

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 patents 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 your acceptance of the following assurances:

1. 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; Tel. (631) 344-7134.
2. 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 as listed below, and any derivatives you may make of them.

By keeping or using the enclosed materials, you accept the above assurances and agree to be bound by the terms of this license. Should you not accept the above assurances, you must return the enclosed materials unused.

<i>E. coli</i> B834(DE3)	<i>E. coli</i> Rosetta(DE3)pLysS
<i>E. coli</i> B834(DE3)pLysS	<i>E. coli</i> Rosetta(DE3)pLacI
<i>E. coli</i> BL21(DE3)	<i>E. coli</i> Rosetta 2(DE3)
<i>E. coli</i> BL21(DE3)pLysS	<i>E. coli</i> Rosetta 2(DE3)pLysS
<i>E. coli</i> BL21(DE3)pLysE	<i>E. coli</i> Rosetta 2(DE3)pLacI
<i>E. coli</i> BL26(DE3)pLysE	<i>E. coli</i> RosettaBlue™(DE3)
<i>E. coli</i> BLR(DE3)	<i>E. coli</i> RosettaBlue(DE3)pLysS
<i>E. coli</i> BLR(DE3)pLysS	<i>E. coli</i> RosettaBlue(DE3)pLacI
<i>E. coli</i> HMS174(DE3)	<i>E. coli</i> Rosetta-gami™ 2(DE3)
<i>E. coli</i> HMS174(DE3)pLysS	<i>E. coli</i> Rosetta-gami 2(DE3)pLysS
<i>E. coli</i> HMS174(DE3)pLysE	<i>E. coli</i> Rosetta-gami 2(DE3)pLacI
<i>E. coli</i> NovaBlue(DE3)	<i>E. coli</i> Rosetta-gami B(DE3)
<i>E. coli</i> Origami™ 2(DE3)	<i>E. coli</i> Rosetta-gami B(DE3)pLysS
<i>E. coli</i> Origami 2(DE3)pLysS	<i>E. coli</i> Rosetta-gami B(DE3)pLacI
<i>E. coli</i> Origami 2(DE3)pLacI	<i>E. coli</i> Tuner™(DE3)
<i>E. coli</i> Origami B(DE3)	<i>E. coli</i> Tuner(DE3)pLysS
<i>E. coli</i> Origami B(DE3)pLysS	<i>E. coli</i> Tuner(DE3)pLacI
<i>E. coli</i> Origami B(DE3)pLacI	Bacteriophage λCE6
<i>E. coli</i> Rosetta™(DE3)	Bacteriophage λDE3

© 2009 EMD Chemicals Inc., an affiliate of Merck KGaA, Darmstadt, Germany. All rights reserved. The Novagen® logo and name, Origami™, Rosetta™, RosettaBlue™, Rosetta-gami™, and Tuner™ are trademarks of EMD Chemicals Inc.

USA and Canada

Tel (800) 628-8470
bioscienceshelp@
emdchemicals.com

Europe

France
Freephone
0800 126 461

Germany
Freecall
0800 100 3496

Ireland
Toll Free
1800 409 445

United Kingdom
Freephone
0800 622 935

**All other
European Countries**
+44 115 943 0840

All Other Countries

Contact Your Local Distributor
www.novagen.com
bioscienceshelp@
emdchemicals.com

techservice@merckbio.eu

www.novagen.com

FOR RESEARCH USE ONLY. NOT FOR HUMAN OR DIAGNOSTIC USE.