

Duet Vectors

Table of Contents

| | |
|--|----|
| About the Kits..... | 2 |
| Description | 2 |
| Storage | 2 |
| Multiple Cloning Sites and Fusion Tags | 3 |
| Vector and Host Strain Compatibility | 3 |
| Vector compatibility | 3 |
| Host strain compatibility | 5 |
| Procedures..... | 7 |
| Cloning | 7 |
| Analysis of Duet recombinants | 7 |
| Transformation into expression host strains | 8 |
| Induction | 8 |
| Induction analysis and protein detection, purification, and quantification | 9 |
| References..... | 11 |
| Bacterial Strain Non-distribution Agreement | 12 |

© 2011 EMD Chemicals, Inc., an affiliate of Merck KGaA, Darmstadt, Germany. All rights reserved. Benzonase®, BugBuster®, His•Bind®, PopCulture®, and Novagen® are registered trademarks of Merck KGaA, Darmstadt, Germany. FRETworks™, GST•Tag™, His•Mag™, His•Tag™, LumiBlot™, Lysonase™, Overnight Express™, pACYCDuet™, pCDFDuet™, pCOLADuet™, Perfect Protein™, pETcoco™, pETDuet™, pRSFDuet™, rLysozyme™, RoboPop™, Rosetta™, RosettaBlue™, Rosetta-gami™, Singles™, S•Tag™, Trail Mix™, and Tuner™ are trademarks of Merck KGaA, Darmstadt, Germany.

KOD DNA Polymerases are sold under licensing agreements with F. Hoffman-La Roche, manufactured by TOYOBO, and distributed by EMD Chemicals Inc. KOD XL DNA Polymerase is licensed under U.S. Patent No. 5,436,149 owned by Takara Shuzo, Co., Ltd. Purchase of KOD DNA Polymerases includes an immunity from suit under patents specified in the product inserts to use only the amount purchased for the purchaser's own internal research. No other patents rights are conveyed expressly, by implication, or by estoppel. Further information on purchasing licenses may be obtained by contacting the Director of Licensing, Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California, 94404, USA.

The T7 expression system host strains (DE3 lysogens) are covered by US Patent 5,693,489. For academic and non-profit laboratories, a non-distribution agreement is included with the product. Commercial customers must obtain a license agreement from Brookhaven Science Associates before purchase.

Origami™ 2, Origami™ B, Rosetta™ 2, Rosetta-gami™, Rosetta-gami™ 2, and Rosetta-gami™ B strains are proprietary strains sold under license by EMD Chemicals, Inc.

USA and Canada

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

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.merck4biosciences.com
bioscienceshelp@
emdchemicals.com

techservice@merckbio.eu

www.merck4biosciences.com

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

About the Kits

| | | |
|------------------|-------|---------|
| pACYCDuet™-1 DNA | 10 µg | 71147-3 |
| pETDuet™-1 DNA | 10 µg | 71146-3 |
| pCDFDuet™-1 DNA | 10 µg | 71340-3 |
| pRSFDuet™-1 DNA | 10 µg | 71341-3 |
| pCOLADuet™-1 DNA | 10 µg | 71406-3 |

Description

Coexpression of multiple target genes in *E. coli* is advantageous for studying protein complexes. Coexpression often achieves optimal yield, solubility, and activity and may protect individual subunits from degradation (1–7). The Duet vectors are T7 promoter expression vectors, each designed to coexpress two target proteins in *E. coli*. The Duet vectors carry compatible replicons and antibiotic resistance markers and may be used together in appropriate host strains to coexpress up to eight proteins. Certain combinations of Duet vectors and pET or pETcoco™ vectors are also compatible for coexpression. The capability of Duet vectors to be cotransformed, propagated, and induced for robust target protein coexpression makes them ideal for the analysis of protein complexes (8, 9).

The Duet vectors are designed with compatible replicons (8–11) and drug resistance genes for effective propagation and maintenance of four plasmids in a single cell. pETDuet-1 carries the ColE1 replicon and *bla* gene (ampicillin resistance), pACYCDuet-1 carries the P15A replicon and *cat* gene (chloramphenicol resistance), and pCDFDuet-1 carries the CloDF13 replicon (12) and *aadA* gene (streptomycin/spectinomycin resistance). Two kanamycin-resistant Duet vectors are available; pRSFDuet-1 carries the RSF1030 replicon (13, 14) and pCOLADuet-1 carries the ColA replicon (15). Each vector carries two expression units each controlled by a T7lac promoter for high-level protein expression. Each promoter is followed by a ribosome binding site and multiple cloning site (MCS) region. A T7 terminator follows the second MCS. The multiple cloning regions have restriction sites that facilitate the cloning of two genes and the transfer from other Novagen pET constructs. The Duet vectors provide the option of producing native unfused proteins, or fusions to His•Tag® and S•Tag™ sequences for detection and purification of protein complexes.

Storage

Store DNA at –20°C or –70°C.

Multiple Cloning Sites and Fusion Tags

The Duet Vectors have two *T7lac* promoters, two MCS regions, and a single T7 terminator for the cloning and expression of two target open reading frames (ORFs). The plasmids also carry the *lacI* gene to ensure the expression of sufficient *lac* repressor to control basal expression. In all vectors, MCS1 has an *Nco* I restriction site at the ATG (Met) translation initiation site which can be used to produce unfused protein and has several restriction sites common to most pET vectors (*Bam*H I, *Eco*R I, *Sac* I, *Sal* I, *Hind* III, and *Not* I) for easy transfer of clones. MCS1 also encodes an amino-terminal 6-amino acid (aa) His•Tag[®] sequence for detection and purification. In all vectors, MCS2 has an *Nde* I restriction site at the ATG (Met) translation initiation site, which can be used to produce unfused protein and several restriction sites (*Bgl* II, *Mun* I and *Xho* I) that generate overhangs compatible with *Bam*H I, *Eco*R I, and *Sal* I, respectively, for easy transfer of inserts in pET vectors. MCS2 also encodes an optional carboxy-terminal 15-aa S•Tag[™] sequence for detection, purification, and quantification. The design of MCS regions facilitates the generation of two unfused proteins or one fusion protein with an N-terminal His•Tag, and/or one fusion protein with a C-terminal S•Tag, as desired for detection, purification, or quantification of protein complexes. Both MCS regions include 8-base pair (bp) rare cutting restriction enzymes, *Sse* 8387I and *Not* I in MCS1 and *Fse* I and *Sgf* I in MCS2, to facilitate the insertion of two ORFs into each vector.

Vector and Host Strain Compatibility

Vector compatibility

The vectors differ in their antibiotic resistance markers, replicons, and copy numbers.

Table 1 summarizes the antibiotic-resistant markers of the Duet Vectors.

Table 1 Duet Vector antibiotic resistance markers

| Plasmid | Antibiotic resistance | Marker | Concentration* | Antibiotic Cat. No. |
|---|-------------------------------|-------------|----------------|--|
| pETDuet [™] -1 | ampicillin or carbenicillin | <i>bla</i> | 50 µg/ml | Ampicillin; 171254 Carbenicillin; 69101-3 |
| pACYCDuet [™] -1 | chloramphenicol | <i>cat</i> | 34 µg/ml | 220551 |
| pCDFDuet [™] -1 | streptomycin or spectinomycin | <i>aadA</i> | 50 µg/ml | Streptomycin; 5711 Spectinomycin; 567570 |
| pRSFDuet [™] -1 pCOLADuet [™] -1 | kanamycin | | 30 µg/ml | 420311 |

*When cotransforming four Duet plasmids, the antibiotic concentrations should be reduced by half.

The various replicons carried by the Duet vectors, P15A (pACYCDuet-1), ColE1 (pETDuet-1), CloDF13 (pCDFDuet-1) are compatible with each other (9, 10) and with the replicons carried by the two kanamycin-resistant Duets, COLA (pCOLADuet-1), and RSF1030 (pRSFDuet-1). Duet vectors with different drug resistance markers can be used in combination for coexpression in the appropriate host strains. Duet vectors can also be used with the pET vectors and other constructs that have compatible replicons. Table 2 (page 4) summarizes the replicons and compatible replicons used in Novagen *E. coli* expression vectors and strains. The difference in target gene dosage attributed to plasmid copy number between any of the plasmids could be used to influence relative target protein expression levels.

Note: The combination of a “plain” T7 promoter pET plasmid [i.e., pET-3a-d, pET-20b(+), etc.] with a T7lac promoter plasmid is not recommended.

Determining the optimal combination of Duet vectors for the coexpression of any given set of ORFs is typically an empirical process, however, previous expression results may be useful when choosing expression constructs (9, 11). In general, the target ORFs expressed from pETDuet[™]-1 (ColE1 replicon), pACYCDuet[™]-1 (P15A replicon), and pCDFDuet[™]-1 (CloDF13 replicon) vectors were expressed at higher levels, when used in combination with pCOLADuet[™]-1 (COLA replicon) vector, than when used in combination with pRSFDuet[™]-1 (RSF1030 replicon) vector (11). In one experiment, it was found that in all 2-Duet vector combinations, and in three out of four 3-Duet vector combinations, the ORFs cloned on the higher copy pRSFDuet[™]-1 vector were observed to have the highest expression levels (9). Note, however, that expression of ORFs cloned into pRSFDuet-1 were substantially reduced in coexpression

experiments with pETDuet-1 and pACYCDuet-1 in the same cell. A significantly different pattern of expression was obtained when the lower copy number pCOLADuet™-1 was substituted for pRSFDuet-1 in an otherwise identical set of experiments (11). In this experiment, the expression of the ORFs encoded by pCOLADuet-1 were not substantially reduced when used in combination with both pACYCDuet-1 and pETDuet-1.

Table 2 Plasmid replicons and compatibility

| Plasmid(s) | Replicon (source) | Copy number* | Compatible Replicons |
|--|---------------------------------|--------------------|---|
| pET (all), pETDuet-1 | ColE1 (pBR322) | ~40 | P15A, Mini-F/RK2, CloDF13, RSF1030, ColA |
| pACYCDuet-1, pLysS, pLysE, pLacI, pRARE, pRARE-2 | P15A (pACYC184) | 10–12 | ColE1, Mini-F/RK2, CloDF13, RSF1030, ColA |
| pCDFDuet™-1, pCDF | CloDF13 | 20–40 | ColE1, P15A, RSF1030, ColA |
| pRSFDuet-1, pRSF | RSF1030 | > 100 | ColE1, P15A, CloDF13 |
| pCOLADuet™-1 | COLA (ColA) | 20–40 | ColE1, P15A, CloDF13, Mini-F/RK2 |
| pETcoco™ (all) | Mini-F/RK2 (pBeloBAC11, RK2) | amplifiable to ~40 | ColE1, P15A, ColA |

* Copy number was estimated based on gel analysis (9, 16)

Host strain compatibility

For protein production, the Duet recombinants are transferred to an *E. coli* expression host (DE3) containing a chromosomal copy of the gene for T7 RNA polymerase. The choice of expression host strain is based on strain characteristics and expression vector compatibility. Review the Competent Cells Protocol (User Protocol TB009) for complete descriptions of the host strain characteristics. Use the following tables to determine compatibility. Compatible vectors and host strains are listed in Table 3 below. For expression host strain group, see Table 4 (page 6). For compatible combinations with pETcoco™ plasmid, please consult User Protocol TB333.

Note: The pETcoco vectors are not compatible with pCDFDuet-1 or pRSFDuet-1.

| Table 3 Vector and host strain compatibility | | | | | |
|---|--|---|--------------------------------|--|---|
| Compatible Vector Combinations | | | | Number of coexpressed target proteins | Compatible expression host strains |
| Vector 1 | Vector 2 | Vector 3 | Vector 4* | | |
| pETDuet™-1 (Amp ^R) | pACYCDuet™-1 (Cam ^R) | pRSFDuet™-1 or pCOLADuet™-1 (Kan ^R) | pCDFDuet™-1 (Sm ^R) | 8 | Group A |
| pETDuet-1 (Amp ^R) | pRSFDuet-1 or pCOLADuet-1 (Kan ^R) | pCDFDuet-1 (Sm ^R) | | 6 | Group C |
| pETDuet-1 (Amp ^R) | pACYCDuet-1 (Cam ^R) | pRSFDuet-1 or pCOLADuet-1 (Kan ^R) | | 6 | Group A |
| pETDuet-1 (Amp ^R) | pACYCDuet-1 (Cam ^R) | pCDFDuet-1 (Sm ^R) | | 6 | Group B |
| pRSFDuet-1 or pCOLADuet-1 (Kan ^R) | pCDFDuet-1 (Sm ^R) | pACYCDuet-1 (Cam ^R) | | 6 | Group A |
| pETDuet-1 (Amp ^R) | pRSFDuet-1 or pCOLADuet-1 (Kan ^R) | | | 4 | Group C |
| pETDuet-1 (Amp ^R) | pCDFDuet-1 (Sm ^R) | | | 4 | Group D |
| pETDuet-1 (Amp ^R) | pACYCDuet-1 (Cam ^R) | | | 4 | Group B |
| pRSFDuet™-1 or pCOLADuet-1 (Kan ^R) | pCDFDuet-1 (Sm ^R) | | | 4 | Group C |
| pACYCDuet-1 (Cam ^R) | pRSFDuet™-1 or pCOLADuet-1 (Kan ^R) | | | 4 | Group A |
| pACYCDuet-1 (Cam ^R) | pCDFDuet-1 (Sm ^R) | | | 4 | Group B |

Amp; ampicillin/carbenicillin, 50 µg/ml; Kan; kanamycin, 30 µg/ml; Cam; chloramphenicol, 34 µg/ml;

Sm; streptomycin/spectinomycin, 50 µg/ml

*When cotransforming four Duet plasmids, the antibiotic concentrations should be reduced by half.

Table 4 Vector and Host Strain Compatibility

| Vector | Compatible expression host strains |
|---------------------------|---|
| pETDuet-1 | Group D |
| pACYCDuet-1 | Group B |
| pCDFDuet-1 | Group D |
| pRSFDuet-1 or pCOLADuet-1 | Group C |

Table 4 Strain groups

| Group A | Group B | Group C | Group D |
|---|-----------------------------|---|---------------------------------------|
| B834(DE3) | B834(DE3) | B834(DE3) | B834(DE3) |
| BL21(DE3) | BL21(DE3) | B834(DE3)pLysS ² | B834(DE3)pLysS |
| BLR(DE3) | BLR(DE3) | BL21(DE3) | BL21(DE3) |
| HMS174(DE3) | HMS174(DE3) | BL21(DE3)pLysS ² | BL21(DE3)pLysS |
| NovaBlue(DE3) | NovaBlue(DE3) | BLR(DE3) | BLR(DE3) |
| Origami TM 2(DE3) ¹ | Origami(DE3) ¹ | BLR(DE3)pLysS ² | BLR(DE3)pLysS |
| Tuner TM (DE3) | Origami 2(DE3) ¹ | HMS174(DE3) | HMS174(DE3) |
| | Origami B(DE3) | HMS174(DE3)pLysS ² | HMS174(DE3)pLysS |
| | Tuner(DE3) | NovaBlue(DE3) | NovaBlue(DE3) |
| | | Origami 2(DE3) ¹ | Origami(DE3) ¹ |
| | | Origami 2(DE3)pLysS ^{1, 2} | Origami(DE3)pLysS ¹ |
| | | Rosetta TM (DE3) ² | Origami 2(DE3) ¹ |
| | | Rosetta(DE3)pLysS ² | Origami 2(DE3)pLysS ¹ |
| | | Rosetta 2(DE3) ² | Origami B(DE3) |
| | | Rosetta 2(DE3)pLysS ² | Origami B(DE3)pLysS |
| | | RosettaBlue TM (DE3) ² | Rosetta(DE3) |
| | | RosettaBlue(DE3)pLysS ² | Rosetta(DE3)pLysS |
| | | Rosetta-gami TM 2(DE3) ^{1, 2} | Rosetta 2(DE3) |
| | | Rosetta-gami 2(DE3)pLysS ^{1, 2} | Rosetta 2(DE3)pLysS |
| | | Tuner(DE3) | RosettaBlue(DE3) |
| | | Tuner(DE3)pLysS ² | RosettaBlue(DE3)pLysS |
| | | | Rosetta-gami(DE3) ¹ |
| | | | Rosetta-gami(DE3)pLysS ¹ |
| | | | Rosetta-gami 2(DE3) ¹ |
| | | | Rosetta-gami 2(DE3)pLysS ¹ |
| | | | Rosetta-gami B(DE3) |
| | | | Rosetta-gami B(DE3)pLysS ¹ |
| | | | Tuner(DE3) |
| | | | Tuner(DE3)pLysS |

1.) These strains carry a mutation in ribosomal protein (*rpsL*) conferring resistance to streptomycin; therefore streptomycin is not necessary to maintain strain genotype. If using pCDF vectors, spectinomycin must be used for antibiotic selection because *rpsL* mutation confers streptomycin resistance.

2.) These strains carry a plasmid with a P15A replicon; when used with pETDuetTM-1 and pRSFDuetTM-1 for coexpression substantially reduced expression levels of ORFs cloned into pRSFDuet-1 may result. This reduction in coexpression has not been observed when pCOLADuetTM-1 is substituted for pRSFDuet-1. See page 4 for more information.

Procedures

Cloning

Cloning two ORFs into the same plasmid requires some extra planning. When creating two-ORF constructs, the first insert should lack restriction sites that will be used to insert the second ORF. Typically, the first insert is cloned and an intermediate plasmid is isolated and verified. Then the second ORF is inserted into the remaining MCS to generate the final construct that also requires verification. Unique restriction sites can be added to the second ORF by PCR amplification with primers that contain the desired restriction sites (17, 18). We recommend the use of the robust, high-fidelity KOD HiFi, Hot Start, or XL DNA polymerases, which greatly reduce the chance of generating PCR-based mutations. Standard cloning procedures, including vector and insert preparation and ligation reactions, can be found in the pET System Manual (User Protocol TB055). A high efficiency *recA*⁻, *endA*⁻ host strain such as NovaBlue (Cat. No. 70181) should be used for cloning.

Note: For greatest specificity use KOD Hot Start DNA Polymerase (Cat. No. 71086-3). For greatest yield of long complex targets, use KOD XL DNA Polymerase (Cat. No. 71087-3). (See User Protocol TB341 and TB342, respectively).

Analysis of Duet recombinants

Plasmid DNA from candidate recombinants should be verified for the presence of the correct insert and reading frame. Verification should occur prior to cotransformation to isolate and analyze a single plasmid clone. Several methods available for analysis of transformants include colony PCR, plasmid preparation, restriction analysis, sequencing, and *in vitro* transcription and translation. These methods are described in the pET System Manual (User Protocol TB055).

Duet plasmid DNA can be isolated for transformation into expression hosts, restriction mapping, *in vitro* transcription/translation, and sequence analysis. Isolate pETDuetTM-1, pACYCDuetTM-1, pCOLADuetTM-1, and pCDFDuetTM-1 plasmids using low-copy number DNA purification methods, as found in scientific literature. For pRSFDuetTM-1 plasmids, use the high-copy number protocol provided. Purified DNA must be RNase-free. Plasmid DNA isolated using SpinPrepTM Plasmid Kits, or protocols from scientific literature may require an additional phenol:CIAA extraction (1:1; CIAA is 24 parts chloroform, 1 part isoamyl alcohol) to eliminate RNases (described in the pET System Manual). Use the table below to determine an appropriate plasmid preparation kit.

Sequencing primers

The following table lists appropriate primers to use for PCR and sequence analysis. Note that because the Duet vectors have two T7lac promoters each, the T7 Promoter Primer is not appropriate for PCR or sequence analysis.

| MCS | Primer Type | pACYCDuet TM -1, pCDFDuet TM -1, pRSFDuet TM -1, pCOLADuet TM -1 | pETDuet TM -1 |
|-------------|-------------|--|--|
| MCS1 | Sense | pACYCDuetUP1 Primer Cat. No. 71178-3 | pET Upstream Primer Cat. No. 69214-3 |
| | Antisense | DuetDOWN-1 Primer Cat. No. 71179-3 | DuetDOWN-1 Primer Cat. No. 71179-3 |
| MCS2 | Sense | DuetUP2 Primer Cat. No. 71180-3 | DuetUP2 Primer Cat. No. 71180-3 |
| | Antisense | T7 Terminator Primer Cat. No. 69337-3 | T7 Terminator Primer Cat. No. 69337-3 |

Transformation into expression host strains

Follow the protocols provided in User Protocol TB009 (see User Protocol TB333 if cotransforming with pETcoco™ vectors) for the transformation of Duet vectors into competent cells. For transformations into expression strains using supercoiled plasmid, add 1 µl containing 10–40 ng of each plasmid into competent cells. Perform a 1 h outgrowth prior to plating. Plate 10–70 µl of the transformation mixture. When cotransforming four Duet plasmids, plate the entire transformation mixture, using several plates if necessary and reduce antibiotic concentration by half. Note that antibiotics appropriate for all vectors must be included in the plates and media when cotransforming multiple vectors. Use the tables on pages 5–6 to determine which expression hosts are appropriate for any combination of expression vectors.

Induction

After the plasmids are established in a λDE3 lysogen, expression of the target ORF can be induced by using medium prepared with Overnight Express™ Autoinduction System components (19), or by adding IPTG to a conventional medium. Medium produced with Overnight Express components directs high-density cell growth in the absence of induction followed by autoinduction during the overnight incubation (see User Protocol TB383 for more information). If using IPTG for induction, a final concentration of 1 mM IPTG should be added when the cells reach an OD₆₀₀ of 0.6. Induce for 3 h. Follow the induction protocols in the pET System Manual (User Protocol TB055). See User Protocol TB333 if using a pETcoco™ recombinant with pETDuet or pACYCDuet vectors.

Induction analysis and protein detection, purification, and quantification

For recommendations and protocols regarding induction analysis and optimization, and sample preparation, purification, detection, and quantification, review the pET System Manual (User Protocol TB055) and the following Technical Bulletins, as appropriate.

Coexpression experiments may result in different expression levels of target proteins (9, 20). These differences may be due to unique translation rates or unequal copy numbers for the two expression plasmids (21). If dissimilar expression levels were caused by unequal plasmid copy number, cloning the target genes into the same plasmid may alleviate this problem.

| Detection/Assay Tools for Fusion Tags | | | |
|---|-----------------|-------------|--|
| His•Tag® detection | Cat. No. | Size | User Protocol No./Applications |
| His•Tag Monoclonal Antibody | 70796-4 | 3 µg | TB283 immunofluorescence, immunoprecipitation, Western blotting |
| | 70796-3 | 100 µg | |
| His•Tag AP Western Reagents | 70972-3 | 25 blots | TB283 colorimetric detection |
| His•Tag AP LumiBlot™ Reagents | 70973-3 | 25 blots | TB283 chemiluminescent detection |
| His•Tag HRP LumiBlot Reagents | 70974-3 | 25 blots | TB283 chemiluminescent detection |
| S•Tag™ detection | Cat. No. | Size | User Protocol No./Applications |
| S-protein AP Conjugate | 69598-3 | 50 µl | TB097 Western blotting |
| S-protein HRP Conjugate | 69047-3 | 50 µl | TB136 Western blotting |
| Biotinylated S-protein | 69218-3 | 250 µl | Western blotting |
| S-protein FITC Conjugate | 69060-3 | 200 µl | TB143 immunofluorescence |
| S•Tag AP Western Blot Kit | 69213-3 | 25 blots | TB082 colorimetric detection |
| S•Tag AP LumiBlot Kit | 69099-3 | 25 blots | TB164 chemiluminescent detection |
| S•Tag HRP LumiBlot Kit | 69058-3 | 25 blots | TB145 chemiluminescent detection |
| Quantitative assay | Cat. No. | Size | User Protocol No./Sensitivity |
| FRETWorks™ S•Tag Assay Kit | 70724-3 | 100 assays | TB251 fluorescent assay, Limit < 1 fmol |
| | 70724-4 | 1000 assays | |
| S•Tag Rapid Assay Kit | 69212-3 | 100 assays | TB082 Limit 20 fmol |
| Western blot protein markers | Cat. No. | Size | User Protocol No./Size standards |
| Perfect Protein™ Western Markers | 69959-3 | 25 lanes | TB102; 15, 25, 35, 50, 75, 100 and 150 kDa |
| Trail Mix™ Western Markers | 70982-3 | 25 lanes | TB310; 15, 25, 35, 50, 75, 100 and 150 kDa, and 15, 16, 100 kDa prestained markers |
| Extraction reagents | Cat. No. | Size | User Protocol No./Capacity and features |
| BugBuster® Protein Extraction Reagent | 70584-3 | 100 ml | TB245 Use 5 ml/g wet cell paste. Tris-buffered |
| | 70584-4 | 500 ml | |
| BugBuster HT Protein Extraction Reagent | 70922-3 | 100 ml | TB245 Use 5 ml/g wet cell paste. Tris-buffered and premixed with Benzonase® Nuclease |
| | 70922-4 | 500 ml | |
| BugBuster Master Mix | 71456-3 | 100 ml | TB245 Use 5 ml/g wet cell paste. Tris-buffered and premixed with Benzonase and rLysozyme™ Solution |
| | 71456-4 | 500 ml | |
| BugBuster 10X Protein Extraction Reagent | 70921-3 | 10 ml | TB245 Dilute to 1X with choice of buffer and use 5 ml/g wet cell paste |
| | 70921-4 | 50 ml | |
| | 70921-5 | 100 ml | |
| BugBuster (primary amine-free) Extraction Reagent | 70923-3 | 100 ml | TB245 Use 5 ml/g wet cell paste. PIPPS-buffered |
| | 70923-4 | 500 ml | |
| PopCulture® Reagent | 71092-3 | 15 ml | TB323 Use 0.1 volume per ml of culture |
| | 71092-4 | 75 ml | |
| | 71092-5 | 250 ml | |
| rLysozyme Solution | 71110-3 | 300 KU | TB334 and TB323 Use 40 U per ml of culture volume with PopCulture Reagent and 1 KU per ml of BugBuster Reagent |
| | 71110-4 | 1200 KU | |
| | 71110-5 | 6000 KU | |

| Extraction Reagents | Cat. No. | Size | User Protocol No./Capacity and Features |
|---|-------------------------------|----------------------------|--|
| Benzonase [®] Nuclease, Purity > 90% | 70746-3 70746-4 | 10,000 U 2,500 U | TB245, 323, 261; Use 25 U per ml original culture volume with PopCulture [®] and BugBuster [®] Reagent |
| Lysenase [™] Bioprocessing Reagent | 71230-3 71230-4 71230-5 | 0.2 ml 1 ml 5 × 1 ml | TB361 Optimized blend of rLysozyme [™] Solution and Benzonase Nuclease. Use 3 µl per ml lysis buffer |
| His•Tag[®] purification | Cat. No. | Size | User Protocol No./Capacity and Features |
| Ni-NTA His•Bind [®] Resin | 70666-3 70666-4 70666-5 | 10 ml 25 ml 100 ml | TB273 Capacity is 5–10 mg/ml settled resin |
| Ni-NTA Superflow | 70691-3 70691-4 70691-5 | 10 ml 25 ml 100 ml | TB273 Capacity is 5–10 mg/ml settled resin; high flow rates and pressures |
| Ni-NTA Buffer Kit | 70899-3 | | TB273 All buffers for native purification using Ni-NTA His•Bind and Ni-NTA Superflow resins |
| His•Bind Resin | 69670-3 69670-4 69670-5 | 10ml 50 ml 100ml | TB054 Capacity is 8 mg/ml settled resin |
| His•Bind Buffer Kit | 69755-3 | | TB054 All buffers for native purification using His•Bind Resin |
| His•Bind Columns | 70971-3 70971-4 | pkg/5 pkg/25 | TB054 pre-packed, pre-charged; Capacity is 10 mg per column |
| His•Mag [™] Agarose Beads | 71002-3 71002-4 | 2 ml 10 ml | TB054 magnetic agarose beads, pre-charged; Capacity is 5 mg per ml settled beads |
| His•Bind Purification Kit | 70239-3 | | TB054 10 ml His•Bind Resin, Buffers and Chromatography Columns |
| BugBuster Ni-NTA His•Bind Purification Kit | 70751-3 | | TB273 10 ml Ni-NTA His•Bind Resin, BugBuster, Benzonase, and Chromatography Columns |
| BugBuster His•Bind Purification Kit | 70793-3 | | TB054 10 ml His•Bind Resin and Buffer, BugBuster, Benzonase, and Chromatography Columns |
| PopCulture His•Mag Purification Kit | 71114-3 | | TB054 Process 40 × 3 ml cultures purifying up to 375 µg per 3 ml culture |
| RoboPop [™] His•Mag Purification Kit | 71103-3 | | TB327 Purify up to 12 mg per 96 wells |
| RoboPop Ni-NTA His•Bind Kit | 71188-3 | | TB346 Purify up to 96 mg per 96 wells |
| S•Tag[™] purification | Cat. No. | Size | User Protocol No./Capacity and Features |
| S-protein Agarose | 69704-3 69704-4 | 2 ml 5 × 2 ml | TB087, TB160; Purify up to 1 mg per 2 ml settled resin |
| S•Tag Thrombin Purification Kit | 69232-3 | | TB087 Purify and cleave up to 1 mg target protein per kit (2 ml settled resin) |
| S•Tag rEK Purification Kit | 69065-3 | | TB160 Purify and cleave up to 1 mg target protein per kit (2 ml settled resin) |

References

1. Sanchez, A., Trappier, S.G., Stroher, U., Nichol, S.T., Bowen, M.D., and Feldmann, H. (1998) *Virology* **240**, 138.
2. Stebbins, C.E., Kaelin, W.G., Jr., and Pavletich, N.P. (1999) *Science* **284**, 455–461.
3. Shuman, S. (1990) *J. Biol. Chem.* **265**, 11960–11966.
4. Li, C., Schwabe, J.W., Banayo, E., and Evans, R.M. (1997) *Proc. Natl. Acad. Sci. USA* **94**, 2278–2283.
5. Ishiai, M., Sanchez, J.P., Amin, A.A., Murakami, Y., and Hurwitz, J. (1996) *J. Biol. Chem.* **271**, 20868–20878.
6. Henriksen, L.A., Umbricht, C.B., and Wold, M.S. (1994) *J. Biol. Chem.* **269**, 11121–11132.
7. Tan, S. (2001) *Protein Expr. Purif.* **21**, 224–234.
8. Novy, R., Yaeger, K., Held, D., and Mierendorf, R. (2002) *inNovations* **15**, 2–6.
9. Held, D., Yaeger, K., and Novy, R. (2003) *inNovations* **18**, 4–6.
10. Selzer, G., Som, T., Itoh, T., and Tomizawa, J. (1983) *Cell* **32**, 119–129.
11. Held, D., Yaeger, K., and Novy, R. (2004) *inNovations* **19**, 17–19.
12. Nijkamp, H.J., de Lang, R., Stuitje, A.R., van den Elzen, P.J., Veltkamp, E., and van Putten, A.J. (1986) *Plasmid* **16**, 135–160.
13. Som, T. and Tomizawa, J. (1982) *Mol. Gen. Genet.* **187**, 375–383.
14. Cannon, P.M. and Strike, P. (1992) *Plasmid* **27**, 220–230.
15. Zverev, V.V. and Khmel, I.A. (1985) *Plasmid* **14**, 192–199.
16. Sektas, M. and Szybalski, W. (2002) *inNovations* **14**, 6–8.
17. Novy, R. (1999) *inNovations* **10**, 13.
18. Novy, R., Kolb, K., and Mierendorf, R. (1999) *inNovations* **9**, 6–8.
19. Rucker, P., Torti, F.M., and Torti, S.V. (1997) *Protein Eng.* **10**, 967–973.
20. Grabski, A., Drott, D., and Mehler, M. (2003) *inNovations* **16**, 11–13.
21. Tsao, K.L. and Waugh, D.S. (1997) *Protein Expr. Purif.* **11**, 233–240.

Bacterial Strain Non-distribution Agreement

By purchase of the Origami™ 2, Origami™ B, Rosetta™ 2, RosettaBlue™, Rosetta-gami™, Rosetta-Gami™ 2, or Rosetta-Gami™ B host strains and acceptance of the following terms, Novagen grants a limited license to use the Origami 2, Origami B, Rosetta 2, RosettaBlue, Rosetta-Gami 2, or Rosetta-Gami B host strains for the cloning and expression of genes. The intent of this license is not to limit the research use of these materials, but to protect against unauthorized commercial distribution of the strains by third parties.

1. The Origami 2, Origami B, Rosetta 2, RosettaBlue, Rosetta-Gami 2, or Rosetta-Gami B host strains or any derivative therefrom is not to be offered for resale or distributed outside your laboratory.
2. Gene clones and libraries in the Origami 2, Origami B, Rosetta 2, RosettaBlue, Rosetta-Gami 2, or Rosetta-Gami B host strains may be distributed for research purposes only, provided that the recipient acknowledge the foregoing condition.
3. Commercial customers must obtain a research license agreement from Brookhaven Science Associates before purchasing and using DE3 lysogens of host strains Origami 2, Origami B, Rosetta 2, RosettaBlue, Rosetta-Gami 2, or Rosetta-Gami B.

The initial purchaser may refuse to accept the above conditions by returning the kit unopened and the enclosed materials unused. By accepting or using the kit or the enclosed materials, you agree to be bound by the foregoing conditions.