fluorescent Protein Products

Any use of this product by a commercial entity requires a separate license from either GE Healthcare or Invitrogen Corporation. For information on obtaining a commercial license to use this product, please refer to the contact information located at the bottom of this statement. No rights are conveyed to modify or clone the gene encoding GFP contained in this product. For information on licensing, contact Out Licensing, Life Technologies Corporation, 5791 Van Allen Way, Carlsbad, California 92008. Phone (760) 603-7200.

Limited Use Label License No. 306: Baculovirus Vectors

This product is for research use only by those researchers in laboratories of academic, government, industrial and/or clinical institutions engaged in the investigation of biological or biochemical processes, or research and development of biological products. This product is not to be used in the manufacture, use or sale of human or animal diagnostic, therapeutic or prophylactic products.

Limited Use Label License No. 332: BacMam Virus 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 solely in accordance with the accompanying product literature or manual. Purchase of this product does not convey a license to expand, amplify, or otherwise propagate the provided viral particles or to otherwise modify or alter the virus by any means.

The trademarks mentioned herein are the property of Life Technologies Corporation or their respective owners.

©2012 Life Technologies Corporation. All rights reserved.

