

Product Information

pT7-FLAG™-SBP-1 Expression Vector

Catalog Number **P3871**
Storage Temperature $-20\text{ }^{\circ}\text{C}$

TECHNICAL BULLETIN

Product Description

The pT7-FLAG-SBP-1 Expression Vector is a 4931 bp *Escherichia coli* expression vector used for cloning and cytoplasmic expression of a properly inserted open reading frame. The protein is expressed as an N-terminal Met-FLAG®, C-terminal SBP fusion protein containing the FLAG epitope (DYKDDDDK)¹ and the 38 amino acid Streptavidin Binding Peptide (SBP) tag.² This vector requires the use of *E. coli* cells containing the DE3 lysogen as a source of the T7 polymerase,³ such as BL21(DE3)T1^R cells, Catalog Number B2935.

pT7-FLAG-SBP-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 PCR products. The MCS has been optimized for use of the well-validated *Hind* III/*Bgl* II pair of 5' overhang producing restriction enzymes often used in the Director Universal PCR System. The N-terminal FLAG fusion protein may be detected using the ANTI-FLAG® M2 monoclonal antibody, Catalog Number F3165, and purified using the ANTI-FLAG M2 Affinity Gel, Catalog Number A2220. The SBP tag may be detected using Streptavidin-Alkaline Phosphatase, Catalog Number S2890, or Streptavidin-Peroxidase, Catalog Number S5512. The fusion protein may also be purified using Streptavidin-Agarose, Catalog Number S1638. Sigma offers a wide selection of ANTI-FLAG and streptavidin products; please visit www.sigma-aldrich.com for a complete listing of antibody conjugates, resins, and affinity capture plates.

The following table provides map positions to key features in the pT7-FLAG-SBP-1 vector. Sequence verification of the MCS and C-terminal junction 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.

pT7- FLAG-SBP-1 Features

Feature	Map Position
pT7 Promoter	79-98
<i>lacO</i>	99-118
Recommended 5' primer sequence binding site	31-53
Ribosomal Binding Site	150-155
FLAG epitope	165-188
SBP tag	218-337
MCS	186-217
C-24 Sequencing Primer Binding Site	360-383
T1/T2 terminator	391-761
β -lactamase (<i>amp</i> ^r)	860-1717
pBR322 ori	1925-2044
f1 ori	2708-3171
<i>lacI</i>	3849-4931

Reagents

- pT7-FLAG-SBP-1 Expression Vector
Catalog Number E1030
10 μg , 0.5 mg/ml in 10 mM Tris-HCl, pH 8.0, 1 mM EDTA.
- pT7-SBP-1+ BAP Control Vector
Catalog Number C0364
1 μg , 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. Please consult the Material Safety Data Sheet for information regarding hazards and safe handling practices.

Storage/Stability

This product ships on dry ice and storage at $-20\text{ }^{\circ}\text{C}$ is recommended.

References

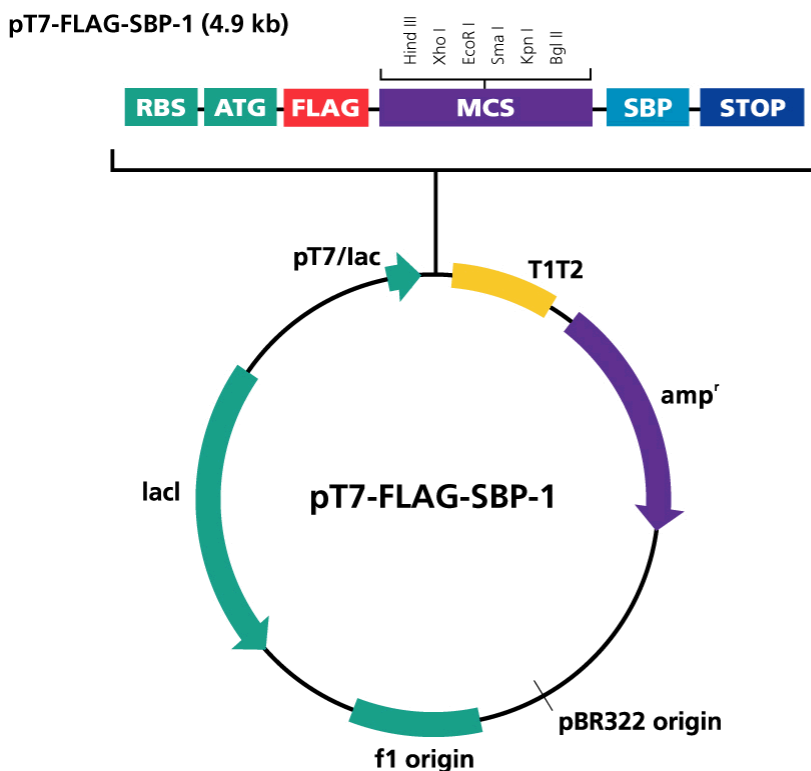
1. Hopp, T. V., et al., *Bio/Technology*, **6**, 1204-1210 (1988).
2. Keefe, A. D., et al., *Protein Expr. Purif.*, **23**, 440-446 (2001).
3. Studier, F. W., et al., *Methods Enzymol.* **185**, 60-89 (1990).

Director and pT7-FLAG are trademarks, and ANTI-FLAG and FLAG are registered trademarks, of Sigma-Aldrich Biotechnology LP and Sigma-Aldrich Co.

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.



Multiple Cloning Site (pT7-FLAG-SBP-1)

FLAG Peptide Sequence																		
Met	Asp	Tyr	Lys	Asp	Asp	Asp	Asp	Lys		Xho I	EcoR I	Sma I	Kpn I	Bgl II				
ATG	GAC	TAC	AAG	GAC	GAT	GAC	GAC	AAG	CTT	CTC	GAG	AAT	TCC	CGG	GTA	CCA	GAT	CT*
TAC	CTG	ATG	TTC	CTG	CTA	CTG	CTG	TTC	GAA	GAG	CTC	TTA	AGG	GCC	CAT	GGT	CTA	GA
										Hind III								

SBP Sequence																		
Met	Asp	Glu	Lys	Thr	Thr	Gly	Trp	Arg	Gly	Gly	His	Val	Val	Glu	Gly	Leu	Ala	Gly
ATG	GAC	GAA	AAA	ACC	ACC	GGT	TGG	CGT	GGT	GGT	CAC	GTT	GTT	GAA	GGT	CTG	GCT	GGT
TAC	CTG	CTT	TTT	TGG	TGG	CCA	ACC	GCA	CCA	CCA	GTG	CAA	CAA	CTT	CCA	GAC	CGA	CCA

SBP Sequence																			
Glu	Leu	Glu	Gln	Leu	Arg	Ala	Arg	Leu	Glu	His	His	Pro	Gln	Gly	Gln	Arg	Glu	Pro	STOP
GAA	CTG	GAA	CAG	CTG	CGT	GCT	CGT	CTG	GAA	CAC	CAC	CCG	CAG	GGT	CAG	CGT	GAA	CCG	TAA
CTT	GAC	CTT	GTC	GAC	GCA	CGA	GCA	GAC	CTT	GTG	GTG	GGC	GTC	CCA	GTC	GCA	CTT	GGC	ATT

*For presentation purposes, the reading frame changes here. Please note that cloning into the *Hind III* and *Bgl II* sites produces an in-frame dual tagged fusion protein.

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.