

Dharmacon™ Yeast Genomic Tiling Collection

Cat #YSC4613

Product Description

The Yeast Genomic Tiling Collection consists of over 1,500 glycerol stock cultures each containing a unique clone of a segment of the yeast *S. cerevisiae* genome in an *E. coli*-yeast shuttle vector, designed by Dr. Greg Prelich of Albert Einstein College of Medicine. These plasmid clones represent a unique and virtually complete overlapping clone collection of the entire *S. cerevisiae* genome. This clone collection constitutes a "minimal functional pathway" through the yeast genome and represents ~ 97% of the length of the yeast genome at the physical level, and ~ 95% of the genome at the functional level. The genomic location represented by each insert was identified in collaboration with the Sanger Institute by sequencing the insert-vector junctions and comparing the obtained sequences to the complete yeast genome sequence in the public database. Analysis revealed only ~ 284 kb of the genome missing in 92 gaps (for a 3 kb average gap size). The genome coordinates (such as the known yeast genes) and well position on the microtiter plate of each clone is precisely known and supplied to the customer. This collection was designed specifically for systematic overexpression screens.

The library was constructed by partially digesting prototrophic yeast genomic DNA with *Mbo*I and subcloning into the *Bam*HI sites of the *E. coli*-yeast shuttle vector, pGP564. Since the library was not made by PCR, it contains no PCR-induced mutations. The proteins are untagged and expressed from their endogenous wild-type promoter. The inserts are flanked by attL sites, allowing transfer of inserts to appropriate destination vectors with attR sites by the Gateway reaction.

The pGP564 shuttle vector contains the *LEU2* selectable marker and 2-micron plasmid sequences necessary for maintenance of a high copy number in yeast. The average insert size in this library is approximately 10 kb, with each insert containing an average of 4-5 genes.

The collection is provided as bacterial cultures of *E. coli* (DH10B) in LB broth with an inert growth indicator + 8% glycerol + kanamycin at a concentration of 50 µg/mL.

Open Labs Products

We provide clone resources developed by leading academic laboratories. Many of these resources address the needs of specialized research communities not served by other vendors. In order to provide these as a public resource, we depend on the contributing laboratories for quality control. Therefore, these are distributed in the format provided by the contributing institution with no additional product validation or guarantee. Additional information can be found in the product manual as well as in associated published articles (if available).

Shipping and Storage

Plates are shipped on dry ice and should be stored at -80 °C.

Source Of Yeast Genomic DNA

The parental yeast strain for isolation of genomic DNA was FY4, a MATa prototroph, originated from Dr. Fred Winston's lab (Harvard Medical School Department of Genetics).



Plate Replication Protocol

To allow any CO₂ that may have dissolved into the medium from the dry ice in shipping to dissipate, store plates at -80 °C for at least 48 hours before thawing.

Procedure

Replication of Plates

1. Dispense ~ 160 µL of sterile (low salt) LB medium into 96-well microtiter plates. The LB should be supplemented with 8% glycerol and the appropriate antibiotic.

Prepare Source Plates

1. Remove the foil seals from the source plates. Removing the seals while the source plates are frozen will minimize cross-contamination.
2. Thaw the source plates with the lids on. Wipe away any condensation under the lid.

Replicate

1. Gently place a disposable replicator into the thawed source plate and lightly move the replicator around inside the well to mix the culture. Make sure to scrape the bottom of the plate of the well.
2. Carefully remove the replicator from the source plate and gently place the replicator into the target plate. Lightly move the replicator back and forth in the target plate to transfer cells.
3. Discard the replicator.
4. Place the lids back on the source plates and target plates.
5. Seal the source plates, being mindful to avoid cross contamination.
6. Repeat this process until all plates have been replicated.
7. Return the source plates to the -80 °C freezer.
8. Place the inoculated target plates in a 37 °C incubator. Incubate the plates for 12-24 hours.

Obtaining Clone Information

Individual clone information and plate locations can be found on the USB drive that accompanies the collection.

Alternatively, the datasheet can be downloaded from our website: gelifesciences.com/dharmacon

Reference

Jones, G. M., J. Stalker, *et al.* (2008). A systematic library for comprehensive overexpression screens in *Saccharomyces cerevisiae*. *Nat Methods* 5(3): 239-41.

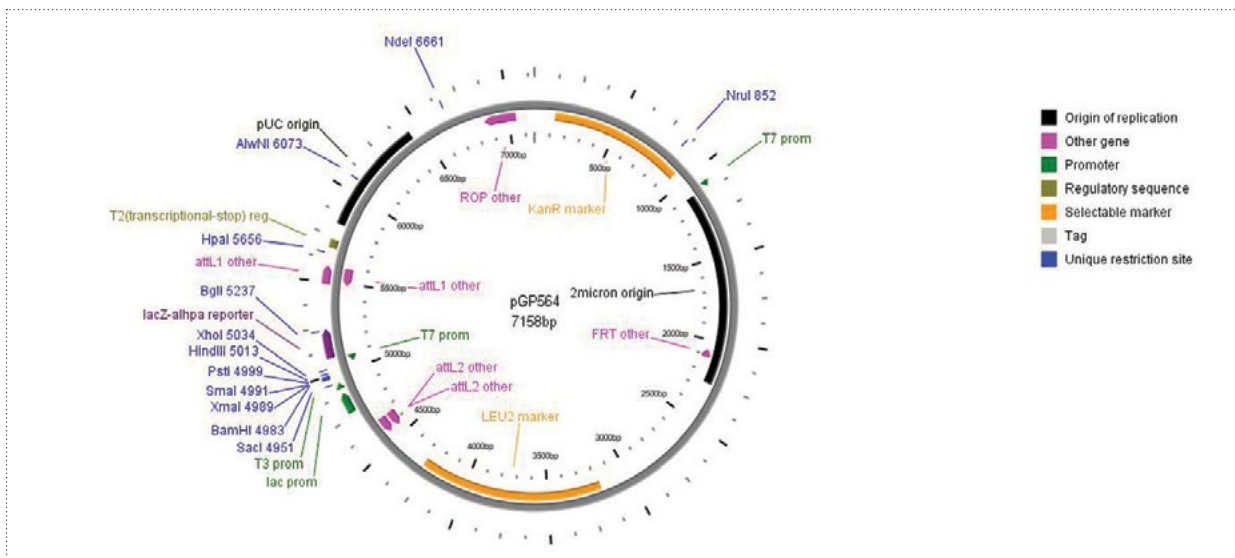


Figure 1. pGP564 Map. The vector is pBR322-based (not pUC-based), so the DNA yield may be a bit lower than a pUC-based vector yield.

Appendix I

Vector Sequence pGP564

LOCUS	7158 bp
FEATURES	<i>Location/Qualifiers</i>
Marker	125..934 /gene="KanR marker"
ORF	125..934 /sequence="ORF_2 rf(6)"
misc_binding	849..854 /dbxref="REBASE:Nrul"
Promoter	1068..1087 /gene="T7 prom"
Rep_Origin	1102..2262 /gene="2micron origin"
Other Gene	2065..2112 /gene="FRT other"
Marker	3175..4281 /gene="LEU2 marker"
ORF	3175..4281 /sequence="ORF_1 rf(4)"
Other Gene	4559..4658 /gene="attL2 other"
Promoter	4774..4889 /gene="lac prom"
Promoter	4915..4934 /gene="T3 prom"
misc_binding	4946..4951 /dbxref="REBASE:Sacl"
misc_binding	4982..4987 /dbxref="REBASE:BamHI"
misc_binding	4988..4993 /dbxref="REBASE:SmaI"
misc_binding	4988..4993 /dbxref="REBASE:XmaI"
misc_binding	4994..4999 /dbxref="REBASE:PstI"
misc_binding	5012..5017 /dbxref="REBASE:HindIII"
misc_binding	5033..5038 /dbxref="REBASE:XhoI"
Promoter	5061..5080 /gene="T7 prom"
Reporter	5088..5246 /gene="lacZ- α reporter"
misc_binding	5230..5240 /dbxref="REBASE:BglI"
Other Gene	5489..5588 /gene="attL1 other"
misc_binding	5653..5658 /dbxref="REBASE:HpaI"
Regulatory_Seq	5687..5730 /gene="T2(transcriptional-stop) reg"
Rep_Origin	5821..6440 /gene="pUC origin"
misc_binding	6067..6075 /dbxref="REBASE:AlwNI"
misc_binding	6659..6664 /dbxref="REBASE:NdeI"
Other Gene	6853..7044 /gene="ROP other"

BASE COUNT

1888 a 1655 c 1613 g 2002 t 0 others

ORIGIN

1	agatctctga	cgctcagtg	aacgacgct	aactcacgtt	aagggat	ggtcatgagc
61	ttgcgccgtc	cogtcaagtc	agcgtaatgc	tctgccagt	ttacaaccaa	ttaaccaatt
121	ctgattagaa	aaactcatcg	agcatcaaat	gaaactgcaa	tttattcata	tcaggattat
181	caataccata	tttttgaaaa	agccgtttct	gtaatgaagg	agaaaactca	ccgaggcagt
241	tccataggat	ggcaagatcc	tggtatcgg	ctgcgattcc	gactcgtcca	acatcaatac
301	aacctattaa	tttccctcg	tcaaaaataa	ggttatcaag	tgagaaatca	ccatgagtga
361	cgactgaatc	cggtgagaat	ggcaaaagtt	tatgcatttc	tttccagact	tgttcaacag
421	gccagccatt	acgctcgtca	tcaaatcac	tcgcatcaac	caaaccgtta	ttcattcgtg
481	attgcgctg	agcgagacga	aatacgcgat	cgctgttaa	aggacaatta	caaacaggaa
541	tcgaatgcaa	ccggcgcagg	aacactgcca	gcgcatcaac	aatattttca	cctgaatcag
601	gatattcttc	taatacctgg	aatgctgttt	ttccgggat	cgcagtggtg	agtaaccatg
661	catcatcagg	agtacggata	aaatgcttga	tggtcggaag	aggcataaat	tccgtcagcc
721	agtttagtct	gaccatctca	tctgtaacat	cattggcaac	gctaccttg	ccatgtttca
781	gaacaactc	tggcgcatcg	ggcttcccat	acaagcgata	gattgtcgca	cctgattgcc
841	cgacattatc	gcgagcccat	ttatacccat	ataaatcagc	atccatgttg	gaatthaatc
901	gcggcctcga	cgtttccctg	tgaatatggc	tcataacacc	ccttgatta	ctgtttatgt
961	aagcagacag	ttttattggt	catgatgata	tatttttata	ttgtgcaatg	taacatcaga
1021	gattttgaga	cacgggcccag	agctgccagg	aaacagctat	gaccatgtaa	tacgactcac
1081	tataggggat	cgatcctcct	gaacgaagca	tctgtgcttc	attttgtaga	acaaaaatgc
1141	aacgcgagag	cgtaat	tcaaaaaag	aatctgagct	gcatttttac	agaacagaaa
1201	tgcaacgcga	aagcgtat	ttaccaacga	agaatctgtg	cttcattttt	gtaaaaacaaa
1261	aatgcaacgc	gagagccta	atttttcaaa	caaagaatct	gagctgcatt	tttacagaac
1321	agaaatgcaa	cgcgagagcg	ctat	aaacaaagat	ctatacttct	ttttgttct
1381	acaaaaatgc	atcccagag	cgctat	ctaacaagc	atcttagatt	acttttttc
1441	tcctttgtgc	gctctataat	gcagtctct	gataactttt	tgcaactgtg	gtccgttaag
1501	gttagaagaa	ggctactttg	gtgtctat	tctcttccat	aaaaaaagcc	tgactccact
1561	tcccgcgttt	actgattact	agcgaagctg	cggtgcatt	ttttcaagat	aaaggcatcc
1621	ccgattatat	tctataccga	tgtggattgc	gcatactttg	tgaacagaaa	gtgatagcgt
1681	tgatgattct	tcattggtca	gaaattatg	aacggtttct	tctat	ctctatatac
1741	tacgtatagg	aaatgtttac	at	tg	t	aatagttctt
1801	actacaattt	ttttgtctaa	agagtaatac	tagagataaa	cataaaaaat	gtagaggtcg
1861	agtttagatg	caagttcaag	gagcgaag	tgatgggta	ggttatatag	ggatatagca
1921	cagagatata	tagcaaagag	atacttttga	gcaatgtttg	tggaagcgg	attcgcaata
1981	tttttagtagc	tcgttacagt	ccggtgcgtt	tttggttttt	tgaagtgcg	tcttcagagc
2041	gcttttggtt	ttcaaaagcg	ctctgaagtt	cctatacttt	ctagagaata	ggacttcgg
2101	aataggaact	tcaaaagcgtt	tccgaaaacg	agcgttccg	aaaatgcaac	gcgagctgcg
2161	cacatacagc	tcaactgttca	cgctgcacct	atatctgcgt	g	tatatata
2221	catgagaaga	acggcatagt	gogt	g	gtacttatat	gcgtctat
2281	atgtaggatg	aaaggtagtc	tagtacctcc	tgatgata	tccattcca	tgccgggtat
2341	cgatgcttc	cttcagcact	acccttagc	tg	g	ctcaattgga
2401	ttagtctcat	ccttcaatgc	tatcatttcc	ttgatattg	gatcacta	agaaccatt
2461	attatcatga	cattaaccta	taaaaatag	cgatcacga	ggcccttcg	tctgcgcgt
2521	ttcggtagt	acggtgaaaa	cctctgacac	atgcagctcc	cgagagcgg	cacagcttgt
2581	ctgtaagcgg	atgccgggag	cagacaagcc	cgtcagggcg	cgtcagcggg	tg
2641	tg	ggcttaacta	tgccgcatca	gagcagattg	tactgagagt	gcaccatatac
2701	gactacgtcg	taaggccgtt	tctgacagag	t	gagggaaactt	tcaccattat
2761	gggaaatgct	tcaagaaggt	atgacttaa	actccatcaa	atggtcaggt	cattgagtgt
2821	ttttat	ttgtat	tttttttag	agaaaatcct	ccaatatcaa	attaggaatc
2881	gtagtttcat	gattttctgt	tacacctaac	ttttgtgtg	gtgccctcct	cctgtcaat
2941	attaatgtta	aagtgaatt	cttttctct	atcacgttga	gccattagta	tcaatttct

3001	tacctgtatt	cctttactat	cctccttttt	ctccttcttg	ataaatgtat	gtagattgcg
3061	tatatagttt	cgctacccct	atgaacatat	tccattttgt	aatttcgtgt	cgtttctatt
3121	atgaatttca	tttataaagt	ttatgtacaa	atatcataaa	aaaagagaat	ctttttaagc
3181	aaggattttc	ttaacttctt	cgcgacagc	atcaccgact	tcggtggtac	tgttggaacc
3241	acctaaatca	ccagttctga	tacctgcatc	caaaccttt	ttaactgcat	cttcaatggc
3301	cttaccttct	tcaggcaagt	tcaatgacaa	tttcaacatc	attgcagcag	acaagatagt
3361	ggcगतtaggg	tcaaccttat	tctttggcaa	atctggagca	gaaccgtggc	atggttcgta
3421	caaaccaaat	gcggtgttct	tgtctggcaa	agaggccaag	gacgcagatg	gcaacaaacc
3481	caaggaacct	gggataacgg	aggcttcatc	ggagatgata	tcaccaaaca	tgttgctggt
3541	gattataata	ccatttaggt	gggttgggtt	cttaactagg	atcatggcgg	cagaatcaat
3541	gattataata	ccatttaggt	gggttgggtt	cttaactagg	atcatggcgg	cagaatcaat
3601	caattgatgt	tgaaccttca	atgtagggaa	ttcgttcttg	atggtttctt	ccacagtttt
3661	tctccataat	cttgaagagg	ccaaaagatt	agctttatcc	aaggaccaaa	taggcaatgg
3721	tggctcatgt	tgtagggcca	tgaagcggc	cattcttgtg	attctttgca	cttctggaac
3781	gggtgattgt	tcactatccc	aagcgacacc	atcaccatcg	tcttcctttc	tcttaccaaa
3841	gtaaatacct	ccactaatt	ctctgacaac	aacgaagtca	gtacctttag	caaattgtgg
3901	cttgattgga	gataagtcta	aaagagagtc	ggatgcaaag	ttacatggtc	ttaagttggc
3961	gtacaattga	agttctttac	ggatttttag	taaaccttgt	tcaggtctaa	cactaccggg
4021	acccattta	ggaccagcca	cagcacctaa	caaaccgca	tcaaccttct	tggaggcttc
4081	cagcgcctca	tctggaagtg	ggacacctgt	agcatcgata	gcagcaccac	caattaaatg
4141	attttcgaaa	tcgaacttga	cattggaacg	aacatcagaa	atagctttaa	gaaccttaat
4201	ggcttcggct	gtgatttctt	gaccaacgtg	gtcacctggc	aaaacgacga	tcttcttagg
4261	ggcagacata	ggggcagaca	ttagaatggt	atatccttga	aatatatata	tatattgctg
4321	aaatgtaaaa	ggtaagaaaa	gttagaaaagt	aagacgattg	ctaaccacct	attggaaaaa
4381	acaataggtc	cttaaataat	attgtcaact	tcaagtattg	tgatgcaagc	atttagtcat
4441	gaacgcttct	ctattctata	tgaaaagccg	gttccggcct	ctcacctttc	ctttttctcc
4501	caatttttca	gttgaaaaag	gtatatgcgt	cagggcacct	ctaggatcag	ctggatggca
4561	aataatgatt	ttattttgac	tgatagtgac	ctgttcgttg	caacaaattg	ataagcaatg
4621	ctttcttata	atgccaaact	tgtacaagaa	agctgggtcg	gcgcgcccac	ccttgcgcc
4681	aatacgc aaa	ccgcctctcc	ccgcgcgttg	gccgattcat	taatgcagct	ggcacgcagc
4741	gtttcccgc	tggaaagcgg	gcagtgcagc	caacgcaatt	aatgtgagtt	acctactca
4801	ttaggcaccc	caggttttac	actttatgct	tcgggctcct	atgttgtgtg	gaattgtgag
4861	cggataacaa	tttcacacag	gaaacagcta	tgaccatgat	tacgccaagc	gcgcaattaa
4921	ccctactaa	agggaacaaa	agctggagct	ccaccgcggt	ggcggccgct	ctagaactag
4981	tggatccccc	gggctgcagg	aattcgatat	caagcttata	gataccgtcg	acctcgaggg
5041	ggggcccggg	acccaattcg	ccctatagtg	agctgtatta	cgcgcgctca	ctggccgtcg
5101	ttttacaacg	tcgtgactgg	gaaaaccctg	gcgttaccca	acttaatcgc	cttgcagcac
5161	atcccccttt	cgccagctgg	cgtaatagcg	aagaggcccg	caccgatcgc	ccttcccaac
5221	agttgcgcag	cctgaatggc	gaatggcgcg	acgcgcctcg	tagcggcgca	ttaagcgcgg
5281	cgggtgtggg	ggttacgcgc	agcgtgacgg	ctacacttgc	cagcgcctca	gcgcccgtc
5341	ctttcgcttt	cttcccttcc	ttctcgcoca	cgttcgcggg	ctttcccogt	caagctctaa
5401	atcgggggct	ccctttaggg	ttccgattta	gtgctttacg	gcacctcgac	ccccaaaaac
5461	ttgattaggg	tgaagggggc	ggccgcggag	cctgcttttt	tgtacaaaagt	tggcattata
5521	aaaagcatt	gctcatcaat	ttgttgcaac	gaacaggtca	ctatcagtca	aaataaaatc
5581	attatttggg	gcccgagctt	aagactggcc	gtcgttttac	aacgtcgtga	ctgggaaaaac
5641	atccatgcta	gcgttaacgc	gagagttagg	aactgccagg	catcaaataa	aacgaaaggc
5701	tcagtccgaa	gactgggcct	ttcgttttat	ctgttgtttg	tcggtgaacg	ctctcctgag
5761	taggacaaat	ccgccgggag	cggatttgaa	cgttgtgaag	caagatctac	cccgtagaaa
5821	agatcaaagg	atcttcttga	gatccttttt	ttctgcgcgt	aatctgctgc	ttgcaacaaa
5881	aaaaccacc	gctaccagcg	gtggtttgtt	tgccggatca	agagctacca	actctttttc
5941	cgaaggtaac	tggcttcagc	agagcgcaga	taccaaatac	tgtccttcta	gtgtagccgt
6001	agttagggcca	ccacttcaag	aactctgtag	caccgcctac	atacctcgct	ctgctaatcc
6061	tgttaccagt	ggctgctgcc	agtggcgata	agtcgtgtct	taccgggttg	gactcaagac

6121	gatagttacc	ggataaggcg	cagcggtcgg	gctgaacggg	gggttcgtgc	acacagccca
6181	gcttgagcgc	aacgacctac	accgaactga	gataacctaca	gcgtgagcta	tgagaaagcg
6241	ccacgcttcc	cgaagggaga	aaggcggaca	ggtatccggt	aagcggcagg	gtcgggaacag
6301	gagagcgcac	gagggagcctt	ccagggggaa	acgcctggta	tctttatagt	cctgtcgggt
6361	ttcgccacct	ctgacttgag	cgctgatttt	tgtgatgctc	gtcagggggg	cgagacctat
6421	ggaaaaacgc	cagcaacgcg	gcctttttac	ggttcctggc	cttttgctgg	ccttttgctc
6481	acatgttctt	tctcgcgtta	tcccctgatt	ctgtggataa	ccgtattacc	gcctttgagt
6541	gagctgatac	cgctcgccgc	agccgaacga	ccgagcgcag	cgagtcagtg	agcgagggaag
6601	cggaagagcg	cctgatgcgg	tattttctcc	ttacgcatct	gtgcggtatt	tcacaccgca
6661	tatggtgcac	tctcagtaca	atctgctctg	atgccgcata	gttaagccag	tataactcc
6721	gctatcgcta	cgtagctggg	tcattgctgc	gccccgacac	ccgccaacac	ccgctgacgc
6781	gccctgacgg	gcttgtctgc	tcccggcatc	cgcttacaga	caagctgtga	ccgtctccgg
6841	gagctgcatg	tgtcagaggt	tttcaccgtc	atcaccgaaa	cgcgcgaggc	agctgcggta
6901	aagctcatca	gcgtggtcgt	gaagcgattc	acagatgtct	gcctgttcat	ccgctccag
6961	ctcgttgagt	ttctccagaa	gcgttaatgt	ctggcttctg	ataaagcggg	ccatgttaag
7021	ggcggttttt	tctgttttgg	tcaactgatgc	ctccgtgtaa	gggggatttc	tgttcatggg
7081	ggtaatgata	ccgatgaac	gagagaggat	gctcacgata	cgggttactg	atgatgaaca
7141	tgcccggtta	ctggaacg				
	//					

FAQS/Troubleshooting

For answers to questions that are not addressed here, please email technical support ts.dharmacon@ge.com

Label licenses:

The gene editing Products, use and applications, are covered by pending and issued patents. Certain Label licenses govern the use of the products, these can be found at gelifesciences.com/webapp/wcs/stores/servlet/catalog/en/GELifeSciences-us/about-us/licensing-statements. It is each Buyer's responsibility to determine which intellectual property rights held by third parties may restrict the use of Products for a particular application.

GE, imagination at work and GE monogram are trademarks of General Electric Company. Dharmacon, Inc., a General Electric company doing business as GE Healthcare. All other trademarks are the property of General Electric Company or one of its subsidiaries. ©2016 General Electric Company—All rights reserved. Version published June 2016. GE Healthcare UK Limited, Amersham Place, Little Chalfont, Buckinghamshire, HP7 9NA, UK



Orders can be placed at:
gelifesciences.com/dharmacon

Customer Support: cs.dharmacon@ge.com
 Technical Support: ts.dharmacon@ge.com or
 1.800.235.9880; 303.604.9499 if you have any questions.