

## Certificate of Analysis

### pNL3.2.NF- $\kappa$ B-RE[*NlucP*/NF- $\kappa$ B-RE/Hygro] Vector:

Part No. N111A  
Size 20 $\mu$ g

Part# 9PIN111  
Revised 9/16



Instructions for use of this product can be found in the Nano-Glo<sup>®</sup> Luciferase Assay System Technical Manual #TM369, available online at:  
[www.promega.com/protocols](http://www.promega.com/protocols)

**Description:** The pNL3.2.NF- $\kappa$ B-RE[*NlucP*/NF- $\kappa$ B-RE/Hygro] Vector<sup>(a,b)</sup> contains 5 copies of an NF- $\kappa$ B response element (NF- $\kappa$ B-RE) that drives transcription of a destabilized form of NanoLuc<sup>®</sup> luciferase, an engineered 23.3kDa luciferase fusion protein. The *NlucP* reporter consists of NanoLuc<sup>®</sup> luciferase with a C-terminal fusion to PEST, a protein destabilization domain, which responds more quickly and with greater magnitude to changes in transcriptional activity than unmodified NanoLuc<sup>®</sup> luciferase. The *NlucP* gene is codon optimized for expression in mammalian cells, and all pNL vectors and *Nluc* genes have minimal consensus transcription factor-binding sites to reduce anomalous expression. The vector contains an ampicillin-resistance gene for selection in *E. coli* and a selectable marker for hygromycin resistance in mammalian cells. All forms of NanoLuc<sup>®</sup> luciferases should be used with the optimized substrate, furimazine, found in the Nano-Glo<sup>®</sup> Assay Reagents.

**Concentration:** 1 $\mu$ g/ $\mu$ l.

**GenBank<sup>®</sup> Accession Number:** JQ513377.

**Storage Buffer:** pNL3.2.NF- $\kappa$ B-RE[*NlucP*/NF- $\kappa$ B-RE/Hygro] Vector is supplied in 10mM Tris-HCl (pH 7.4), 1mM EDTA.

**Storage Conditions:** See the product information label for storage recommendations and expiration date.

## Quality Control Assays

### Contaminant Assays

**Contaminating Nucleic Acids:** RNA, single-stranded DNA and chromosomal DNA are not evident in specified quantities of the vector as determined by agarose gel electrophoresis.

**Nuclease Assay:** Following incubation of 1 $\mu$ g of the 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:**  $A_{260}/A_{280} \geq 1.80$ ,  $A_{260}/A_{250} \geq 1.05$  at pH 7.4.

### Functional Assays

**Identity Assay:** The vector was sequenced completely and has 100% identity with the published sequence available at:  
[www.promega.com/vectors](http://www.promega.com/vectors)

**Restriction Digestion:** The functional purity of vector DNA is verified by successful digestion with restriction enzymes at the optimal temperature for one hour. Samples are examined by agarose gel electrophoresis to compare cut and uncut vector DNA with marker DNA.

PLEASE SEE THE BACK OF THIS PAGE  
FOR PROTOCOL INFORMATION.

Signed by:

R. Wheeler, Quality Assurance

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

Researchers may use this product for research use only; no transfer or commercial use of this product is allowed. Commercial use means any and all uses of this product or derivatives by a party in exchange for consideration, including, but not limited to (1) use in further product manufacture; (2) use in provision of services, information or data; and (3) 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 researcher 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 of this product or derivatives is authorized without the prior express written consent of Promega.

#### In addition, researchers must:

(1a) use Nano-Glo<sup>®</sup>-branded luminescent assay reagents (LARs) for all determinations of luminescence activity of this product and its derivatives; or

(1b) contact Promega to obtain a license for use of the product and its derivatives with LARs not manufactured by Promega.

**For users of Nano-Glo<sup>®</sup>-branded LARs intended for energy transfer (such as bioluminescence resonance energy transfer) to acceptors other than a genetically encoded autofluorescent protein, researchers must:**

(2a) use NanoBRET<sup>™</sup>-branded energy acceptors (e.g., BRET-optimized HaloTag<sup>®</sup> ligands) for all determinations of energy transfer activity by this product and its derivatives; or

(2b) contact Promega to obtain a license for use of the product and its derivatives for energy transfer assays to energy acceptors not manufactured by Promega.

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, prophylactic or commercial 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 REGARD TO THE PRODUCT. The terms of this agreement shall be governed under the laws of the State of Wisconsin, USA.

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



AF9PIN111 0916N111



# 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>

### 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.

HaloTag, Nano-Glo and NanoLuc are registered trademarks of Promega Corporation. NanoBRET is a trademark of Promega Corporation.

FuGENE is a registered trademark of Fugent, L.L.C., USA. GenBank is a registered trademark of the U.S. Department of Health and Human Services.

© 2012, 2013, 2016 Promega Corporation. All Rights Reserved.

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

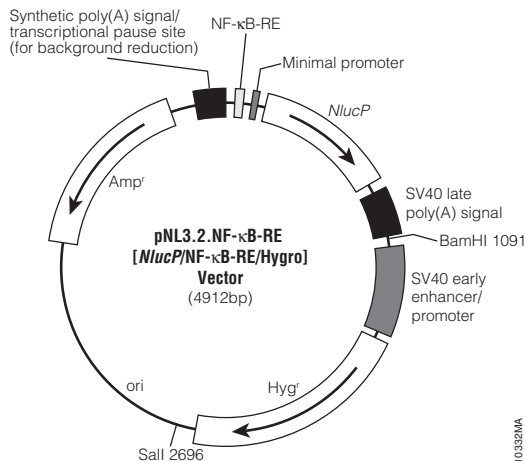
All specifications are subject to change without prior notice.

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# 9PIN111  
Printed in USA. Revised 9/16

## pNL3.2.NF-κB-RE[*NlucP*/NF-κB-RE/Hygro] Vector Features List and Circle Map

NF-κB response elements	33–84
Minimal promoter	117–147
<i>NlucP</i> (NanoLuc®-PEST) reporter gene	180–818
SV40 late poly(A) signal	858–1079
SV40 early enhancer/promoter	1127–1545
Synthetic hygromycin (Hyg <sup>r</sup> ) coding region	1570–2607
Synthetic poly(A) signal	2631–2679
Reporter Vector primer 4 (RVprimer4) binding region	2746–2765
<i>ColE1</i> -derived plasmid replication origin	3003
Synthetic β-lactamase (Amp <sup>r</sup> ) coding region	3794–4654
Synthetic poly(A) signal/transcriptional pause site	4759–4912
Reporter Vector primer 3 (RVprimer3) binding region	4861–4880



**Figure 1.** pNL3.2.NF-κB-RE[*NlucP*/NF-κB-RE/Hygro] Vector map and sequence reference points.

Sequence information and vector maps for the NanoLuc® Vectors are available online at: [www.promega.com/vectors](http://www.promega.com/vectors)

Further information on the use of NanoLuc® Vectors is available in Technical Manual #TM369, which is available online at: [www.promega.com/protocols](http://www.promega.com/protocols)

## Sample Protocol to Determine Induction of Luciferase by TNFα in HEK 293 Cells Transfected with the pNL3.2.NF-κB-RE[*NlucP*/NF-κB-RE/Hygro] Vector

### Materials to be Supplied by User

- Dulbecco's PBS (DPBS)
- 0.05% (w/v) trypsin in DPBS
- DMEM
- DMEM supplemented with 10% fetal bovine serum (DMEM/FBS)
- Tumor necrosis factor-α (Sigma T0157), 10μg/ml solution in PBS containing 1mg/ml BSA
- Nano-Glo® Luciferase Assay System (Cat.# N1110)
- HEK 293 cells
- Transfection reagent (e.g., FuGENE® HD Transfection Reagent, Cat.# E2311)

### Day 1: Plate Cells

1. Grow HEK 293 cells in DMEM/FBS to approximately 75% confluency.
2. Harvest cells via trypsinization: Remove the DMEM/FBS, wash the cells with DPBS and add the trypsin/DPBS (1X volume). After 2 minutes, add a 4X volume of DMEM/FBS, collect the cell suspension and pellet the cells by centrifugation. Aspirate the supernatant and resuspend in DMEM/FBS. We have routinely used a concentration of 15,000 viable cells/100μl DMEM/FBS.
3. Dispense 100μl of the cell suspension into the wells of a 96-well plate. Plate enough wells to perform each test condition in triplicate.
4. Cover the plate and place it in a tissue culture incubator at 37°C overnight (or for 24 hours).

### Day 2: Transfect Cells

1. Transfect the cells using a high-efficiency transfection reagent (e.g., FuGENE® HD Transfection Reagent, Cat.# E2311). Transfection conditions may require optimization.
2. Cover the plate and place it in a tissue culture incubator at 37°C overnight or as needed for cell recovery depending on the transfection method used. We have used 24 hours recovery time for lipid-mediated transfections.

### Day 3: Induce Transfected Cells

1. Prepare 1X induction and 1X control solutions. Calculate the volume of 1X induction and 1X control solution by multiplying the number of wells needed for each solution by 100μl and prepare 110% of this amount.
  - 1X induction solution: Dilute 10μg/ml TNFα solution to 20ng/ml in DMEM/FBS. Final TNFα concentration will be 20ng/ml.
  - 1X control solution: DMEM/FBS.
2. Remove media from wells that will be treated with either 1X induction solution or 1X control solution.
3. Add 100μl of 1X induction solution to the cells to be induced and 100μl of 1X control solution to the control noninduced cells.
4. Return the plate to the tissue culture incubator and induce for 5 hours.
5. Analyze luciferase activity using the Nano-Glo® Luciferase Assay System (Cat.# N1110, Technical Manual #TM369).
6. Calculate the fold induction as follows:

$$\text{Fold Induction} = \frac{\text{Average relative light units of induced cells}}{\text{Average relative light units of control cells}}$$