

Certificate of Analysis

pFN6K (HQ) Flexi® Vector:

| Part No. | Size |
|----------|------|
| C852A | 20µg |

Description: The pFN6K (HQ) Flexi® Vector^(a,b,c) is designed for use with the Flexi® System, Entry/Transfer (Cat.# C8640), and the Flexi® System, Transfer (Cat.# C8820). The vector contains a T7 promoter for bacterial or in vitro expression of a protein-coding region. The vector appends an N-terminal MKHQHQQAIA 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 PmeI sites, which allow easy insertion or transfer of the sequence of interest. Inserts containing a protein-coding region can be easily transferred from the pFN6K (HQ) Flexi® Vector to other Flexi® Vectors with different expression options. For more information, see the *Flexi® Vector Systems Technical Manual #TM254*.

Usage Information

Concentration: 100ng/µl.

GenBank® Accession Number: DQ132630.

Storage Buffer: The pFN6K (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 pFN6K (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 PmeI 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 PmeI, SgfI and Bgl II.

Part# 9PIC852

Revised 4/18



AF9PIC852 0418C852



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.

© 2005–2018
Promega Corporation. All Rights Reserved.

Flexi and HaloTag are registered 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# 9PIC852
Printed in USA. Revised 4/18.

^(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

pFN6K (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 |
| MKHQHQAIA coding region (HQ tag) | 70-102 |
| Sgfl site | 94-101 |
| barnase coding region | 125-460 |
| PmeI site | 462-469 |
| T7 terminator | 589-636 |
| kanamycin-resistance coding region | 1017-1811 |
| Co/E1-derived plasmid origin of replication | 1980-2016 |
| cer site (site for <i>E. coli</i> XerCD recombinase) | 2687-2972 |
| rrnB transcription terminator | 3023-3424 |

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 (Sgfl & PmeI) | 25µl | R1851 |
| | 100µl | R1852 |
| Carboxy Flexi Enzyme Blend (Sgfl & EcolCRI) | 50µl | R1901 |
| HaloTag® Flexi® Vectors-CMV Dilution Series Sample Pack | 9 × 2µg | G3780 |

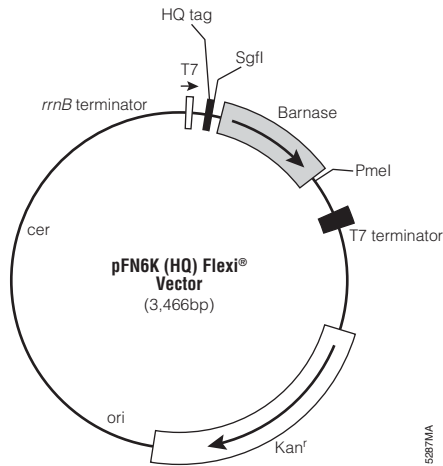


Figure 1. pFN6K (HQ) Flexi® Vector circle map and sequence reference points.

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