

Genetic Markers in E. coli and EndA+ versus EndA- Strains

Symbol	Description	Effect of Mutation
<i>ara-</i> 14	Mutation in arabinose metabolism	Blocks arabinose catabolism.
araD	L-ribulose phosphate 4-epimerase mutation; part of an inducible operon araBAD repressed by L-arabinose	Blocks arabinose catabolism.
argA	N-Acetylglutamate synthase mutation; inhibited by the presence of arginine	Arginine required from growth in minimal media.
cycA	Mutation of a gene product involved in D-alanine, glycine, D-serine and D-cycloserine transport, and an L-alanine carrier	Mutants cannot use D-alanine as a carbon source.
dam	DNA adenine methylase mutation	Blocks methylation of adenine residues in the sequence 5'G"ATC3'.
dapD	Succinyl-diaminopimelate aminotransferase mutation	Reflects impaired synthesis of succinyl CoA. Mutants need to be supplemented with succinate or lysine + methionine.
dcm	DNA cytosine methylase mutation	Blocks methylation of cytosine in the sequence 5'C"CAGG3' or 5'C"CTGG3'.
deoC	Deoxyribose-phosphate aldolase mutation	
deoR	Regulatory gene mutation allowing constitutive expression of genes for deoxyribose synthesis	Allows efficient propagation of large plasmids.
dut1	Mutation of deoxyuridine triphosphatase, which catalyzes the conversion of dUTP to dUMP and PPi	Mutants are impaired in conversion of dUTP to dUMP, leading to higher dUTP pools, which can lead to misincorporation of uracil instead of thymidine. Stable incorporation of dUTP needs mutation in <i>ung</i> gene.
endA1	DNA-specific endonuclease I mutation	Improves quality of plasmid DNA isolations.
galE	Part of the galETK operon that encodes UDP galactose- 4-epimerase	Mutant is more resistant to bacteriophage P1 infection.
galK	Galactokinase mutation	Blocks catabolism of galactose.
galT	Galactose-1-phosphate uridylyltransferase mutation	Blocks catabolism of galactose.
gyrA96	DNA gyrase mutation	Confers resistance to nalidixic acid.
hflA150	Protease mutation that leads to stabilization of cll gene products	Leads to high frequency of lysogeny by λ phages (1).
hflB	Gene encodes a possible protease component	Leads to high frequency of bacteriophage lambda lysogenization.
hsdR	Host DNA restriction and methylation system mutation.	Allows cloning without cleavage of transformed DNA by endogenous restriction
(r _K ⁻ , m _K +)	Restriction minus, modification positive for the <i>E. coli</i> K strain methylation system	endonucleases. DNA prepared from this strain can be used to transform $r_{\mbox{\tiny K}}$ + <i>E. coli</i> strains.
hsdS20	Mutation of specificity determinant for host DNA	Allows cloning without cleavage of transformed DNA by endogenous restriction
$(r_{\rm B}^{-}, m_{\rm B}^{-})$	restriction and methylation system. Restriction minus, modification minus for the <i>E. coli</i> B strain methylation system	endonucleases. DNA prepared from this strain is unmethylated by the <i>hsdS</i> 20 methylases.
lac1 ^q	Overproduction of the <i>lac</i> repressor protein	Leads to high levels of the <i>lac</i> repressor protein, inhibiting transcription from the <i>lac</i> promoter.
lacY	Galactoside permease mutation	Blocks lactose utilization.
<i>lacZ</i> ∆M15	Partial deletion of β-p-galactosidase gene	Allows complementation of β -galactosidase activity by α -complementation sequence in pGEM®-Z Vectors. Allows blue/white selection for recombinant colonies when plated on X-Gal.
leuB	β-isopropylmalate dehydrogenase mutation	Requires leucine for growth on minimal media.
Δ (lon)	Deletion of <i>lon</i> protease	Reduces proteolysis of expressed proteins.
LysS	pLysS plasmid is integrated into the host genome	Strains carrying this plasmid will be tet resistant and produce T7 lysozyme, a natural inhibitor of T7 RNA polymerase, thus lowering background transcription of sequences under the control of the T7 RNA polymerase promoter (2).
mcrA	Mutation in methylcytosine restriction system	Blocks restriction of DNA methylated at the sequence 5'GmCGC3'.
mcrB	Mutation in methylcytosine restriction system	Blocks restriction of DNA methylated at the sequence 5'AGmCT3'.
metB	Cystathionine γ-synthase mutation	Requires methionine for growth on minimal media.
metC	Cystathionine beta-lyase mutation; involved in methionine biosynthesis	Requires methionine for growth on minimal media.
mtl	Mutation in mannitol metabolism	Blocks catabolism of mannitol.
mutS	Methyl-directed mismatch repair mutation	Prevents repair of the newly synthesized, unmethylated strand.
отрТ	Mutation of protease VII, an outer membrane protein	Reduces proteolysis of expressed proteins.
P2	P2 bacteriophage lysogen present in host	λ phages containing the <i>red</i> and <i>gam</i> genes of λ are growth inhibited by P2 lysogens (3).
proA	γ-glutamyl phoshate reductase mutation	proA/argD mutant will not block proline synthesis but will be repressed by arginine. Mutants excrete proline on minimal media and are resistant to proline analogs. proA/ argD/argR triple mutant grows slowly on minimal media + arginine.
proAB	Mutations in proline metabolism	Requires proline for growth in minimal media.
recA1, recA13	Mutation in recombination	Minimizes recombination of introduced DNA with host DNA, increasing stability of inserts. Inserts are more stable in <i>recA</i> 1 than <i>recA</i> 13 hosts.
recB, recC, recD	Exonuclease V mutations. The Rec BCD trimer (exonuclease V) progressively degrades ssDNA and dsDNA in an ATP-dependent manner to form oligonucleotides; implicated in homologous recombination	Reduces general recombination and affects repair of radiation damage. Allows easier propagation of sequences with inverted repeats.



Genetic Markers in E. coli and EndA+ versus EndA- Strains

Genetic Markers in E. coli (continued).

Symbol	Description	Effect of Mutation
recF	Recombination and repair mutation	Mutant cannot repair daughter strand gaps (post-replicational repair).
relA	ppGpp synthetase I mutation.	Allows RNA synthesis in the absence of protein synthesis. The ribosome-associated ppGpp synthetase I can sense ribosomes with an uncharged tRNA in the A site, a sign of starvation.
rha	Use L-rhamnose, a methylpentose	Blocks rhamnose catabolism.
rpsL	Mutation in subunit S12 of 30S ribosome	Confers resistance to streptomycin.
sbcB	Exonuclease I mutation	Allows general recombination in <i>rec</i> BC mutant strains.
strA	Mutant alters ribosome protein S12	Confers resistance to streptomycin.
supB, supC, supG, supL, supM, supN, supO	Suppressor mutations	Suppresses ochre (UAA) and amber (UAG) mutations.
supD, supE, supF	Suppressor mutations	Suppresses amber (UAG) mutations.
thi-1	Mutation in thiamine metabolism	Mutants require thiamine for growth in minimal media.
thr	Threonine biosynthesis mutation	Mutants are obligate threonine auxotrophs.
thyA	Thymidylate synthase; dTTP biosynthesis	Mutants are obligate thymidine auxotrophs.
Tn5	Transposon	Encodes resistance to kanamycin.
Tn10	Transposon	Encodes resistance to tetracycline.
tonA	Mutation in outer membrane protein	Confers resistance to bacteriophage T1.
traD36	Transfer factor mutation	Prevents transfer of F' episome.
trpC	Phosphoribosyl anthranilate isomerase mutation; part of tryptophan biosynthesis pathway	
trpR	trpR aporepressor; regulates the biosynthesis of tryptophan and its transport	
tsx	T6 and colicin K phage receptor; outer membrane protein involved in specific diffusion of nucleosides; transports the antibotic albicidin	Confers resistance to bacteriophage T6 and colicin K.
ung1	Uracil-DNA N-glycosylase mutation	Allows uracil to exist in plasmid DNA.
xyl-5	Mutation in xylose metabolism	Blocks catabolism of xylose.

References

- 1. Hoyt, A. et al. (1982) Cell 31, 565.
- 2. Studier, F.W. (1991) J. Mol. Biol. 219, 37.
- 3. Kaiser, K. and Murray, N. (1985) In: DNA Cloning, Vol. 1, Glover, D., ed., IRL Press Ltd., Oxford, UK.

Importance of EndA+ versus EndA- E. coli Strains.

Endonuclease I is a 12kDa periplasmic protein encoded by the gene endA that degrades double-stranded DNA. The $E.\ coli$ genotype endA1 refers to a mutation in the wildtype endA gene, which produces an inactive form of the nuclease. $E.\ coli$ strains with this mutation are referred to as EndA negative (EndA—); the wildtype is indicated as EndA+. The table below contains a list of EndA— and EndA+ $E.\ coli$ strains. High-quality DNA is easily obtained from both EndA+ and EndA— strains using Promega PureYieldTM and Wizard[®] Plus SV Plasmid Purification Systems. However, the level of endonuclease I produced is strain-dependent, and these systems may not totally exclude endonuclease I from plasmid DNA prepared from very high endonuclease I-producing strains. In general, we recommend the use of EndA— strains whenever possible.

DIE102luoVI 10 Cold®

ABLE® Cins are EndA+)	KW251
ABLE® K	LE392
BL21 (DE3)	MC1061
BMH 71-18	NM522 (all NM series strains are EndA+)
CJ236	P2392
C600	PR700 (all PR series strains are EndA+)
DH12S™	Q358
ES1301	RR1
HB101	TB1
HMS174	TG1
JM83	TKB1
JM101	Y1088 (all Y10 series strains are EndA+)
JM110	

BJ5183lueXL10-Gold®	MM294	
DH1	SK1590	
DH10B	SK1592	
DH20	SK2267	
DH21	SRB	
$DH5\alpha^{TM}$	Stbl2™	
JM103	Stbl4™	
JM105	SURE®	
JM106	TOP10	
JM107	XLO	
JM108	XL1-Blue	
JM109	XL10-Gold®	
KRX		

NANAOA