**GIBCOBRL** 

## Plasmid pCMV•SPORT-βgal

Page 1 of 2

Cat. No. 10586-014 Lot No.  $25 \ \mu g; 0.5 \ \mu g/\mu l$ Store at 4°C

**Description:** Plasmid pCMV•SPORT- $\beta$ gal is used as a positive control for monitoring expression in eukaryotic cells. The plasmid contains the reporter gene  $\beta$ -galactosidase ( $\beta$ -gal) from *E. coli* cloned as a *Not* I fragment into plasmid pCMV•SPORT1. The plasmid also contains a CMV promoter upstream of the  $\beta$ -gal gene, followed by the SV40 t-intron and polyadenylation signal. The  $\beta$ -lactamase gene allows selection for ampicillin resistance in *E. coli*. The sequence and the location of sites for restriction endonucleases, that cleave up to 10 location of sites for restriction endonucleases, that cleave up to 10 times, can be found in the TECH-ONLINE<sup>SM</sup> section of Life Technologies' web page, http://www.lifetech.com.

Storage Buffer: 10 mM Tris-HCl (pH 7.4) 5 mM NaCl 0.1 mM EDTA

<u>**Ouality Control Assay:**</u> This product has passed the following quality control assays: Purity, DNA structure and selected restriction sites are verified by agarose gel analysis; Expression of β-gal activity is verified by transfection into BHK-21 cells using LIPOFECTAMINE<sup>TM</sup> Reagent and *in situ* staining with X-gal.

Doc. Rev. 091197

This product is distributed for laboratory research use only. CAUTION: Not for diagnostic use. The safety and efficacy of this product in diagnostic or other clinical uses has not been established.

For technical questions about this product, call the Life Technologies TECH-LINE® (800) 828-6686]

Page 2 of 2



MAP of pCMV•SPORT-βgal

Cat No. 10586-014