

Product Information

pT7-FLAG™-1 Expression Vector

Catalog Number **P1118**

Storage Temperature -20 °C

TECHNICAL BULLETIN

Product Description

The pT7-FLAG-1 Expression Vector is a 4889 bp *E. coli* expression vector used for cloning and cytoplasmic expression of a properly inserted open reading frame as a N-terminal Met-FLAG fusion protein containing the FLAG® epitope (DYKDDDDK).¹ This vector requires the use of *E. coli* cells containing the DE3 lysogen as a source of the T7 polymerase,² such as BL21(DE3)T1^R cells, Catalog Number B2935.

The N-terminal FLAG fusion protein may be detected using the Monoclonal ANTI-FLAG® M2, Catalog Number F3165, and purified using the ANTI-FLAG M2 Affinity Gel, Catalog Number A2220.

See the chart under pT7-FLAG-1 Features for map positions of the features in the pT7-FLAG-1 vector.

Reagents Provided

- pT7-FLAG-1 Expression Vector
Catalog Number P6742
- pT7-FLAG-1-BAP Control Vector
Catalog Number P6992

Precautions and Disclaimer

This product is for R&D use only, not for drug, household, or other uses. Please consult the Material Safety Data Sheet for information regarding hazards and safe handling practices.

Storage

This product ships on dry ice and storage at -20 °C is recommended.

References

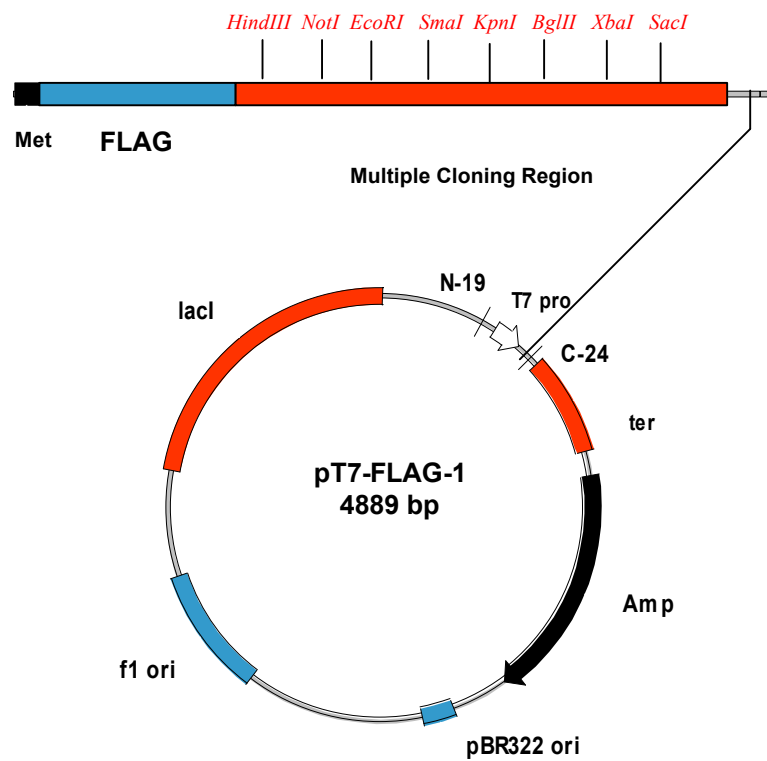
1. Hopp, T.V., et al., *Bio/Technology*, **6**, 1204-1210 (1988).
2. Studier, F.W. et.al., *Journal of Molecular Biology*, **79**, 115-131 (1986).

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. BSA will grant a non-exclusive 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 Economic Development and Technology Transfer, Brookhaven National Laboratory, Bldg. 475D, P.O. Box 5000, Upton, New York 11973-5000, telephone (516)-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(DE3pLysE, and any derivatives you may make of them.

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.

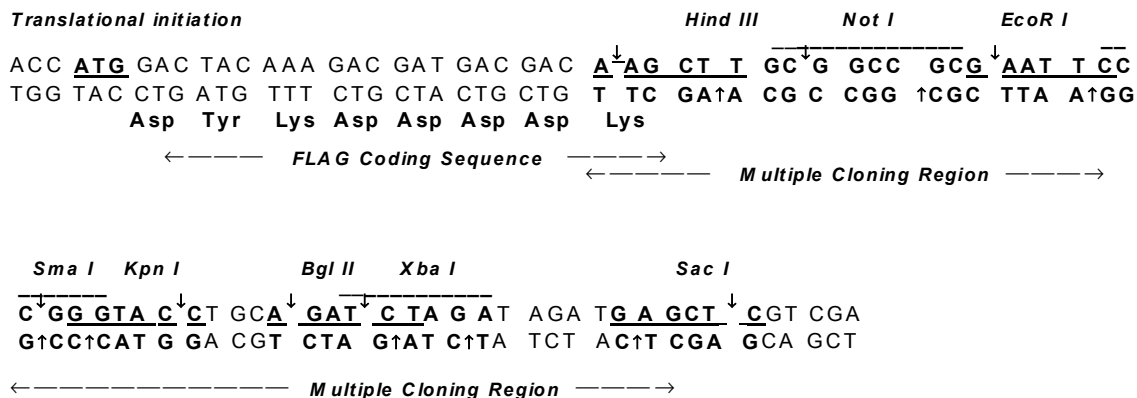


pT7-FLAG-1 Features

Feature	Map Position
pT7 Promoter	436-452
FLAG	521-544
MCS	542-593
T1/T2 ter.	634-1004
Amp	1096-1956
pBR322 ori	2168-2287
f1 ori	2951-3414
lacI	3807-4889

Sequence of the Multiple Cloning Region

Sequence Range: 515 to 598 bp



PT7-FLAG is a trademark, and ANTI-FLAG and FLAG are registered trademarks, of Sigma-Aldrich Biotechnology LP and Sigma-Aldrich Co.

AH,PHC 11/10-1

Sigma brand products are sold through Sigma-Aldrich, Inc.
 Sigma-Aldrich, Inc. warrants that its products conform to the information contained in this and other Sigma-Aldrich publications.
 Purchaser must determine the suitability of the product(s) for their particular use. Additional terms and conditions may apply.
 Please see reverse side of the invoice or packing slip.