

BacMam Aequorin Kit

Catalog Numbers: A13740 and A13741

Literature Lot Number: V1

Literature Part Number: A13741.PIS

Revision date: 2 June 2011

FAST FACTS

BacMam Aequorin utilizes BacMam technology to transiently deliver aequorin, a luminescent calcium sensor, into mammalian cells.

Aequorin has been used extensively to monitor calcium changes in cultured, living mammalian cells and has become a standard method for interrogating calcium second messenger pathways in a high-throughput screening formats.

Upon binding of intracellular calcium ions, aequorin displays a "flash" style luminescence signal while consuming coelenterazine h as a substrate. The luminescence signal resolves any problems with autofluorescent compounds as well as providing an exceptional signal-to-noise ratio.

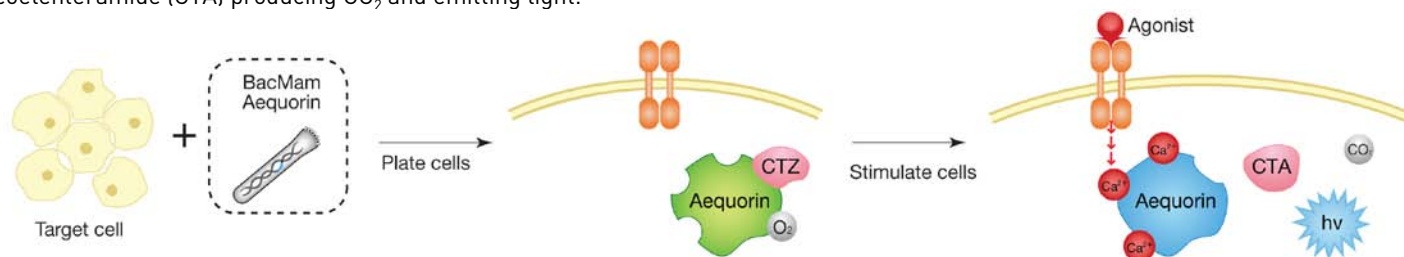
Due to the variation of cell types in which this product may be used, assay optimization may be required for best results (see BacMam Aequorin Transduction Guidelines, next page)

Refer to Kost, T.A., *et al. Drug Disc Today* 2007, 12, 396–403 for examples of BacMam gene expression in cells. For additional information on BacMam visit www.invitrogen.com/bacmam.

Kit components	Amount		Storage	Handling
	A13740	A13741		
BacMam Aequorin Reagent	10 mL	100 mL	4°C	<ul style="list-style-type: none"> Do not freeze Minimize exposure to ambient light Use aseptic technique Aliquot to minimize handling, if necessary
Enhancer Solution	400 µL	4 mL	–20°C	<ul style="list-style-type: none"> Protect from light

Additional materials required, but not provided	Source	Part number
Coelenterazine h	Invitrogen	C6780
Opti-MEM® I Reduced Serum Medium	Invitrogen	11058-021
Cell line of interest	Various	Various

Figure1 Aequorin Mechanism of Action. Aequorin-transduced cell lines are loaded with coelenterazine h (CTZ). Upon ligand-binding, intracellular calcium concentrations rise. The free calcium binds to the aequorin molecule leading to oxidation of CTZ to coelenteramide (CTA) producing CO₂ and emitting light.



BacMam Aequorin Transduction Guidelines

Due to the variation of cell types in which this product may be used, assay optimization may be required for best results.

- We recommend performing a virus titration to determine the optimal percentage of virus for transducing your cell line. Test a range of v/v dilutions of the BacMam reagent such as 50%, 25%, 12.5%, 6.25%, and 3.13% (v/v) for initial optimization.
- For difficult-to-transduce cells, such as primary and stem cell lines, we recommend following the transduction protocol described for BacMam GFP control (Catalog no. B10383)
- During the optimization process, we recommend incubating the cells after BacMam transduction in the presence and absence of the 1X BacMam Enhancer Solution. For cell lines such as CHO, HEK293T and HeLa, using the BacMam Enhancer Solution greatly increases the target expression. For certain cell types such as U2-OS, the Enhancer Solution may be toxic or non-beneficial. Add the enhancer solution to the cells during the 16–24 hours incubation period.
- For the best possible assay performance, we recommend optimizing the number of cells per well. For most cell lines, 2,000–10,000 cells per well in 25 μ L of medium in 384-well format is optimal.
- We recommend a clear-bottom plate when using instruments with bottom-read capabilities and a solid-bottom plate when using instruments with top-read capabilities.
- To optimize coelenterazine h loading time and temperature, we recommend testing 2, 4, 6, 8, and 24-hour loading times at room temperature and 37°C. Some cell lines can tolerate an overnight incubation at 37°C before significant signal intensity is lost.

Transduction Procedure

1. Add the following to a T-75 or 10-cm culture dish (both have approximately 75 cm² of surface area): 4 million cells, 9 mL of culture medium (with or without serum) and 1 mL (10%) BacMam Aequorin reagent.
2. Incubate the cells at 37°C and 5% CO₂ for 16–24 hours to allow for expression of aequorin protein.
3. Dissociate the cells from flask or dish by trypsinization or other method.
4. Count the cells using a hemacytometer or the Countess[®] Automated Cell Counter.
5. Centrifuge the cells and resuspend them at 1 million cells per 2.5 mL of Opti-MEM[®] I Reduced Serum Medium supplemented with 1% charcoal-dextran stripped FBS.
Note: (Optional) At this stage, cells can be cryopreserved and stored in liquid nitrogen for future use.
6. Add coelenterazine h to a final concentration of 5 μ M (1X). Mix well.
7. Plate the cells in the desired format (i.e., 96-, 384-, or 1,536-well plates) at one-half the working volume of the plate type (i.e., 25 μ L in a standard 384-well plate with a working volume of 50 μ L) using 10,000 cells/well at 25- μ L volume in a 384-well format.
8. Incubate the cells at 37°C or room temperature for 2–4 hours.

Calcium Mobilization Assay

Due to the “flash” style luminescence of aequorin, a luminometer with liquid handling injectors is needed. Examples include Hamamatsu FDSS6000 with luminescence option, Hamamatsu FDSS7000, MDS Analytical Technologies FLIPR Tetra[®] with luminescence option, Perkin Elmer’s LumiLux[®] and LumiLux[®] CS, CyBio CyBi[®]-Lumax or the BMG Labtech FLUOstar OPTIMA.

1. Make dilutions of test compounds in Opti-MEM[®] I Reduced Serum Medium at a 2X concentration.
Note: A calcium mobilization experiment is typically performed by first collecting a baseline reading for approximately 10 seconds before initiating calcium mobilization by compound addition.
2. Add an equal volume of compound to the cells in the plate (i.e., 25 μ L of compound to 25 μ L of cells in a 384-well plate).
Note: The aequorin signal should appear immediately after compound addition and only last for a few seconds
3. Monitor the luminescent signal for 45 seconds after adding the compound.
4. Process the data by computing the area under the curve and plotting that value versus compound concentration.

Technical Support

For additional assistance in using BacMam Aequorin Reagent, contact Technical Support. Email: probestech@invitrogen.com or phone (800) 438-2209 or (541) 335-0353.

Frequently Asked Questions

Visit www.invitrogen.com and search for A13740 or A13741 to download answers to Frequently Asked Questions (FAQs) for this product. The FAQ sheet is located under the “How to Use” tab on the product page.

Purchaser Notification

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 applications of any kind, including, without limitation, quality control and commercial services such as 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, 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. 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. 308: WPRE Element

This product contains the Woodchuck Post-transcriptional Regulatory Element ("WPRE") which is the subject of intellectual property owned by The Salk Institute for Biological Studies, and licensed to Life Technologies 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). 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) not to 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; and/or (4) resale of the product or its components, whether or not such product or its components are resold for use in research. In addition, any use of WPRE outside of this product or the product's authorized use requires a separate license from the Salk Institute. Life Technologies will not assert a claim against the buyer of infringement of patents owned by Life Technologies and claiming 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 or for a Commercial Purpose. If the purchaser is not willing to accept the limitations of this limited use statement, Life Technologies 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 Licensing Department, Life Technologies Corporation, 5791 Van Allen Way, Carlsbad, California 92008, Phone (760) 603-7200. Fax (760) 602-6500, or The Salk Institute for Biological Studies, 10010 North Torrey Pines Road, La Jolla, CA 92037, Attn.: Office of Technology Management, Phone: (858) 453-4100 extension 1275, Fax: (858) 546-8093.

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.

FLIPR Tetra is a trademark of Molecular Devices, Inc.

LumiLux is a trademark of PerkinElmer LAS, Inc.

CyBi is a trademark of CyBio AG.

Lumax is a trademark of Monobind, Inc.

©2011 Life Technologies Corporation. All rights reserved.