

NovaBright™ β-galactosidase Enzyme Reporter Gene Chemiluminescent Detection System for Yeast or Mammalian Cells

Catalog no. N10563, N10564, N10565, N10566

Table 1. Contents and storage information.

Material	N10563	N10564	N10565	N10566	Storage*	Stability
Galacton-Star® substrate concentrate (Component A)	1 mL	4.2 mL	1 mL	4.2 mL	<ul style="list-style-type: none">• 2–6°C• Protect from light• DO NOT FREEZE	When stored as directed this kit is stable for 1 year.
Assay buffer *for mammalian cells* (Component B)	20 mL	100 mL	--	--		
Assay buffer *for yeast cells* (Component B)	--	--	20 mL	100 mL		
Number of assays: Sufficient material is supplied for 200 (Cat. no. N10563, N10565) or 1,000 (Cat. no. N10564, N10566) microplate assays based on the protocol below.						

Introduction

The NovaBright™ β-galactosidase Enzyme Reporter Gene Chemiluminescent Detection System is designed for the rapid, simple, and sensitive detection of β-galactosidase reporter enzyme directly in microwell cultures of mammalian or yeast cells. The assay system incorporates the Galacton-Star® substrate with a luminescence enhancer to generate glow light emission kinetics. A single reagent, which provides cell lysis components and luminescent substrate and enhancer, is added directly to cells in culture medium. Light emission typically reaches maximum intensity within 60–90 minutes and remains constant for 45–90 minutes or longer (depending upon assay temperature, see Figure 1). Light signal is measured in a luminometer, without the need for automatic injection. The simple assay format and glow light emission achieved with the NovaBright™ β-galactosidase Enzyme Reporter Gene Chemiluminescent Detection System reagents provide an ideal assay system for automated high-throughput screening applications, enabling simple processing and measurement of multiple microplates. The dynamic range of the assay spans five orders of magnitude, from picogram to nanogram levels, enabling detection of a wide range of reporter enzyme concentration in cells.

The NovaBright™ β-galactosidase Enzyme Reporter Gene Chemiluminescent Detection System provides a chemiluminescent assay format that is sensitive, simple to use, and designed specifically for optimal performance in automated, high-throughput screening applications. The protocol is adaptable for use in multiple microplate formats with mammalian or yeast cells.

Applications

The bacterial β-galactosidase gene is widely used as a reporter enzyme for the study of gene regulation, for identification of protein-protein interactions and in assays for cell fusion.

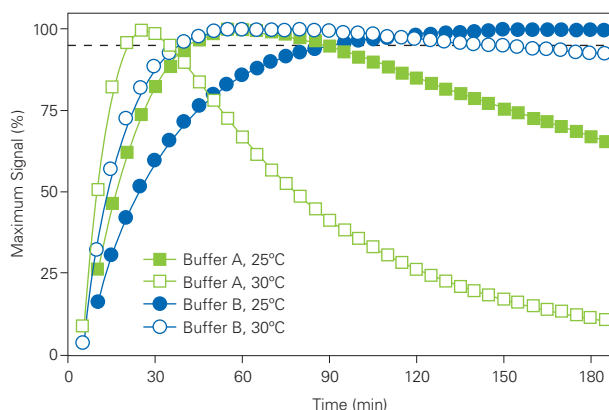


Figure 1. NovaBright™ β -galactosidase Enzyme Reporter Gene Chemiluminescent Detection assays were performed with purified β -galactosidase (500 pg) with assay buffer for mammalian cells (Buffer A) or assay buffer for yeast cells (Buffer B). Repeated measurements were made in the Applied Biosystems TR717™ microplate luminometer at room temperature (25°C) or 30°C. Data is presented as percentage of maximum signal, and dotted line indicates 95% of maximum signal.

Chemiluminescent 1,2-dioxetane substrates for β -galactosidase, including Galacton-Plus® and Galacton-Star® substrates, provide highly sensitive enzyme detection¹⁻⁴ and have been utilized extensively in reporter assays in mammalian cell and tissue extracts, and in a combined assay for luciferase and β -galactosidase activities^{5,6} (NovaBright™ β -galactosidase and Firefly Luciferase Dual Enzyme Reporter Gene Chemiluminescent Detection System, Invitrogen Cat. no. N10561 and N10562).

The NovaBright™ β -galactosidase Enzyme Reporter Gene Chemiluminescent Detection System is used widely for traditional reporter gene assays in:

- Transiently and stably-transfected mammalian cells^{7,8} including assays for studying viral infectivity and function⁹⁻¹²
- Yeast cells,¹³⁻¹⁵ including quantitative yeast two hybrid¹⁷⁻¹⁹ and one hybrid analysis²⁰
- Fish cells²¹
- Bacterial cells²²
- β -galactosidase complementation assays used for intracellular monitoring of protein-protein interactions,²³ protein translocation,²⁴ cell fusion,²⁵ and receptor dimerization/activation,^{26,27} including for high-throughput compound screening for receptor activation²⁸

Before Starting

Materials Required but Not Provided

- β -galactosidase reporter expressing mammalian or yeast cells
- 96- or 384-well tissue culture-treated microplates (solid white or clear-bottom white)
- Microplate luminometer
- *Optional:* β -galactosidase (Sigma Cat. no. G-5635)

Caution

Component A is highly flammable and irritating to eyes. Vapors may cause drowsiness and dizziness. Keep the container tightly closed and away from sources of ignition. No smoking. Avoid contact with eyes and skin. In case of contact with eyes, rinse immediately with plenty of water and seek medical advice.

General Guidelines

Assay Plates

The use of white, tissue culture-treated microplates is recommended for optimal assay sensitivity. Clear-bottom white plates can be used to allow microscopic examination of cultures. White backing sheets may be applied to the plate bottom prior to signal measurement, if desired. The absolute signal will be higher (approximately 2-fold), since the white backing reflects light toward the photomultiplier tube detector and eliminates light absorption by the black plate platform, but relative signal levels are unaffected. Black plates can be used but are not recommended, since they cause reduced light yield due to absorbance of the luminescent signal.

When using 384-well plates, reduce the culture and assay reagent volume to 25 μ L (or as desired, maintaining a 1:1 ratio of culture:reagent volume without overfilling the wells).

Culture Medium

Mammalian cell culture medium containing phenol red indicator can be used without affecting assay sensitivity. The inclusion of phenol red can result in a reduction of signal due to absorbance of some of the emitted light, but the signal/noise is not affected. The type of culture medium and presence/absence of serum may contribute to some slight variability in light signal and assay kinetics. Initially perform a kinetic analysis of the particular cell line/culture medium system to determine optimum signal measurement time. It is not necessary to measure light signal at the time of peak light emission, as long as all samples are measured in a consistent time frame.

Endogenous β -galactosidase Activity

Chemiluminescent β -galactosidase reporter assays may be conducted in mammalian cells that have endogenous β -galactosidase activity. The reaction is performed at a pH which is more favorable for bacterial enzyme activity, which has a higher pH optimum than the endogenous mammalian enzyme. Assay the level of endogenous enzyme in non-transfected cells to determine the assay background signal.

Assay Temperature

Assays are ideally performed at 26–28°C. If the incubation temperature is significantly different, the kinetics of light emission vary, since the rate of an enzyme reaction is dependent on temperature (see Figure 1).

Luminometers

We recommend using a single-mode luminometer or a multi-mode detection instrument set for luminescence measurement to measure light emission from 96- or 384-well microplates. The linear range of detection varies according to cell type and on the reporter enzyme expression level. Optimize the number of cells used per well to prevent a measurement signal that is outside the linear range of the luminometer. Extremely high light signals can saturate the detector (very unlikely for experimental samples), resulting in erroneous measurements. Refer to your luminometer user's manual to determine the upper limit for your specific luminometer.

Controls

The recommended positive and negative controls are described below:

β -Galactosidase Positive Control

Prepare stock enzyme by reconstituting lyophilized β -galactosidase (Sigma Cat. no. G-5635) to 1 mg/mL in 0.1 M sodium phosphate pH 7.0, 0.1% BSA. Store at 4°C. Generate a standard curve by serially diluting the stock enzyme in cell culture medium. For the high end of the dilution curve, use 2–20 ng of enzyme. Purified enzyme provides a positive control for the assay reagents, as well as a means to determine the range of detection of the luminometer

instrumentation, if desired. The purified enzyme standard curve is not intended (or accurate) for absolute quantitation of reporter enzyme concentrations, as the specific activity of the purified enzyme preparation and the reporter enzyme may differ significantly. Additional positive controls can include use of control β -galactosidase constructs that provide constitutive expression of reporter enzyme as a positive control for cell transfection.

Negative Control

Assay a volume of mock-transfected extract equivalent to that of experimental extract to determine endogenous cellular background. In experiments involving induction of reporter expression, assay uninduced cells as a negative control for total assay background.

Amount of Cells

For mammalian cell cultures, a typical cell density is 10,000–50,000 cells/well in 100 μ L for 96-well plates or 1,000–10,000 cells/well in 25 μ L for 384-well plates.

For yeast cells, a typical cell density is 10,000–75,000 cells/well in 100 μ L for 96-well plates or 5,000–40,000 cells/well in 25 μ L for 384-well plates.

Experimental Protocol

β -Galactosidase Detection Protocol

Perform all assays in duplicate or triplicate. See guidelines above for preparing positive and negative controls.

Assay buffer for mammalian cells is optimized for mammalian cell cultures, and has been used with transfected adherent and non-adherent cell lines, including NIH/3T3, CHO-K1, and K562. Assay buffer for yeast cells is optimized for yeast cells, and is also been shown to work with mammalian cells. The alternative buffer formulations result in differing light emission kinetics. For mammalian cells, the assay buffer for mammalian cells provides faster kinetics, which may be advantageous at lower incubation temperatures. The choice of reagents depends on the particular assay requirements and desired kinetic performance.

- 1.1 Prepare diluted substrate fresh each time. Dilute the substrate 1:25 with assay buffer (i.e., mix 40 μ L substrate with 960 μ L assay buffer, 100 μ L/well required) to make reaction assay buffer.
- 1.2 Add 100 μ L/well of reaction assay buffer to a 96-well plate containing 100 μ L/well of cells in culture medium.
- 1.3 Incubate the samples at 26–28°C for approximately 60–90 minutes or until constant light emission is reached.
- 1.4 Place the plate in a luminometer and measure for 0.1–1 second/well.

References

1. Bioluminescence and Chemiluminescence: Molecular Reporting with Photons (Hastings, JW, Kricka, LJ, and Stanley, PE, eds.), John Wiley, Chichester, England (1997) p. 525-528; 2. Bioluminescence and Chemiluminescence: Fundamentals and Applied Aspects (Campbell, AK, Kricka, LJ, and Stanley, PE, eds), John Wiley, Chichester, England (1994) p. 20-23; 3. Clin Chem 42, 1542 (1996); 4. Anal Biochem 199, 119 (1991); 5. Bioluminescence and Chemiluminescence: Molecular Reporting with Photons (Hastings, JW, Kricka, LJ, and Stanley, PE, eds), John Wiley, Chichester, England (1997), p. 451-457; 6. BioTechniques 21, 520 (1996); 7. Genes & Devel 18, 2120 (2004); 8. BMC Genomics 8,390 (2007); 9. J Biol Chem 280, 9345

(2005); **10.** Antimicrob Agents Chemother 49, 1139 (2005); **11.** J Virology 79, 1713 (2005); **12.** J Virol 80, 130 (2006). 137; **13.** J Biol Chem 280, 8906 (2005); **14.** Genetics 159, 159 (2001); **15.** Proc Natl Acad Sci USA 97, 12985 (2000); **16.** Nature Cell Biol 7, 1179 (2005); **17.** Toxicological Sciences 57, 54 (2000); **18.** Curr Genet 44, 184 (2003); **19.** J Virol 77, 4113 (2003); **20.** Blood 110, 3039 (2007); **21.** Biotechnol Letters 23, 481 (2001); **22.** J Bacteriol 186, 4694 (2004); **23.** Mol Cell Biol 24, 7456 (2004); **24.** Nature Methods 2, 521 (2005); **25.** J Cell Biol 180, 1005 (2008); **26.** Nature Biotech 18, 218 (2000); **27.** J Biol Chem 278, 15217 (2003); **28.** J Biomol Screening 7, 451 (2002).

Product List

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

Cat. no.	Product Name	Unit Size
N10563	NovaBright™ β -galactosidase Enzyme Reporter Gene Chemiluminescent Detection Kit *for mammalian cells* *200 assays*	1 kit
N10564	NovaBright™ β -galactosidase Enzyme Reporter Gene Chemiluminescent Detection Kit *for mammalian cells* *1000 assays*	1 kit
N10565	NovaBright™ β -galactosidase Enzyme Reporter Gene Chemiluminescent Detection Kit *for yeast cells* *200 assays*	1 kit
N10566	NovaBright™ β -galactosidase Enzyme Reporter Gene Chemiluminescent Detection Kit *for yeast cells* *1000 assays*	1 kit
Related Products		
N10561	NovaBright™ β -galactosidase and Firefly Luciferase Dual Enzyme Reporter Gene Chemiluminescent Detection Kit *200 assays*	1 kit
N10562	NovaBright™ β -galactosidase and Firefly Luciferase Dual Enzyme Reporter Gene Chemiluminescent Detection Kit *600 assays*	1 kit

Contact Information

Corporate Headquarters

5791 Van Allen Way
Carlsbad, CA 92008
Phone: (760) 603-7200
Fax: (760) 602-6500

European Headquarters

Inchinnan Business Park.
3 Fountain Drive
Paisley PA4 9RF, UK
Phone: +44 (0) 141 814 6100
Fax: +44 (0) 141 814 6117
Email: euinfo@invitrogen.com
Technical Services: euotech@invitrogen.com

Japanese Headquarters

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

For country-specific contact information, visit www.invitrogen.com.

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

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 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. 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) 335-0354.

Limited Warranty

Invitrogen (a part of Life Technologies Corporation) is committed to providing our customers with high-quality goods and services. Our goal is to ensure that every customer is 100% satisfied with our products and our service. If you should have any questions or concerns about an Invitrogen product or service, contact our Technical Support Representatives. All Invitrogen products are warranted to perform according to specifications stated on the certificate of analysis. The Company will replace, free of charge, any product that does not meet those specifications. This warranty limits the Company's liability to only the price of the product. No warranty is granted for products beyond their listed expiration date. No warranty is applicable unless all product components are stored in accordance with instructions. The Company reserves the right to select the method(s) used to analyze a product unless the Company agrees to a specified method in writing prior to acceptance of the order. Invitrogen makes every effort to ensure the accuracy of its publications, but realizes that the occasional typographical or other error is inevitable. Therefore the Company makes no warranty of any kind regarding the contents of any publications or documentation. If you discover an error in any of our publications, please report it to our Technical Support Representatives. Life Technologies Corporation shall have no responsibility or liability for any special, incidental, indirect or consequential loss or damage whatsoever. The above limited warranty is sole and exclusive. No other warranty is made, whether expressed or implied, including any warranty of merchantability or fitness for a particular purpose.

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

Copyright 2009, Life Technologies Corporation. All rights reserved. For research use only. Not intended for any animal or human therapeutic or diagnostic use.