



pMC1neo and pMC1neo Poly A Vectors

Instruction Manual

Catalog #213201

Revision C0

For Research Use Only. Not for use in diagnostic procedures.

213201-12



Agilent Technologies

LIMITED PRODUCT WARRANTY

This warranty limits our liability to replacement of this product. No other warranties of any kind, express or implied, including without limitation, implied warranties of merchantability or fitness for a particular purpose, are provided by Agilent. Agilent shall have no liability for any direct, indirect, consequential, or incidental damages arising out of the use, the results of use, or the inability to use this product.

ORDERING INFORMATION AND TECHNICAL SERVICES

Email

techservices@agilent.com

World Wide Web

www.genomics.agilent.com

Telephone

Location	Telephone
United States and Canada	800 227 9770
Austria	01 25125 6800
Benelux	02 404 92 22
Denmark	45 70 13 00 30
Finland	010 802 220
France	0810 446 446
Germany	0800 603 1000
Italy	800 012575
Netherlands	020 547 2600
Spain	901 11 68 90
Sweden	08 506 4 8960
Switzerland	0848 8035 60
UK/Ireland	0845 712 5292
All Other Countries	Please visit www.agilent.com/genomics/contactus

CONTENTS

Materials Provided	1
Storage Conditions	1
Vector Sequences	1
Preparation Of Host Cells	1
Preparation of a -80°C Glycerol Stock	2
Vector Maps	3
Preparation of Media and Reagents	4
Reference	4
MSDS Information	4

pMC1neo and pMC1neo Poly A Vectors

MATERIALS PROVIDED

Material Provided	Quantity
pMC1neo	(25 µg)
pMC1neo poly A	(25 µg)
AG1 Strain: <i>recA1, endA1, gyrA96, thi-1, hsdR17, (r_k⁻, m_k⁺), supE44, relA1</i> , (uncharacterized mutation improves transformation efficiency)	0.5 ml

STORAGE CONDITIONS

Vectors: –20°C

AG1 Strain (Bacterial Glycerol Stock): –80°C

VECTOR SEQUENCES

The complete sequence and list of restriction sites for the pMC1neo and pMC1neo Poly A vectors are available at www.genomics.agilent.com.

PREPARATION OF HOST CELLS

The host strain has been sent as a glycerol stock. For the appropriate media and plates, please refer to the following table:

Bacterial strain	Plates for bacterial streak	Media for glycerol stock
AG-1	LB agar	LB agar

On arrival, prepare the following from the glycerol stock:

Note *Do not allow the contents of the vial to thaw. The vials can be stored at –20 or –80°C, but most strains remain viable longer if stored at –80°C.*

1. Revive the stored cells by scraping off splinters of solid ice with a sterile wire loop.
2. Streak the splinters onto an LB agar plate. Restreak the cells fresh each week.

PREPARATION OF A -80°C GLYCEROL STOCK

1. In a sterile 50-ml conical tube, inoculate 10 ml of the appropriate broth with one or two colonies from the plate. Grow the cells to late log phase.
2. Add 4.5 ml of a sterile glycerol–broth solution (5 ml of glycerol + 5 ml of broth) to the bacterial culture from step 1. Mix well.
3. Aliquot into sterile centrifuge tubes (1 ml/ tube).

This preparation may be stored at -20°C for 1–2 years or at -80°C for more than 2 years.

Vector Maps

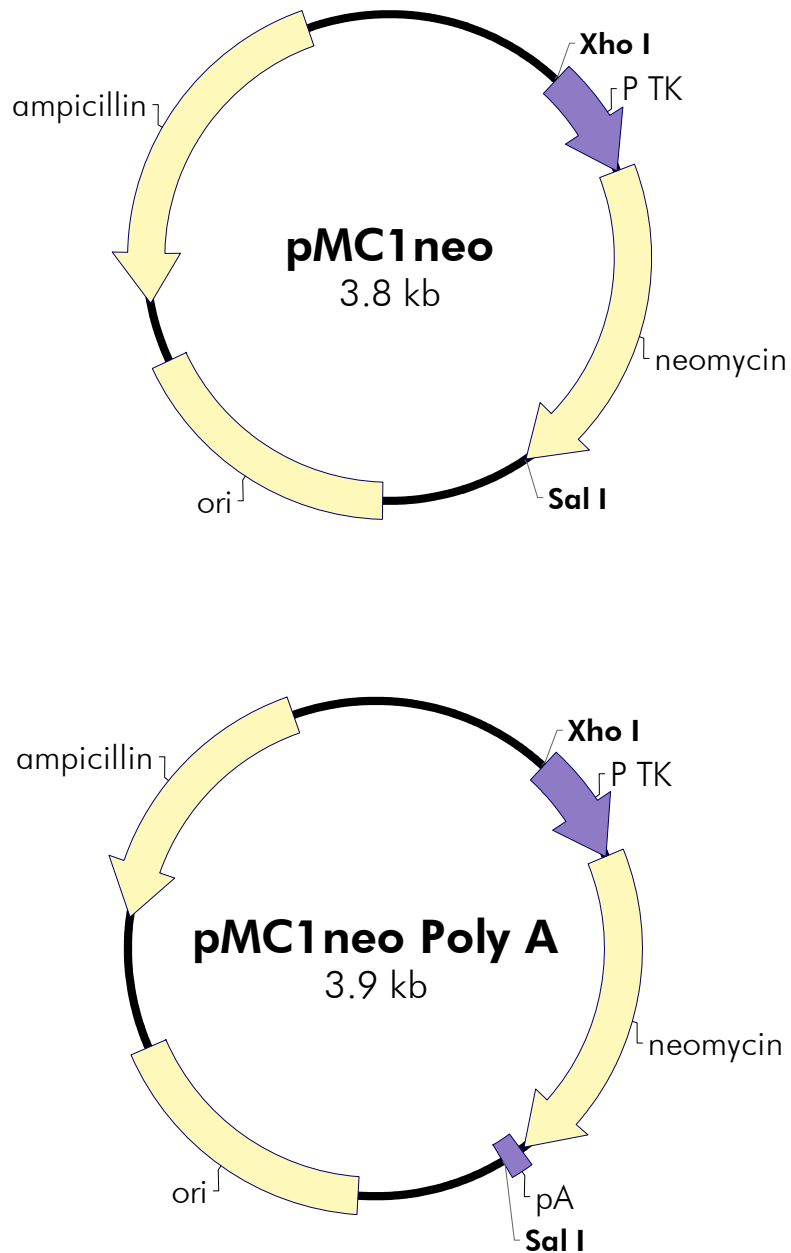


FIGURE 1 Map of the pMC1neo vector (top) and pMC1neo Poly A vector (bottom). The complete sequence and list of restriction sites for the vectors are available at www.genomics.agilent.com.

PREPARATION OF MEDIA AND REAGENTS

LB Broth (per Liter)	LB Agar (per Liter)
10 g of NaCl 10 g of tryptone 5 g of yeast extract Add deionized H ₂ O to a final volume of 1 liter Adjust to pH 7.0 with 5 N NaOH Autoclave	10 g of NaCl 10 g of tryptone 5 g of yeast extract 20 g of agar Add deionized H ₂ O to a final volume of 1 liter Adjust pH to 7.0 with 5 N NaOH Autoclave Pour into petri dishes (~25 ml/100-mm plate)

REFERENCE

1. Thomas, K. R. and Capecchi, M. R. (1987) *Cell* 51(3):503-12.

MSDS INFORMATION

Material Safety Data Sheets (MSDSs) are provided online at <http://www.chem.agilent.com/en-US/search/library/Pages/MSDSSearch.aspx>. MSDS documents are not included with product shipments.