



Deciphering the pGL4 Vector Code

The pGL4 Vectors^(a-e) are the most advanced luciferase reporter gene vectors that maximize expression, minimize off-target responses such as anomalous expression, and respond rapidly to transcriptional dynamics. pGL4 Vectors that express firefly luciferase (Table 1) or *Renilla* luciferase (Table 2) are available. The pGL4 Vectors offer a variety of choices for the three key components shown in Figure 1: The luciferase reporter gene, the upstream elements, including any promoters and response elements, and the selectable marker. The vector nomenclature for the pGL4 Vectors logically reflects these features. As an example, we will decipher the name of the pGL4.31[*luc2P*/CRE/Hygro] Vector.

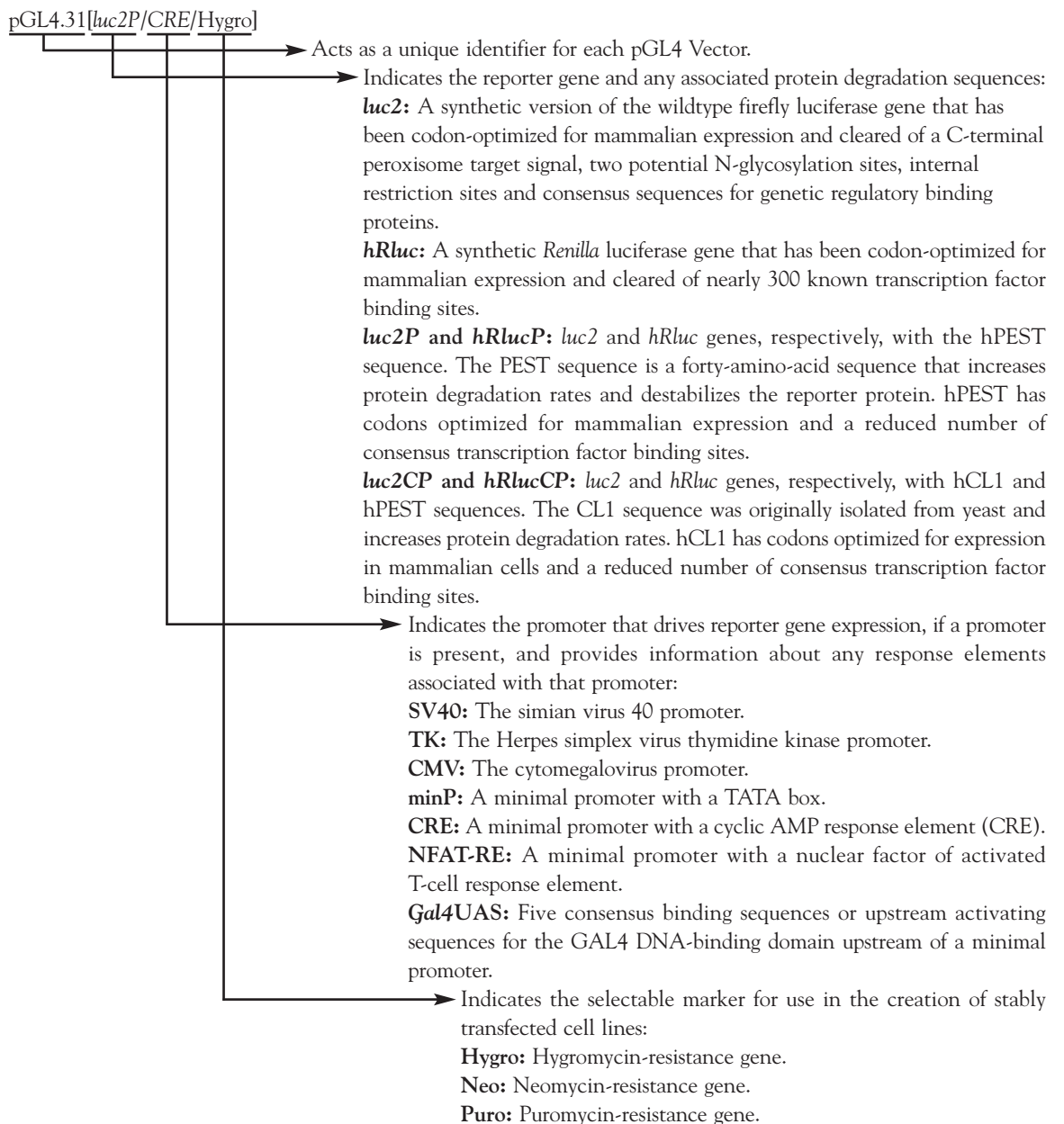


Table 1. pGL4 Vectors With the Synthetic Firefly Luciferase *luc2* gene.

Vector Name	Multiple Cloning Region	Catalog Number
pGL4.10[<i>luc2</i>]	Yes	E6651
pGL4.11[<i>luc2P</i>]	Yes	E6661
pGL4.12[<i>luc2CP</i>]	Yes	E6671
pGL4.13[<i>luc2/SV40</i>]	No	E6681
pGL4.14[<i>luc2/Hygro</i>]	Yes	E6691
pGL4.15[<i>luc2P/Hygro</i>]	Yes	E6701
pGL4.16[<i>luc2CP/Hygro</i>]	Yes	E6711
pGL4.17[<i>luc2/Neo</i>]	Yes	E6721
pGL4.18[<i>luc2P/Neo</i>]	Yes	E6731
pGL4.19[<i>luc2CP/Neo</i>]	Yes	E6741
pGL4.20[<i>luc2/Puro</i>]	Yes	E6751
pGL4.21[<i>luc2P/Puro</i>]	Yes	E6761
pGL4.22[<i>luc2CP/Puro</i>]	Yes	E6771
pGL4.23[<i>luc2/minP</i>]	Yes	E8411
pGL4.24[<i>luc2P/minP</i>]	Yes	E8421
pGL4.25[<i>luc2CP/minP</i>]	Yes	E8431
pGL4.26[<i>luc2/minP/Hygro</i>]	Yes	E8441
pGL4.27[<i>luc2P/minP/Hygro</i>]	Yes	E8451
pGL4.28[<i>luc2CP/minP/Hygro</i>]	Yes	E8461
pGL4.29[<i>luc2P/CRE/Hygro</i>]	No	E8471
pGL4.30[<i>luc2P/NFAT-RE/Hygro</i>]	No	E8481
pGL4.31[<i>luc2P/Gal4UAS/Hygro</i>]	No	C9351

Table 2. pGL4 Vectors with the Synthetic *Renilla* Luciferase *hRluc* gene.

Vector Name	Multiple Cloning Region	Catalog Number
pGL4.70[<i>hRluc</i>]	Yes	E6881
pGL4.71[<i>hRlucP</i>]	Yes	E6891
pGL4.72[<i>hRlucCP</i>]	Yes	E6901
pGL4.73[<i>hRluc/SV40</i>]	No	E6911
pGL4.74[<i>hRluc/TK</i>]	No	E6921
pGL4.75[<i>hRluc/CMV</i>]	No	E6931
pGL4.76[<i>hRluc/Hygro</i>]	Yes	E6941
pGL4.77[<i>hRlucP/Hygro</i>]	Yes	E6951
pGL4.78[<i>hRlucCP/Hygro</i>]	Yes	E6961
pGL4.79[<i>hRluc/Neo</i>]	Yes	E6971
pGL4.80[<i>hRlucP/Neo</i>]	Yes	E6981
pGL4.81[<i>hRlucCP/Neo</i>]	Yes	E6991
pGL4.82[<i>hRluc/Puro</i>]	Yes	E7501
pGL4.83[<i>hRlucP/Puro</i>]	Yes	E7511
pGL4.84[<i>hRlucCP/Puro</i>]	Yes	E7521

(a) For research use only. This product and/or its use is subject to one or more of Promega's U.S. Pat. Nos. 6,387,675 and 6,552,179, Australian Pat. No. 698424 and other pending and issued patents. Researchers must do one of the following: (1) use luminescent assay reagents purchased from Promega for all determinations of luminescence activity resulting from the research use of this product and its derivatives; or (2) contact Promega to obtain a license for the use of the product and its derivatives in conjunction with luminescent assay reagents not purchased from Promega. For all other uses, please contact Promega.

(b) Cat.# C9351, E6651, E6661, E6671, E6681, E6691, E6701, E6711, E6721, E6731, E6741, E6751, E6761, E6771, E8411, E8421, E8431, E8441, E8451, E8461, E8471, E8481: 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.

(c) Cat.# E6881, E6891, E6901, E6911, E6921, E6931, E6941, E6951, E6961, E6971, E6981, E6991, E7501, E7511, E7521: Licensed from University of Georgia Research Foundation, Inc., under U.S. Pat. Nos. 5,292,658, 5,418,155, Canadian Pat. No. 2,105,984 and related patents.

(d) Cat.# E6931: The CMV promoter and its use are covered under U.S. Pat. Nos. 5,168,062 and 5,385,839 owned by the University of Iowa Research Foundation, Iowa City, Iowa, and licensed FOR RESEARCH USE ONLY. Research Use includes contract research for which monetary or other consideration may be received. Other commercial users must obtain a license to these patents directly from the University of Iowa Research Foundation.

(e) Cat.# E8481: The NFAT response element and its use are licensed under one or more of the following patents: U.S. Pat. Nos. 5,837,840, 6,197,925, 5,989,810, 6,096,515, 6,388,052, 6,150,099, 6,171,781, 6,352,830, 6,312,899 and 6,875,571.

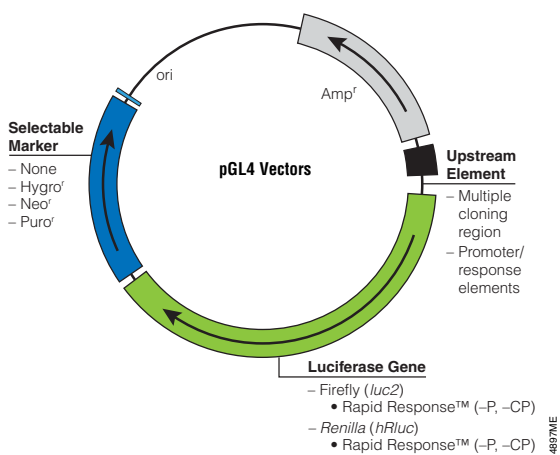


Figure 1. Schematic diagram of the pGL4 Vectors.