pMYs-GFP Retroviral Vector

CATALOG NUMBER: RTV-052 STORAGE: -20°C

QUANTITY AND CONCENTRATION: 10 μ g at 0.25 μ g/ μ L in TE

Background

Retroviruses are efficient tools for delivering heritable genes into the genome of dividing cells. Most retrovirus vectors including pBABE and pMXs are based on Moloney murine leukemia virus (MMLV). MMLV-based vectors usually are silenced in immature cells including embryonic carcinoma (EC) cells and embryonic stem (ES) cells, and possibly hematopoietic stem cells. Myeloproliferative sarcoma virus (MPSV) and PCC4-cell-passaged myeloproliferative sarcoma virus (PCMV) are mutants of MMLV and can stably express genes in immature cells including ES cells.

Cell Biolabs' pMYs-GFP retroviral vector includes hybrid LTRs containing elements from both MMLV and MPSV/PCMV and is capable of expressing genes in hematopoietic stem cells.

The vector contains the ampicillin-resistance gene, LTRs, package signal and GFP insert (Figure 1).

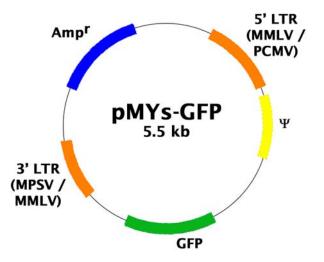


Figure 1. Schematic representation of pMYs-GFP retroviral vector.

Safety Consideration

Remember that you will be working with samples containing infectious virus. Follow the recommended NIH guidelines for all materials containing BSL-2 organisms. Always wear gloves, use filtered tips and work under a biosafety hood.



References

1. Kitamura T., et al., (2003) Exp. Hematol. 31, 1007-1014.

License Information

This product is licensed from the University of Tokyo.

Warranty

These products are warranted to perform as described in their labeling and in Cell Biolabs literature when used in accordance with their instructions. THERE ARE NO WARRANTIES THAT EXTEND BEYOND THIS EXPRESSED WARRANTY AND CELL BIOLABS DISCLAIMS ANY IMPLIED WARRANTY OF MERCHANTABILITY OR WARRANTY OF FITNESS FOR PARTICULAR PURPOSE. CELL BIOLABS's sole obligation and purchaser's exclusive remedy for breach of this warranty shall be, at the option of CELL BIOLABS, to repair or replace the products. In no event shall CELL BIOLABS be liable for any proximate, incidental or consequential damages in connection with the products.

This product is for RESEARCH USE ONLY; not for use in diagnostic procedures.

Contact Information

Cell Biolabs, Inc. 7758 Arjons Drive San Diego, CA 92126 Worldwide: +1 858-271-6500 USA Toll-Free: 1-888-CBL-0505 E-mail: <u>tech@cellbiolabs.com</u> www.cellbiolabs.com

©2008: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.

