

## pMLuc-3 CMV Positive Control

	<b>Cat. No.</b>
pMLuc-3 Positive Control	71219-3

### pMLuc-3 CMV Positive Control sequence landmarks

CMV promoter	1-738
IL-2 signal peptide coding sequence	816-902
Rluc coding sequence	903-1832
Rabbit $\beta$ -globin polyadenylation signal	1842-2124
pUC origin	2288
<i>bla</i> (Ap) coding sequence	2891-3748
Synthetic polyadenylation signal	3822-3977

The pMLuc vectors are designed for cloning and analysis of promoter and enhancer elements. pMLuc-3 CMV Positive Control contains the CMV promoter followed by the IL-2 signal peptide coding sequence and an optimized *Renilla* luciferase (Rluc) cDNA. Secreted Rluc activity is conveniently measured in culture supernatants using the MightyLight™ Rluc Assay Kit (Cat. No. 71220-3). Unique sites are shown on the circle map.

