

Product Information

pT7-MAT-Tag™-1 Expression Vector

Catalog Number **E5780**
Storage Temperature $-20\text{ }^{\circ}\text{C}$

TECHNICAL BULLETIN

Product Description

pT7-MAT-Tag-1 is a 4823 bp *Escherichia coli* expression vector used for cytoplasmic expression of a properly inserted open reading frame as a N-terminal MAT-Tag (Metal Affinity Tag) fusion protein. The MAT-Tag (HNHRHKH) is a transition metal binding, e.g. Ni^{+2} and Co^{+2} , sequence useful for high quality purification. The promoter region of the very strong phage T7 promoter^{1,2} drives transcription of MAT-Tag ORF fusion constructs. This vector requires the use of *E. coli* cells containing a source of the T7 polymerase, such as BL21(DE3) cells. Transcription is regulated in these cells by having the T7 polymerase gene under the control of the *lacUV5* promoter. Tighter repression of basal level transcription is provided by the inclusion of *lacO* sequences immediately downstream of the pT7 promoter and having the *lac* repressor gene (*lacI*) on the plasmid.

pT7-MAT-Tag-1 may be used in conjunction with the Director™ Universal PCR System, Catalog Number RDC1, for a simple, rapid and universal method to directionally clone and express PCR products. The MCS has been optimized for use with the *Hind* III and *Bgl* II restriction enzymes often used in the Director system.

N-terminal MAT-Tagged fusion proteins may be purified utilizing the metal affinity properties of the MAT-Tag by using HIS-Select® Nickel Affinity Gel, Catalog Number P6611. Sigma-Aldrich offers a wide selection of related HIS-Select products. Please visit www.sigma-aldrich.com for a complete listing of resins and affinity capture plates.

Reagents

- pT7-MAT-Tag-1 Expression Vector, 10 μg , Catalog Number E4530, 0.5 mg/ml in 10 mM Tris-HCl, pH 8.0, 1 mM EDTA.
- pT7-MAT-Tag-2-BAP Control Vector, 1 μg , Catalog Number C7989, 0.05 mg/ml in 10 mM Tris-HCl, pH 8.0, 1 mM EDTA.

Precautions and Disclaimer

This product is for R&D use only. Not for drug, household or other uses. Consult the MSDS for information regarding hazards and safe handling practices.

Storage/Stability

Product ships on dry ice. Store at $-20\text{ }^{\circ}\text{C}$.

Vector Features

The following table provides map positions to key features in the pT7-MAT-Tag-1 Expression Vector. Sequence verification of the MCS can be performed using the C-24 Sequencing Primer, Catalog Number P7957. The sequence 5'-CTATCATGCCATACCGCGAAAGG-3', available from Sigma-Genosys, is recommended for sequencing through the N-terminal junction.

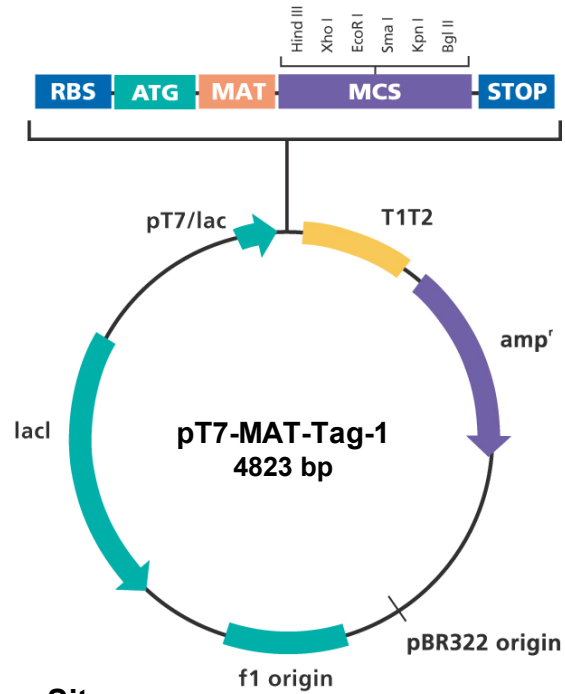
Feature	Map Position
Recommended 5' primer sequence binding site	31-53
pT7 Promoter	72-91
<i>lacO</i>	92-111
Ribosomal Binding Site	143-148
MAT-Tag	161-181
MCS	182-217
C-24 Sequencing Primer Binding Site	252-275
T1/T2 terminator	283-653
β -lactamase (amp^r)	752-1609
pBR322 ori	1817-1936
f1 ori	2600-3063
<i>lacI</i>	3741-4823

References

1. Rosenberg, A. H, et al., Vectors for selective expression of cloned DNAs by T7 RNA polymerase. *Gene*, **56**, 125-135 (1987).

2. Studier, F. W., and Moffatt, B. A., Use of bacteriophage T7 RNA polymerase to direct selective high-level expression of cloned genes. *J. Mol. Biol.*, **189**, 113-130 (1986).

MAT-Tag and HIS-Select are registered trademarks, and Director is a trademark, of Sigma-Aldrich Biotechnology LP and Sigma-Aldrich Co.



Multiple Cloning Site

MAT Sequence									
Met	Ala	His	Asn	His	Arg	His	Lys	His	
ATG	GCT	CAC	AAC	CAC	CGT	CAC	AAA	CAC	
TAC	CGA	GTG	TTG	GTG	GCA	GTG	TTT	GTG	

Hind III	Xho I	EcoR I	Sma I	Kpn I	Bgl II	Stop Codons									
AAG	CTT	CCT	CGA	GTG	AAT	TCC	CCG	GGT	ACC	AGA	TCT	TGA	ATG	AAA	TGA
TTC	GAA	GGA	GCT	CAC	TTA	AGG	GGC	CCA	TGG	TCT	AGA	ACT	TAC	TTT	ACT

Academic and Non-Profit Laboratory Assurance Letter

The T7 system is based on technology developed at Brookhaven National Laboratory under contract with the U.S. Department of Energy and is the subject of U.S. Patent No. 5,693,489 (expiration date, December 2, 2014) assigned to Brookhaven Science Associates, LLC. (BSA). BSA will grant a nonexclusive license for the use of this technology, including the enclosed material, based upon 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 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 Technology Commercialization and Partnerships, Brookhaven National Laboratory, Bldg. 490-C, P.O. Box 5000, Upton, New York 11973-5000, telephone (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 of BL21(DE3), BL21(DE3)pLysS, and BL21(DE3)pLysE, and any derivatives.
3. You may refuse this license by returning the enclosed materials unused. By keeping or using the enclosed materials, you agree to be bound to the terms of this license.