

BL21(DE3)pLysS Competent Cells



BL21(DE3)pLysS Competent Cells allow high-efficiency protein expression of any gene that is under the control of a T7 promoter and has a ribosome binding site. BL21(DE3)pLysS is lysogenic for λ -DE3, which contains the T7 bacteriophage gene I, encoding T7 RNA polymerase under the control of the *lac* UV5 promoter. BL21(DE3)pLysS also contains a plasmid, pLysS, which carries the gene encoding T7 lysozyme. T7 lysozyme lowers the background expression level of target genes under the control of the T7 promoter but does not interfere with the level of expression achieved following induction by IPTG. For researchers doi...

Expand to Read More »

Products
Specifications
Product Resources
Patents & Disclaimers



Prices valid for US customers only

PRODUCT	SIZE	CONC.	CATALOG #	*LIST PRICE	ORDER QTY	
Single-Use BL21(DE3)pLysS Competent Cells Protocol MSDS Components	1ml (20 × 50µl)	-	L1195	\$ 190.00	<input type="text"/>	Add to cart 
BL21(DE3)pLysS Competent Cells, >10 ⁶ cfu/µg Protocol MSDS Components	1ml (5 × 200µl)	-	L1191	\$ 155.00	<input type="text"/>	Add to cart 

Storage Conditions

Store at -70°C.

For product intended use please see Patents & Disclaimers tab.

Protocols

[E. coli Competent Cells Technical Bulletin](#)

Related Citations

ATP modulation of Ca²⁺ release by type-2 and type-3 inositol (1, 4, 5)-triphosphate receptors. Differing ATP sensitivities and molecular determinants of action. 2008 *J. Biol. Chem.* **283**, 21579–87.

ATP binding to a unique site in the type-1 S2- inositol 1,4,5-triphosphate receptor defines susceptibility to phosphorylation by protein kinase A. 2006 *J. Biol. Chem.* **281**, 17410-17419.

Effect of species differences of stromelysin-1 (MMP-3) inhibitor potency: An explanation of inhibitor selectivity using homology modeling and chimeric proteins. 1999 *J. Biol. Chem.* **274**, 24881-24887.

Hyperphenylalaninemia with high levels of 7-biopterin is associated with mutations in the PCBD gene encoding the bifunctional protein pterin-4a-carbinolamine dehydratase and transcriptional coactivator (DCoH). 1998 *Am. J. Hum. Genet.* **62**, 1302-1311.

Use Restrictions

L1195, L1191 For Research Use Only. Not for Use in Diagnostic Procedures.

Patents - Disclaimers

L1195, L1191 Usage Restrictions for the T7 Expression System

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 patents assigned to Brookhaven Science Associates, LLC (BSA). This technology, including bacteria, phage and plasmids that carry the gene for T7 RNA polymerase, is to be used for academic or nonprofit laboratory or licensed commercial research purposes only. By accepting or using the T7 expression technology you agree to be bound by the following conditions set forth by BSA. The initial purchaser may refuse to accept the conditions of this notice by returning this product and the enclosed materials to Promega unused.

Academic and NonProfit Laboratories

No materials that contain the cloned gene for T7 RNA polymerase may be distributed further to third parties outside of your laboratory unless the recipient receives a copy of this assurance notice and agrees to be bound by its terms. This limitation applies to Bacterial Strains JM109(DE3), BL21(DE3)pLysS and KRX and to any derivatives thereof.

Commercial Laboratories

A license is required for any commercial use of the T7 expression system, including use of the T7 system for research purposes or for production purposes by any commercial entity. Information about commercial licenses may be obtained from the Licensing Office, Brookhaven National Laboratory, Upton, NY 11973, Telephone: 631-344-7134, FAX: 631-344-3729.

L1195, L1191 Usage restrictions apply to Bacterial Strains JM109(DE3), BL21(DE3)pLysS and KRX and to any derivatives thereof.

[View All Reviews »](#)

[Review This Product »](#)

by

[View All Reviews »](#)

[Review This Product »](#)

by



It appears that you have Javascript disabled. Our website requires Javascript to function correctly.
For the best browsing experience, please enable Javascript.