

Certificate of Analysis

FastBreak™ Cell Lysis Reagent, 10X:

Part No.	Size (ml)
V857A	15
V857B	60
V857C	100

Description: FastBreak™ Cell Lysis Reagent^(a,b,c) is a proprietary formulation of buffer and detergents designed to lyse *E. coli* cells in culture.

Storage Conditions: Store at 4–25°C. A precipitate may form at low temperature. If this occurs, warm the reagent to room temperature before use.

Part# 9PIV857

Revised 8/16



AF9PIV857 0816V857

Quality Control Assays

Lysis Efficiency

Cells are induced to express luciferase and then grown in LB overnight to an $OD_{600} > 3$. One hundred microliters of FastBreak™ Reagent is added to 900µl of cell culture and mixed for 10 minutes at room temperature. A sample of the crude lysate is removed and centrifuged to remove cell debris. Luciferase activity in the supernatant of the cleared lysate must be at least 50% of the luciferase activity measured in the the same volume of crude lysate.



Promega

Promega Corporation

2800 Woods Hollow Road	
Madison, WI 53711-5399	USA
Telephone	608-274-4330
Toll Free	800-356-9526
Fax	608-277-2516
Internet	www.promega.com

PRODUCT USE LIMITATIONS, WARRANTY, DISCLAIMER

Promega manufactures products for a number of intended uses. Please refer to the product label for the intended use statements for specific products. Promega products contain chemicals which may be harmful if misused. Due care should be exercised with all Promega products to prevent direct human contact.

Each Promega product is shipped with documentation stating specifications and other technical information. Promega products are warranted to meet or exceed the stated specifications. Promega's sole obligation and the customer's sole remedy is limited to replacement of products free of charge in the event products fail to perform as warranted. Promega makes no other warranty of any kind whatsoever, and SPECIFICALLY DISCLAIMS AND EXCLUDES ALL OTHER WARRANTIES OF ANY KIND OR NATURE WHATSOEVER, DIRECTLY OR INDIRECTLY, EXPRESS OR IMPLIED, INCLUDING, WITHOUT LIMITATION, AS TO THE SUITABILITY, PRODUCTIVITY, DURABILITY, FITNESS FOR A PARTICULAR PURPOSE OR USE, MERCHANTABILITY, CONDITION, OR ANY OTHER MATTER WITH RESPECT TO PROMEGA PRODUCTS. In no event shall Promega be liable for claims for any other damages, whether direct, incidental, foreseeable, consequential, or special (including but not limited to loss of use, revenue or profit), whether based upon warranty, contract, tort (including negligence) or strict liability arising in connection with the sale or the failure of Promega products to perform in accordance with the stated specifications.

Signed by:

R. Wheeler, Quality Assurance

© 2003, 2005, 2013, 2016 Promega Corporation. All Rights Reserved.

FastBreak, MagneGST and MagneHis are trademarks of Promega Corporation.

Biomek is a registered trademark of Beckman Coulter, Inc.

Products may be covered by pending or issued patents or may have certain limitations. Please visit our Web site for more information.

All specifications are subject to change without prior notice.

Product claims are subject to change. Please contact Promega Technical Services or access the Promega online catalog for the most up-to-date information on Promega products.

Part# 9PIV857
Printed in USA, Revised 8/16.

^(a)This product is licensed for use under U.S. Pat. No. 6,174,704.

^(b)Patent Pending.

^(c)Certain applications of this product are covered by patents issued and applicable in certain countries. Because purchase of this product does not include a license to perform any patented application, users of this product may be required to obtain a patent license depending upon the particular application and country in which the product is used.

1. Description

FastBreak™ Cell Lysis Reagent is designed for the efficient, gentle lysis of *E. coli* cultures without the need for centrifugation or mechanical cell disruption. This reagent is provided as a 10X concentrate and contains a proprietary nonionic detergent to facilitate lysis. The reagent is added directly to *E. coli* cultures. FastBreak™ Reagent has a different formulation than the other cell lysis buffers used in Promega protein purification systems, which are designed for use with cell pellets. Following a brief incubation, the cells are disrupted, and the protein of interest is released. Recombinant proteins can be directly screened in the resulting extract or purified by the addition of an appropriate affinity matrix such as the MagneHis™ Protein Purification System (Cat.# V8500). This format allows the procedure to be performed manually or on a robotic platform, such as the Beckman Coulter Biomek® 2000 or FX workstations, for high-throughput applications.

2. Manual Protocol

(For robotic procedures refer to: www.promega.com/automethods/)

A. Cell Lysis

Reagents to Be Supplied by the User

- *E. coli* culture
- shaker or rotary platform

1. Add 100µl of FastBreak™ Reagent to 900µl of bacterial culture (see Note 1).
2. Mix for 10–15 minutes at room temperature on a shaker or rotary platform. Adequate mixing is necessary to ensure complete lysis.
3. Proceed with screening or purification. We have observed reduced yields when using the FastBreak™ Buffer with the MagneGST™ Protein Purification System (Cat.# V8600).

Notes:

1. When lysing cultures that have reached a high density (e.g., $OD_{600} > 5$) or when using the *E. coli* strain, BL21(DE3)pLysS, the lysates produced are often viscous due to the release of genomic DNA. This can be resolved by adding DNase (e.g., 20–40 units of RQ1 DNase [Cat.# M6101] or 4 Kunitz units of Sigma DNase [Cat.# D5025] freshly resuspended in 10X FastBreak™ Reagent) and incubating for 10–20 minutes at room temperature on a shaker or rotary platform. The lysate will become less viscous after DNase treatment. If the lysate is still too viscous for easy handling, add more DNase or increase the incubation time.
2. When using MagneHis™ Ni-Particles, it may not be apparent that DNase is needed until after the particles have been added to the cell lysate. In very viscous lysates, the particles will not mix uniformly in the lysate and may not be captured by the magnet. The samples may be treated with DNase after addition of the particles.
3. Lysozyme can be used to enhance cell lysis. Lysozyme (from egg white) is active in the presence of FastBreak™ Reagent. Freshly prepare a 20mg/ml solution of lysozyme in TE buffer (pH 7.4). Add 10µl of lysozyme solution per milliliter of lysate and incubate for 10 minutes. Due to the increased cell lysis when using lysozyme, it is also necessary to add DNase to reduce the viscosity of the lysate. Since lysozyme binds to MagneHis™ Ni-Particles, NaCl (0.5M final concentration) should be added to the lysate following the DNase treatment when using the MagneHis™ Protein Purification System.
4. FastBreak™ Reagent can be used to lyse cells that have been stored at 4°C overnight. However, storage of cells may not be optimal for all expressed proteins.
5. In some cases a precipitate may be seen. This occurs most often with Terrific Broth purification steps.

FastBreak™ Cell Lysis Reagent is a component of the MagneHis™ Protein Purification System. For a detailed protocol using the FastBreak™ Cell Lysis Reagent with the MagneHis™ Protein Purification System, please see the *MagneHis™ Protein Purification System Technical Manual #TM060* (www.promega.com/tbs/tm060/tm060.html).