

## Certificate of Analysis

### pFN11A (BIND) Flexi® Vector:

Part No.                      Size (units)  
C934A                              20µg

Part# 9PIC934

Revised 9/16



Instructions for use of this product can be found in the CheckMate™/Flexi® Vector Mammalian Two-Hybrid System Technical Manual #TM283, available online at: [www.promega.com/protocols](http://www.promega.com/protocols)

**Description:** The pFN11A (BIND) Flexi® Vector<sup>(a-d)</sup> is designed for use with the CheckMate™/Flexi® Vector Mammalian Two-Hybrid System (Cat.# C9360), the Flexi® System, Entry/Transfer (Cat.# C8640) and the Flexi® System, Transfer (Cat.# C8820). The pFN11A (BIND) Flexi® Vector contains the yeast GAL4 DNA-binding domain upstream and in-frame with the cloning site. The vector also expresses the *Renilla reniformis* luciferase gene under the control of the SV40 promoter, allowing normalization for differences in transfection efficiency. This vector can be used to test putative transcriptional activation domains for protein sequences of interest when co-transfected with the pGL4.31[*luc2P*/GAL4UAS/Hygro] Vector (Cat.# C9351). In common with other Flexi® Vectors, the pFN11A (BIND) Flexi® Vector contains a CMV immediate early enhancer/promoter region plus a chimeric intron for mammalian expression, a T7 promoter for in vitro expression of the protein-coding region, a barnase gene for positive selection of the insert, an ampicillin resistance gene for selection of the plasmid, and unique *SgfI* and *PmeI* sites that allow easy insertion and transfer of the insert of interest. **Do not use the pFN11A (BIND) Flexi® Vector without an insert as a negative control** because the presence of the barnase gene decreases the viability of the transfected cells. The pFN11A (BIND) Flexi® Vector containing a cloned protein-coding region can be used to transfer the insert to other Flexi® Vectors with different expression options using the Flexi® Systems. For more information, see the *CheckMate™/Flexi® Vector Mammalian Two-Hybrid System Technical Manual #TM283*, and the *Flexi® Vector Systems Technical Manual #TM254*.

**Concentration:** 100ng/µl.

**GenBank® Accession Number:** DQ487212.

**Storage Buffer:** The pFN11A (BIND) Flexi® Vector is supplied in 10mM Tris-HCl, 1mM EDTA (pH 8.0 at 25°C).

**Storage Conditions:** See the Product Information Label for storage recommendations. Avoid multiple freeze-thaw cycles and exposure to frequent temperature changes.

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



AF9PIC934 0916C934



## 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 pFN11A (BIND) 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:**  $A_{260}/A_{280} \geq 1.80$ .

**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 *BglII*.

Signed by:

R. Wheeler, Quality Assurance

### 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.

© 2006–2016 Promega Corporation. All Rights Reserved.

Flexi is a registered trademark of Promega Corporation. CheckMate is a trademark 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.

<sup>(a)</sup>BY USE OF THIS PRODUCT, RESEARCHER AGREES TO BE BOUND BY THE TERMS OF THIS LIMITED USE LABEL LICENSE. If the researcher is not willing to accept the terms of this label license, and the product is unused, Promega will accept return of the unused product and provide the researcher with a full refund.

Researchers may use this product for research use only, no commercial use is allowed. "Commercial use" means any and all uses of this product and derivatives by a party for money or other consideration and may include but is not limited to use in: (1) product manufacture; and (2) to provide a service, information or data; and/or resale of the product or its derivatives, whether or not such product or derivatives are resold for use in research. Researchers shall have no right to modify or otherwise create variations of the nucleotide sequence of the luciferase gene except that researchers may: (1) create fused gene sequences provided that the coding sequence of the resulting luciferase gene has no more than four deoxynucleotides missing at the affected terminus compared to the intact luciferase gene sequence, and (2) insert and remove nucleic acid sequences in splicing research predicated on the inactivation or reconstitution of the luminescence of the encoded luciferase. No other use or transfer of this product or derivatives is authorized without the prior express written consent of Promega. In addition, researchers must either: (1) use luminescent assay reagents purchased from Promega for all determinations of luminescence activity of this product and its derivatives; or (2) contact Promega to obtain a license for use of the product and its derivatives. Researchers may transfer derivatives to others for research use provided that at the time of transfer a copy of this label license is given to the recipients and recipients agree to be bound by the terms of this label license. With respect to any uses outside this label license, including any diagnostic, therapeutic or prophylactic uses, please contact Promega for supply and licensing information. PROMEGA MAKES NO REPRESENTATIONS OR WARRANTIES OF ANY KIND, EITHER EXPRESSED OR IMPLIED, INCLUDING FOR MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE WITH REGARDS TO THE PRODUCT. The terms of this label license shall be governed under the laws of the State of Wisconsin, USA. This label license relates to Promega patents and/or patent applications on improvements to the luciferase gene.

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

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

<sup>(d)</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.

Part# 9PIC934  
Printed in USA. Revised 9/16

## pFN11A (BIND) Flexi® Vector Features and Circle Map

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

CMV immediate/early enhancer/promoter	1–742
Chimeric intron	857–989
T7 RNA polymerase promoter (–17 to +3)	1033–1052
GAL4 fusion protein	1083–1520
IgA linker	1521–1553
SgfI site	1554–1561
Barnase coding region	1585–1920
PmeI site	1922–1929
SV40 late polyadenylation signal	2081–2302
SV40 enhancer/early promoter	2401–2819
hGH intron	2871–3130
<i>Renilla</i> luciferase coding region	3155–4090
Synthetic polyadenylation signal	4148–4194
$\beta$ -lactamase ( <i>Amp<sup>r</sup></i> ) coding region	4457–5317
ColE1-derived plasmid origin of replication	5472–5508
cer region (site for <i>E. coli</i> XerCD recombinase)	6179–6464

**Note:** The IgA linker encodes the polypeptide AIPSTPPTSPAIA.

## Related Products

Product	Size	Cat.#
Flexi® System, Entry/Transfer	5 entry/20 transfer reactions	C8640
Single Step (KRX) Competent Cells	20 × 50µl	L3002

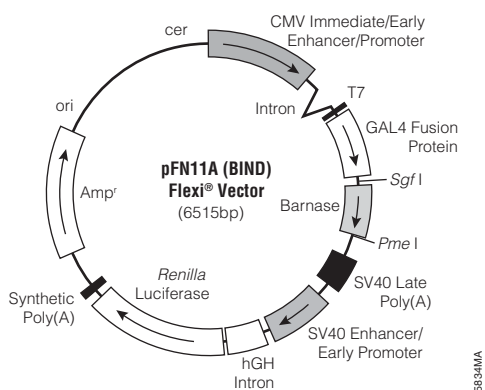


Figure 1. pFN11A (BIND) Flexi® Vector circle map and sequence reference points.

Maps of all the Flexi® Vectors are available at:  
[www.promega.com/vectors/cloning\\_vectors.htm](http://www.promega.com/vectors/cloning_vectors.htm)