

Certificate of Analysis

pNL3.2.CMV Vector:

Part No.	Size
N141A	20µg

Description: The pNL3.2.CMV Vector^(a,b) is designed for constitutive expression of the NanoLuc[®] reporter fused to a PEST destabilization domain (*NlucP*). The NlucP fusion protein naturally accumulates at low intracellular levels due to constitutive proteosomal degradation, thus serving as a negative control for experiments configured to measure regulated changes in NanoLuc[®] luciferase expression levels. This vector also encodes a hygromycin resistance gene for selection in mammalian cells and an ampicillin resistance gene for selection in bacteria.

The pNL3.2.CMV Vector contains the following features:

- A **CMV promoter** for robust expression in mammalian cells
- The **NanoLuc[®] reporter gene**, for high sensitivity detection of the protein of interest.
- **PEST** sequence for maintaining low intracellular NanoLuc[®] luciferase concentration.
- **Ampicillin resistance** gene for selection in bacteria.
- **Hygromycin resistance** gene for selection in mammalian cells.

Concentration: 1µg/µl.

GenBank[®] Accession Number: KF853603.

Storage Buffer: The pNL3.2.CMV Vector is supplied in 10mM Tris-HCl, 1mM EDTA (pH 7.4).

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. See label for expiration date.

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

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

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 Enzyme Digest: The functional purity of the vector DNA is verified by successful digestion with restriction enzymes at the optimal temperature for one hour. Samples are examined by agarose gel electrophoresis, comparing cut and uncut vector DNA with marker DNA.

Part# 9PIN141

Revised 9/16



AF9PIN141 0916N141



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.

© 2013, 2016 Promega Corporation. All Rights Reserved.

FuGENE, Nano-Glo and NanoLuc are registered trademarks of Promega Corporation. NanoBRET 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

Part# 9PIN141
Printed in USA. Revised 9/16

Signed by:

R. Wheeler, Quality Assurance

pNL3.2.CMV Vector Features

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

CMV immediate early enhancer/promoter	14–755
NanoLuc®-PEST protein coding region	859–1497
SV40 late poly(A) region	1537–1758
SV40 early enhancer/promoter	1806–2224
Synthetic hygromycin (Hyg ^r) coding region	2249–3286
Beta-lactamase (Amp ^r) coding region	4473–5333
ColE1-derived plasmid replication origin	3682–3718
Synthetic poly(A) signal	5438–5486

Related Products

Product	Size	Cat. #
Nano-Glo® Luciferase Assay	10ml	N1110
	100ml	N1120
	10 × 10ml	N1130
	10 × 100ml	N1150
FuGENE® HD Transfection Reagent	1ml	E2311
	5 × 1ml	E2312

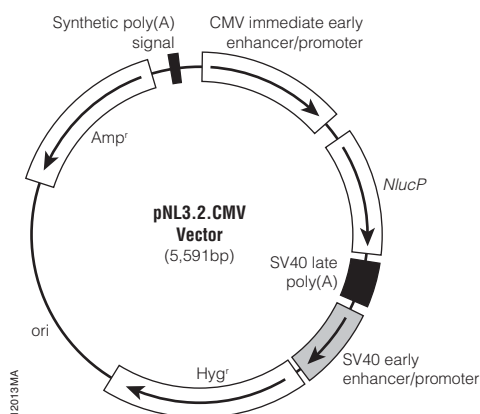


Figure 1. pNL3.2.CMV Vector circle map.

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®-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, researchers 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.

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.

U.S. Pat. No. 8,557,970 and other patents pending.