

Certificate of Analysis

pNLCol2[*luc2*-P2A-*NlucP*/minP/Hygro] Vector:

Part No.	Size
N147A	20µg

Description: The pNLCol2[*luc2*-P2A-*NlucP*/minP/Hygro] Vector^(a,b) is a second generation coincidence reporter vector (Col; 1) that allows expression of both firefly luciferase (Fluc, *luc2* gene) from *Photinus pyralis* and NanoLuc[®] Luciferase fused to a PEST destabilization domain (NanoLuc[®]-PEST, *NlucP* gene) from the same mRNA transcript. The stoichiometric expression of both luciferases is achieved by use of the P2A sequence from porcine teschovirus-1, which promotes a ribosomal skip and expression of the two unfused enzymes with different compound interaction profiles. When used in high-throughput screening, true hits will show a similar response from both reporters, helping to reduce the number of false hits and the workload associated with follow-up screens.

The *luc2* and *NlucP* genes are codon optimized for mammalian expression, and all pNL vectors have minimal consensus transcription factor-binding sites to reduce anomalous expression. The pNLCol2[*luc2*-P2A-*NlucP*/minP/Hygro] Vector contains a multiple cloning region for insertion of promoter elements of choice upstream of a minimal promoter.

The pNLCol2[*luc2*-P2A-*NlucP*/minP/Hygro] Vector contains the following key features:

- Firefly luciferase reporter gene (*luc2*).
- P2A peptide sequence that promotes a ribosomal skip, resulting in the stoichiometric expression of two unfused reporter proteins from the same mRNA transcript.
- NanoLuc[®] Luciferase reporter gene fused to the PEST destabilization domain (*NlucP*). PEST-mediated turnover by the proteasome provides a reporter with a very short intracellular protein half-life, providing a reporter protein that responds more quickly and typically with greater signal-to-background compared to other reporters. NanoLuc[®]-PEST is also substantially brighter than *Renilla* luciferase, providing a dramatic increase in sensitivity for the second generation coincidence reporter system.
- Ampicillin resistance gene for selection in bacteria.
- Hygromycin resistance gene for selection in mammalian cells.
- Multiple cloning region for insertion of promoter elements of choice.
- Minimal promoter

Concentration: 1µg/µl.

GenBank[®] Accession Number: KM359772.

Storage Buffer: The pNLCol2[*luc2*-P2A-*NlucP*/minP/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.

Endotoxin Concentration: Endotoxin Units (EU) are obtained using *Limulus amoebocyte* lysate testing. The specification is <100EU/mg of plasmid DNA.

Nuclease Assay: Following incubation of 1µg 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

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.

Reference

1. Cheng, K.C. and Inglese, J. (2012) A coincidence reporter-gene system for high-throughput screening. *Nat. Methods*, **9**, 937.

Signed by:

R. Wheeler, Quality Assurance

Part# 9PIN147
Revised 9/16



AF9PIN147 0916N147



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.

Dual-Luciferase, HaloTag, NanoBRET, NanoLuc and Nano-Glo are registered trademarks of Promega Corporation.

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

© 2014, 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.

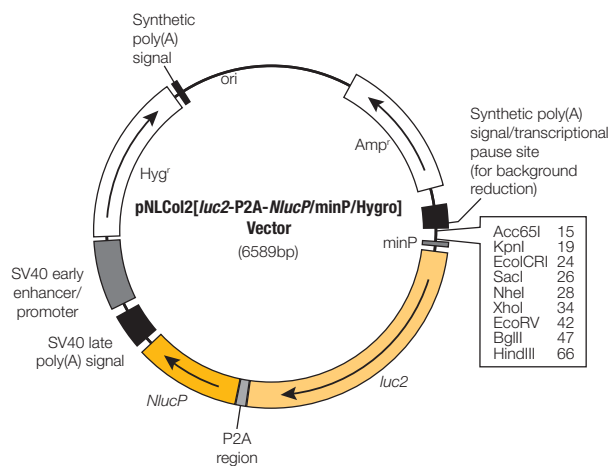
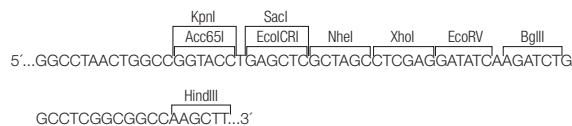
Part# 9PIN147
Printed in USA. Revised 9/16

pNLCol2[*luc2*-P2A-*NlucP*/minP/Hygro] Vector Features and Maps

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

Multiple cloning region	1–70
Minimal promoter (minP)	78–108
<i>luc2</i> reporter gene	141–1790
P2A region	1791–1856
<i>NlucP</i> reporter gene	1857–2495
SV40 late poly(A) region	2535–2756
SV40 early enhancer/promoter	2804–3222
Hygromycin resistance coding region	3247–4284
Synthetic poly(A) sequence	4308–4356
Reporter Vector primer 4 (RVprimer4) binding region	4423–4442
<i>Col</i> /E1-derived plasmid replication origin	4680
Synthetic β-lactamase (<i>Amp^r</i>) coding region	5741–6331
Synthetic poly(A) signal/transcriptional pause region	6436–6589
Reporter Vector primer 3 (RVprimer3) binding region	6538–6557

Multiple cloning region of the pNLCol2[*luc2*-P2A-*NlucP*/minP/Hygro] Vector.



Sequence information and vector maps are available online: www.promega.com/vectors

For information on the Nano-Glo® Dual-Luciferase® Assay System see Technical Manual #TM426, available online: www.promega.com/protocols

^(a) 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 terms of this label license, and the product is unused, Promega will accept return of the unused product and provide researcher with a full refund.

Researcher 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. Researcher shall have no right to modify or otherwise create variations of the nucleotide sequences of the luciferase genes except that researcher may (1) create fused gene sequences provided that the coding sequences of the resulting luciferase genes have no more than four deoxynucleotides missing at the affected terminus compared to the intact luciferase gene sequences; and (2) insert and remove nucleic acid sequences in splicing research predicated on the inactivation or reconstitution of the luminescence of the encoded luciferases. No other use of this product or derivatives is authorized without the prior express written consent of Promega.

In addition, researcher must:

- (1a) use Nano-Glo®-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 uses of Nano-Glo®-branded LARs intended for energy transfer (such as bioluminescence resonance energy transfer) to acceptors other than a genetically encoded autofluorescent protein, researcher must:

- (2a) use NanoBRET®-branded energy acceptors (e.g., BRET-optimized HaloTag® 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.

Researcher 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 label license shall be governed under the laws of the State of Wisconsin, USA.

^(b)The method of recombinant expression of *Coleoptera* luciferase is covered by U.S. Pat. Nos. 5,583,024, 5,674,713 and 5,700,673. A license (from Promega for research reagent products and from The Regents of the University of California for all other fields) is needed for any commercial sale of nucleic acid contained within or derived from this product.