

Certificate of Analysis

pFN26A (BIND) *hRluc*-neo Flexi® Vector

Part No. Size
E138A 20µg

Part# 9PIE138
Revised 12/16



Information on the use of this product can be found in the *Flexi® Vector Systems Technical Manual*, #TM254, available at:
www.promega.com/protocols

Description: The pFN26A (BIND) *hRluc*-neo Flexi® Vector^(a-e) (Cat.# E1380) is designed to functionally express a fusion protein comprised of a DNA-binding domain of the yeast GAL4 gene, a linker segment and an in-frame protein-coding sequence flanked by SgfI and PmeI sites at the 5' and 3' ends, respectively, under the control of the human cytomegalovirus (CMV) immediate early promoter. This vector can be used to test putative transcriptional activation domains for protein sequences of interest, such as the ligand binding domain of many nuclear receptors, when cotransfected with the pGL4.35[*Luc2P/9XGAL4* UAS/Hygro] Vector (Cat.# E1370).

Concentration: 100ng/µl.

GenBank® Accession Number: GQ229578.

Storage Buffer: 10mM Tris-HCl, 1mM EDTA (pH 7.4 at 25°C).

Storage Conditions: See the Product Information Label for storage recommendations and expiration date.



AF9PIE138 1216E138



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

Quality Control Assays

Functional Assays

Identity Assay: The vector has been sequenced completely and has 100% identity with the published sequence available at: www.promega.com/vectors

Restriction Digestion: The functional purity of this vector DNA is verified by complete digestion with selected restriction enzymes at 37°C for 1 hour. Samples are examined by agarose gel electrophoresis, and cut and uncut vector DNA are compared with marker DNA.

Contaminant Assays

Contaminating Nucleic Acid Assay: RNA, single-stranded DNA and chromosomal DNA are not evident in a specified sample of this vector as determined by agarose gel electrophoresis.

Nuclease Assay: Following incubation of 1µg of this vector in Restriction Enzyme Buffer at 37°C for 16–24 hours, no evidence of nuclease activity is detected by agarose gel electrophoresis.

Physical Purity: $A_{260}/A_{280} \geq 1.80$, $A_{260}/A_{250} \geq 1.05$.

Signed by:

R. Wheeler, Quality Assurance

(a) 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.

(b) U.S. Pat. No. 8,008,006 and European Pat. No. 1341808.

(c) U.S. Pat. No. 7,906,282 and European Pat. No. 1341808.

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

(e) Patent Pending.

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.

© 2009, 2016 Promega Corporation. All Rights Reserved.

Flexi is a registered trademark of Promega Corporation.

GenBank is a registered trademark of the U.S. Department of Health and Human Services.

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.

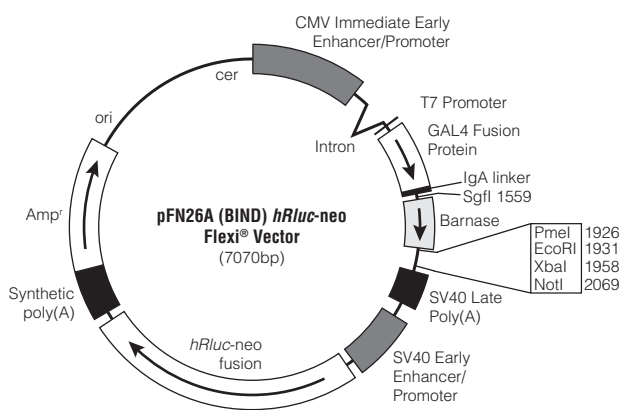
Part# 9PIE138
Printed in USA. Revised 12/16



Promega

pFN26A (BIND) *hRluc*-neo Flexi® Vector Features List and Map:

CMV immediate early enhancer/promoter	1–742
Chimeric intron	857–989
T7 RNA polymerase promoter	1033–1052
GAL4 DNA binding domain fusion protein	1083–1520
IgA linker	1521–1553
Barnase	1585–1920
SV40 late poly(A) region	2081–2302
SV40 early enhancer/promoter	2401–2819
<i>hRluc</i> -neomycin fusion protein	2864–4639
Synthetic poly(A)/transcriptional pause region	4703–5021
Synthetic β-lactamase (Amp ^r) coding region	5012–5872
<i>ColE1</i> -derived plasmid replication origin	6027–6063
Cer	6734–7019



81172MA

pFN26A (BIND) *hRluc*-neo Flexi® Vector Map.

Sequence information is available at: www.promega.com/vectors