

Certificate of Analysis

pFC7K (HQ) Flexi® Vector:

Part No.	Size
C854A	20µg

Description: The pFC7K (HQ) Flexi® Vector^(a,b,c) is designed for use with the Carboxy Flexi® System, Transfer (Cat.# C9320). The vector contains a T7 promoter for bacterial or in vitro expression of a protein coding region. The vector is configured to append a C-terminal VSHQHGHQ coding region, which can be used to purify the expressed protein using the MagneHis™, MagZ™, HisLink™ 96 or HisLink™ Protein Purification Systems (Cat.# V8500 and V8550; V8830; V3680 and V3681; and V8821, respectively). The vector contains the lethal barnase gene for positive selection of the insert, a kanamycin-resistance gene for selection of the plasmid and unique SgfI and EcoICRI sites, which allow easy insertion of the sequence of interest. Inserts containing a protein-coding region can be easily transferred to the pFC7K (HQ) Flexi® Vector from other Flexi® Vectors with different expression options. Once inserted in this vector, the sequence is no longer available for transfer because of the absence of a PmeI site at the 3' end of the open reading frame. For more information, see the *Flexi® Vector Systems Technical Manual #TM254*.

Usage Information

Concentration: 100ng/µl.

GenBank® Accession Number: DQ133907.

Storage Buffer: The pFC7K (HQ) Flexi® Vector is supplied in 10mM Tris-HCl (pH 8.0), 1mM EDTA.

Storage Conditions: Store the vector at -20°C. Avoid multiple freeze-thaw cycles and exposure to frequent temperature changes. These fluctuations can greatly alter product stability. See label for expiration date.

Usage Notes: Concentration gradients may form in frozen products and should be dispersed upon thawing. Mix well prior to use.

Quality Control Assays

Nuclease Assay: Following incubation of 1µg of pFC7K (HQ) Flexi® Vector in Restriction Enzyme Buffer B at 37°C for 16 hours, no evidence of nuclease activity is detected by agarose gel electrophoresis.

Physical Purity: $A_{260}/A_{280} > 1.80$.

Restriction Digestion: The presence of unique restriction sites for EcoICRI and SgfI is confirmed by showing that the vector is linearized and yields the expected fragment sizes after digesting 1µg of vector for 2 hours with 10 units of EcoICRI, SgfI and BglII.

^(a)Patent Pending.

^(b)For research use only. Persons wishing to use this product or its derivatives in other fields of use, including without limitation, commercial sale, diagnostics or therapeutics, should contact Promega Corporation for licensing information.

^(c)U.S. Pat. Nos. 8,293,503 and 8,367,403, European Pat. No. 1685247 and other patents and patents pending.

Signed by:

R. Wheeler, Quality Assurance

Part# 9PIC854

Revised 10/16



AF9PIC854 1016C854



Promega

Promega Corporation

2800 Woods Hollow Road	
Madison, WI 53711-5399	USA
Telephone	608-274-4330
Toll Free	800-356-9526
Fax	608-277-2516
Internet	www.promega.com

PRODUCT USE LIMITATIONS, WARRANTY, DISCLAIMER

Promega manufactures products for a number of intended uses. Please refer to the product label for the intended use statements for specific products. Promega products contain chemicals which may be harmful if misused. Due care should be exercised with all Promega products to prevent direct human contact.

Each Promega product is shipped with documentation stating specifications and other technical information. Promega products are warranted to meet or exceed the stated specifications. Promega's sole obligation and the customer's sole remedy is limited to replacement of products free of charge in the event products fail to perform as warranted. Promega makes no other warranty of any kind whatsoever, and SPECIFICALLY DISCLAIMS AND EXCLUDES ALL OTHER WARRANTIES OF ANY KIND OR NATURE WHATSOEVER, DIRECTLY OR INDIRECTLY, EXPRESS OR IMPLIED, INCLUDING, WITHOUT LIMITATION, AS TO THE SUITABILITY, PRODUCTIVITY, DURABILITY, FITNESS FOR A PARTICULAR PURPOSE OR USE, MERCHANTABILITY, CONDITION, OR ANY OTHER MATTER WITH RESPECT TO PROMEGA PRODUCTS. In no event shall Promega be liable for claims for any other damages, whether direct, incidental, foreseeable, consequential, or special (including but not limited to loss of use, revenue or profit), whether based upon warranty, contract, tort (including negligence) or strict liability arising in connection with the sale or the failure of Promega products to perform in accordance with the stated specifications.

© 2005, 2009, 2013, 2014, 2016
Promega Corporation. All Rights Reserved.

Flexi and HaloTag are registered trademarks of Promega Corporation. HisLink, MagneHis and MagZ are trademarks of Promega Corporation.

GenBank is a registered trademark of US Dept of Health and Human Services.

Products may be covered by pending or issued patents or may have certain limitations. Please visit our Web site for more information.

All specifications are subject to change without prior notice.

Product claims are subject to change. Please contact Promega Technical Services or access the Promega online catalog for the most up-to-date information on Promega products.

Part# 9PIC854
Printed in USA. Revised 10/16.

pFC7K (HQ) Flexi® Vector Features and Circle Map

The following features are present in the vector based on nucleotide sequence.

T7 RNA polymerase promoter (-17 to +2)	21-39
SgfI site	61-68
barnase coding region	70-405
EcoICRI site	425-430
HQHGHQ coding region (HQ tag)	430-447
T7 terminator	543-590
Kanamycin-resistance coding region	971-1765
Co/E1-derived plasmid origin of replication	1934-1970
cer site (site for <i>E. coli</i> XerCD recombinase)	2641-2926
rrnB transcription terminator	2977-3378

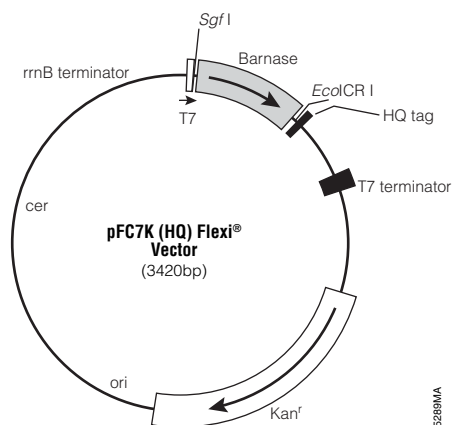


Figure 1. pFC7K (HQ) Flexi® Vector circle map and sequence reference points.

Note: Maps of all the Flexi® Vectors are available at:
www.promega.com/vectors/cloning_vectors.htm

Related Products

Product	Size	Cat. #
Flexi® System, Entry/Transfer	5 entry and 20 transfer reactions	C8640
Flexi® System, Transfer	100 transfer reactions	C8820
Carboxy Flexi® System, Transfer	50 transfer reactions	C9320
10X Flexi® Enzyme Blend (SgfI & PmeI)	25µl	R1851
	100µl	R1852
Carboxy Flexi Enzyme Blend (SgfI & EcoICRI)	50µl	R1901
HaloTag® Flexi® Vectors-CMV Dilution Series Sample Pack	9 × 2µg	G3780
Single Step (KRX) Competent Cells	20 × 50µl	L3002

There are Flexi® Vectors available for many different applications.
 Visit: www.promega.com/applications/cloning to find out more.

Summary of Changes

The following changes were made to the 12/14 revision of this document:
 1. Expired patent or license statements were removed.