

Certificate of Analysis

pBiT3.1-N [CMV/HiBiT/Blast] Vector:

Part No.	Size
N236A	20µg



Instructions for use of this product can be found in the *Nano-Glo® HiBiT Lytic Detection System Technical Manual #TM516* and *Nano-Glo® HiBiT Extracellular Detection System Technical Manual #TM523*, available online at: www.promega.com/protocols

Description: The pBiT3.1-N [CMV/HiBiT/Blast] Vector^(a) is configured to append the 11 amino acid HiBiT peptide tag to the amino terminus of the target protein. The vector contains a multiple cloning region to generate an in-frame HiBiT fusion protein. The vector can be used for both stable and transient gene expression and encodes kanamycin resistance for bacterial selection and blasticidin resistance for mammalian selection.

The pBiT3.1-N [CMV/HiBiT/Blast] Vector contains the following features:

- A **CMV immediate-early enhancer/promoter** for constitutive expression in mammalian cells.
- The **HiBiT peptide tag** for bioluminescent detection of the protein of interest.
- A **multiple cloning region** containing unique restriction sites to facilitate gene insertion into the vector.
- A sequence encoding a flexible **linker** between the protein of interest and the HiBiT tag.
- A **kanamycin-resistance gene** for selection of the plasmid in bacteria and a **blasticidin-resistance gene** for selection in mammalian cells.

Concentration: 1µg/µl.

Storage Buffer: The pBiT3.1-N [CMV/HiBiT/Blast] Vector is supplied in 10mM Tris-HCl, 1mM EDTA (pH 7.4).

Storage Conditions: Store at -30°C to -10°C.

Usage Notes:

- Expression of the HiBiT-tagged protein will only result when the proper reading frame is maintained between the HiBiT tag and the gene of interest.
- The flexible linker will be variable in length depending on the restriction enzyme used.
- The insert should also contain a stop codon at the 3' end for termination of the translation.
- Avoid multiple freeze-thaw cycles.

Expiration Date: See product label for 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.

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

Functional Assays

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

Restriction Digestion: The functional purity of the vector DNA is verified by successful digestion with restriction enzymes at the optimal temperature for 1 hour. Samples are examined by agarose gel electrophoresis, comparing cut and uncut vector DNA with marker DNA.

Part# 9PIN236

Printed 8/17



AF9PIN236 0817N236



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.

^(a)Patents Pending.

© 2017 Promega Corporation. All Rights Reserved.

Nano-Glo is a registered trademark of Promega Corporation.

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.

Signed by:

R. Wheeler, Quality Assurance

Part# 9PIN236
Printed in USA 8/17

pBIT3.1-N [CMV/HiBiT/Blast] Vector Features and Circle Map

The following features are present in the pBIT3.1-N [CMV/HiBiT/Blast] Vector based on nucleotide sequence.

CMV promoter	276–866
Chimeric intron	981–1113
T7 RNA polymerase promoter (–17 to +3)	1157–1176
HiBiT	1190–1225
SV40 late polyadenylation signal	1380–1601
EM7 bacterial promoter	1667–1733
Neo-Kan resistance	1747–2541
Col/E1-derived plasmid origin of replication	2696–2732
Synthetic polyadenylation signal sequence	3413–3461 (Reverse)
Blasticidin resistance (Blast ^r) coding region	3485–3883 (Reverse)
SV40 Min Ori	3945–4010 (Reverse)
SV40 Enhancer	4017–4253

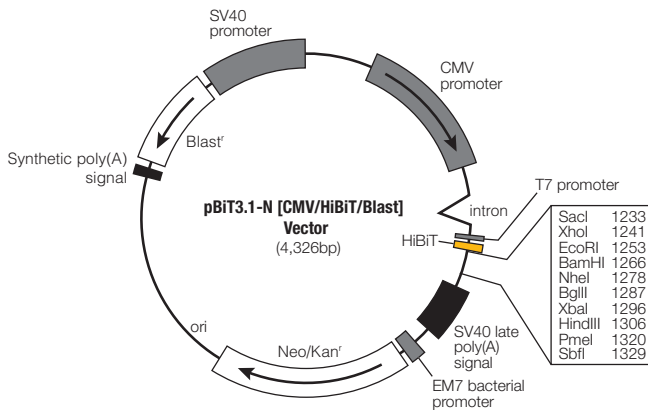


Figure 1. pBIT3.1-N [CMV/HiBiT/Blast] Vector circle map and sequence reference points.



Figure 2. pBIT3.1-N [CMV/HiBiT/Blast] Vector multiple cloning region sequence and unique restriction sites.

Related Products

Product	Size	Cat. #
Nano-Glo [®] HiBiT Lytic Detection System	10ml	N3030
	100ml	N3040
	10 × 100ml	N3050
Nano-Glo [®] HiBiT Extracellular Detection System	10ml	N2420
	100ml	N2421
	10 × 100ml	N2422
Nano-Glo [®] HiBiT Blotting System	100ml	N2410