

BL21(DE3) Chemi-Competent Expression Cells 156-3003

10 x 0.05 ml

For research use only

Store at -70°C

Storage and Stability

Store the BL21 (DE3) competent cells at -70°C in a constant temperature freezer. When stored under these conditions, the cells are stable for 6 months after date of receipt.

Kit Contents

10 x 50 μl BL21 (DE3) chemically competent cells

1 x 10 μl pUC19 control plasmid DNA (10 ng/ μl) in 10 mM Tris-HCl pH 8.0, 1 mM EDTA

1 x 10 ml SOC medium*

* (2% Tryptone, 0.5% yeast extract, 0.4% glucose, 10 mM NaCl, 2.5 mM KCl, 5 mM MgCl_2 , 5 mM MgSO_4)

Description

BL21 (DE3) competent cells are chemically competent *E. coli* cells used for high-level protein expression with T7 RNA polymerase-based expression systems. The strain is a derivative of *E. coli* B, which is deficient in the lon protease as well as the ompT outer membrane protease, facilitating the isolation of intact recombinant proteins. The host is a lysogen of λDE3 and contains the T7 RNA polymerase gene, under the control of the *lacUV5* promoter, integrated into the chromosome. IPTG is used to induce the expression of recombinant proteins cloned into vectors downstream of a T7 RNA promoter and transformed into the BL21 (DE3) cells.

BL21 (DE3) competent cells are packaged in single-use volumes for convenience and maintenance of efficiency. The recommended transformation protocol will yield 10 x 50 μl transformations.

Genotype

E. coli B F⁻ *dcm ompT hsdS*(r_B⁻, m_B⁻) *gal* λ (DE3)

Quality Control

Each batch of BL21 (DE3) competent cells is performance tested to ensure yields of $\geq 1 \times 10^7$ CFU/ μg DNA using nonsaturating amounts of pUC19 control plasmid.

Protocol Notes

1. Thaw cells on wet ice. For best results, use within 15 min of thawing. Refreezing thawed competent cells will result in decreased transformation efficiencies.
2. Mix cells by gently tapping (flicking) the reaction tube. Do not pipet the cells or vortex the tube.
3. For highest transformation efficiencies, follow the recommended time for the duration of the 42°C heat pulse.

Transformation Protocol

1. Thaw on ice one tube of 50 μl competent cells for each transformation. Gently mix the cells by tapping the tube. Once thawed, cells should be used promptly.
2. For determining the transformation efficiency, add 1 μl (10 ng) of pUC19 control DNA to one tube of competent cells. Gently mix by tapping the tube.

- For experimental DNA, add 1 to 50 ng DNA (in $\leq 5 \mu\text{l}$) to one tube of cells. Gently mix by tapping the tube. For best results, the experimental DNA should be free of protein, detergents, organic solvents, and salts.
- Incubate cell/DNA mix on ice for 30 min.
- Heat shock the cells by placing tubes in a 42°C water bath for 30 sec.
- Return the tubes to ice for 2 min.
- Add $250 \mu\text{l}$ room temperature SOC medium to each transformation reaction.
- Secure the tubes in a microcentrifuge tube rack with tape.
- Place the rack on its side in a shaking incubator. Incubate at 37°C for 1 hr while shaking at 225–250 rpm.
- Spread up to $200 \mu\text{l}$ of the transformed cells on LB agar plates containing the appropriate antibiotic (for example $100 \mu\text{g}/\text{ml}$ ampicillin for the pUC19 control plasmid and for Profinity eXact™ tag plasmids). Note: Spreading the liquid until the surface of the plate is dry can significantly reduce the number of transformants.
- Incubate plates 12–16 hr at 37°C .
- Transformation efficiency (CFU/ μg) may be calculated as follows:

$$\frac{\text{CFU on plate}}{\text{ng pUC19 added}} \times \frac{1 \times 10^3 \text{ ng}}{\mu\text{g}} \times \frac{\text{Transformation volume}}{\text{volume plated}} \times \text{dilution factor} = \text{CFU}/\mu\text{g}$$

Example:

If a transformation experiment using 1 ng of pUC19 DNA results in 150 colonies after plating $100 \mu\text{l}$ of a 1:50 dilution, transformation efficiency is:

$$\frac{150 \text{ colonies}}{1 \text{ ng}} \times \frac{1 \times 10^3 \text{ ng}}{\mu\text{g}} \times \frac{0.25 \text{ ml}}{0.1 \text{ ml}} \times 50 = 1.88 \times 10^7 \text{ CFU}/\mu\text{g}$$

Related Products

Catalog#	Description
156-3001	Profinity eXact pPAL RIC-Ready Expression Vector Kit
156-3002	Profinity eXact pPAL Supercoiled Expression Vector Kit
156-3008	Profinity eXact Expression and Purification Starter Kit
156-3006	Profinity eXact Mini Spin Purification Starter Kit, 10 pk
156-3005	Profinity eXact Purification Resin, 10 ml
732-4646	Bio-Scale Mini™ Profinity eXact cartridges, 2 x 1 ml
732-4648	Bio-Scale Mini Profinity eXact cartridge, 1 x 5 ml

Legal Information

Academic and Non-Profit Laboratory Assurance Letter

The T7 expression system is based on technology developed at Brookhaven National Laboratory under contract with the U.S. Department of Energy and is the subject of patent applications assigned to Brookhaven Science Associates, LLC. (BSA). BSA will grant a non-exclusive license for use of this technology, including the enclosed materials, based upon the following assurances:

- These materials are to be used for noncommercial research purposes only. A separate license is required for any commercial use, including the use of these materials for research purposes or production purposes by any commercial entity. Information about commercial licenses may be obtained from the Office of Intellectual Property & Sponsored Research, Brookhaven National Laboratory, Bldg. 185, P.O. Box 5000, Upton, New York 11973-5000; Telephone (631)-344-7134.
- No materials that contain the cloned copy of T7 gene 1, the gene for T7 RNA polymerase, may be distributed further to third parties outside of your laboratory, unless the recipient receives a copy of this license and agrees to be bound by its terms. This limitation applies to strains BL21(DE3), BL21(DE3)pLysS, and BL21(DE3)pLysE, and any derivatives you may make of them.

You may refuse this license by returning the enclosed materials unused. By keeping or using the enclosed materials, you agree to be bound by the terms of this license.

Commercial Entities Doing Business in the United States

A separate license is required for any commercial use, including the use of these materials for research purposes or production purposes by any commercial entity prior to purchasing any T7 products. Please contact your licensing department to confirm that your company already holds a license from BSA or information about commercial licenses may be obtained from the Office of Intellectual Property & Sponsored Research, Brookhaven National Laboratory, Bldg. 185, P.O. Box 5000, Upton, New York 11973-5000; Telephone (631)-344-7134.