

Certificate of Analysis

pGL4.31[*luc2P*/GAL4UAS/Hygro] Vector:

Part No. Size (units)
C935A 20µg

Part# 9PIC935

Revised 10/16



Instructions for use of this product can be found in the CheckMate™/Flexi® Vector Mammalian Two-Hybrid System Technical Manual #TM283, available online at: www.promega.com/protocols

Description: The pGL4.31[*luc2P*/Gal4UAS/Hygro] Vector^(a,b,c) is designed to report transcriptional activation using the firefly luciferase reporter gene. This reporter vector contains five consensus binding sequences, or Upstream Activating Sequences, for the Gal4 DNA-binding domain (*Gal4*UAS) and a minimal adenoviral promoter upstream of the firefly luciferase gene. The vector backbone is based on the pGL4 luciferase reporter vector series, which has been engineered to reduce the occurrence of anomalous transcription. The *luc2P* reporter gene is a synthetic firefly luciferase sequence, which has been codon-optimized for high mammalian expression and contains minimal cryptic DNA regulatory elements. In addition, the *luc2P* gene incorporates a protein degradation sequence, PEST, resulting in a destabilized reporter protein with a faster response to transcriptional regulation. The vector also contains an ampicillin selection marker (Amp) for *E. coli* and a hygromycin selection marker (Hygro) for mammalian cells. The pGL4.31[*luc2P*/Gal4UAS/Hygro] Vector is provided for use with the CheckMate™/Flexi® Vector Mammalian Two-Hybrid System (Cat.# C9360).

Concentration: 1µg/µl.

GenBank® Accession Number: DQ487213.

Storage Buffer: The pGL4.31[*luc2P*/Gal4UAS/Hygro] Vector is supplied in 10mM Tris-HCl (pH 7.4), 1mM EDTA.

Storage Conditions: See the Product Information Label for storage recommendations. Avoid multiple freeze-thaw cycles and exposure to frequent temperature changes.

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



AF9PIC935 1016C935



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

Nuclease Assay: Following incubation of 1µg of the pGL4.31[*luc2P*/Gal4UAS/Hygro] Vector in Restriction Enzyme Buffer B at 37°C for 16 hours, no evidence of nuclease activity is detected by agarose gel electrophoresis.

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

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.

©BY USE OF THIS PRODUCT, RESEARCHER AGREES TO BE BOUND BY THE TERMS OF THIS LIMITED USE LABEL LICENSE. 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, and (2) insert and remove nucleic acid sequences in splicing research. No other use or transfer of this product or derivatives is authorized. 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. For any uses outside this label license, contact Promega for supply and licensing information. This product is for research use only; no commercial use is allowed. For a full copy of this label license, including the definition of "commercial use," go to: www.promega.com/LULL.

©U.S. Pat. No. 8,008,006 and European Pat. No. 1341808.

©U.S. Pat. No. 7,728,118.

© 2006–2015, 2016 Promega Corporation. All Rights Reserved.

Flexi is a registered trademark of Promega Corporation. CheckMate is a trademark of Promega Corporation.

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.

Signed by:

R. Wheeler, Quality Assurance

Part# 9PIC935
Printed in USA. Revised 10/16.

pGL4.31[*luc2P*/GAL4UAS/Hygro] Vector circle map and sequence reference points

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

pGL4.31[*luc2P*/GAL4UAS/Hygro] Vector Sequence Reference Points:

GAL4UAS binding sites	31–133
Adenovirus major late promoter	145–185
<i>luc2P</i> firefly luciferase reporter	238–2013
SV40 late polyadenylation signal	2053–2274
SV40 enhancer/early promoter	2322–2740
Synthetic hygromycin resistance (<i>Hygro^r</i>) coding region	2765–3802
Synthetic polyadenylation signal	3826–3874
ColE1-derived plasmid origin of replication	4198–4234
Synthetic β -lactamase (<i>Amp^r</i>) coding region	4989–5849
Synthetic poly(A) signal/transcriptional pause site	5954–6107

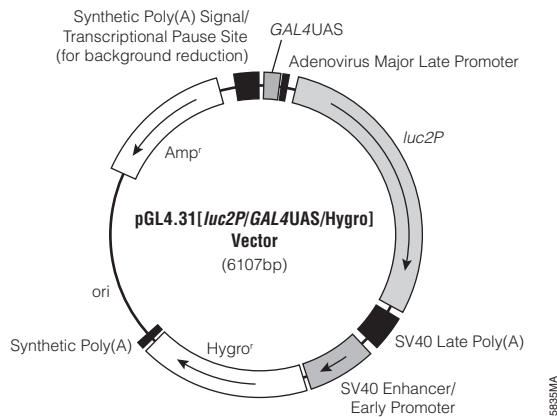


Figure 1. pGL4.31[*luc2P*/GAL4UAS/Hygro] Vector circle map.

Maps of all Promega Vectors are available at:
www.promega.com/vectors