

## Certificate of Analysis

### pFN6A (HQ) Flexi® Vector:

Part No.	Size
C851A	20µg

**Description:** The pFN6A (HQ) Flexi® Vector<sup>(a,b,c)</sup> 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, an ampicillin-resistance gene for selection of the plasmid and unique SgfI and Pml sites, which allow easy insertion or transfer of the sequence of interest. Inserts containing a protein-coding region can be easily transferred from the pFN6A (HQ) Flexi® Vector to other Flexi® Vectors with different expression options. For more information, see the *Flexi® Vector Systems Technical Manual #TM254*.

**Concentration:** 100ng/µl.

**GenBank® Accession Number:** DQ132629.

**Storage Buffer:** The pFN6A (HQ) Flexi® Vector is supplied in 10mM Tris-HCl, 1mM EDTA (pH 8.0).

**Storage Conditions:** See the Product Information Label for storage recommendations and expiration date. Avoid multiple freeze-thaw cycles and exposure to frequent temperature changes. These fluctuations can greatly alter product stability.

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

Part# 9PIC851

Revised 03/18



AF9PIC851 0318C851



## 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	<a href="http://www.promega.com">www.promega.com</a>

## Quality Control Assays

**Nuclease Assay:** Following incubation of 1µg of pFN6A (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 Pml 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 Pml, SgfI and BglII.

### 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, 2018  
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# 9PIC851  
Printed in USA. Revised 03/18.

<sup>(a)</sup>Patent Pending.

<sup>(b)</sup>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.

<sup>(c)</sup>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

## Features List and Map for the pFN6A (HQ) Flexi® Vector

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
β-lactamase (Amp <sup>r</sup> ) coding region	970–1830
Co/E1-derived plasmid origin of replication	1985–2021
cer site (site for <i>E. coli</i> XerCD recombinase)	2692–2977
rrnB transcription terminator	3028–3429

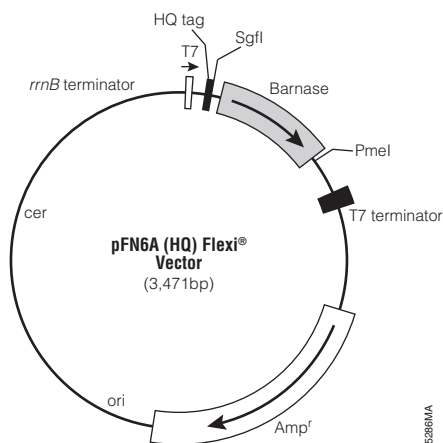


Figure 1. pFN6A (HQ) Flexi® Vector map.

Sequence information and restriction enzyme tables for the Flexi® Vectors are available online at: [www.promega.com/vectors/](http://www.promega.com/vectors/)

## 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 & 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](http://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.