

pCMV-VSV-G Envelope Vector

CATALOG NUMBER: RV-110

STORAGE: -20°C

QUANTITY AND CONCENTRATION: 10 µg at 0.25 µg/µL in TE

Background

Retroviruses are efficient tools for delivering heritable genes into the genome of dividing cells. pCMV-VSV-G expresses the G glycoprotein of the vesicular stomatitis virus (VSV-G) under the control of the CMV immediate-early promoter. VSV-G is used in pseudotyping of Moloney Murine Leukemia Virus (MMLV)-based retroviral vectors by mediating viral entry. VSV-G interacts with phospholipid components of the target cell membrane and fosters the fusion of viral and cellular membranes. VSV-G does not require a cell surface receptor and can serve as a surrogate viral envelope protein. The vector contains the ampicillin-resistance gene for propagation and antibiotic selection in bacteria (Figure 1).

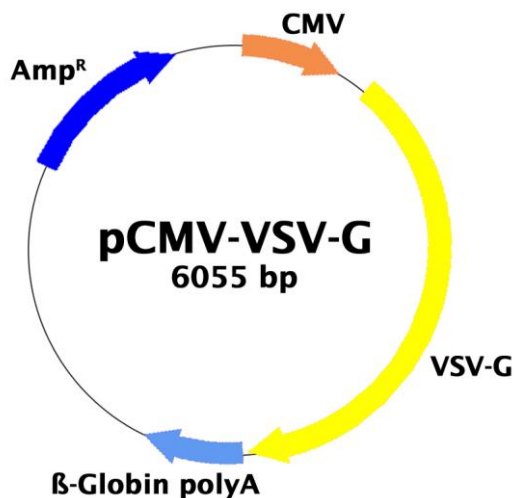


Figure 1. Schematic representation of pCMV-VSV-G envelope vector.

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. Emi, N., *et al.*, (1991) *J. Virol.* **65**, 1202-1207.

Recent Product Citations

1. Jia, R. et al. (2021). The ubiquitin isopeptidase USP10 deubiquitinates LC3B to increase LC3B levels and autophagic activity. *J Biol Chem*. doi: 10.1016/j.jbc.2021.100405.
2. Fu, R.Y. et al (2020). CD4+ T Cells Engineered with FVIII-CAR and Murine Foxp3 Suppress Anti-Factor VIII Immune Responses in Hemophilia A Mice. *Cell Immunol*. doi: 10.1016/j.cellimm.2020.104216.
3. Wąchalska, M. et al. (2020). Palmitoylated mNeonGreen Protein as a Tool for Visualization and Uptake Studies of Extracellular Vesicles. *Membranes (Basel)*. **10**(12):E373. doi: 10.3390/membranes10120373.
4. Simpson, L.M. et al. (2020). Inducible Degradation of Target Proteins through a Tractable Affinity-Directed Protein Missile System. *Cell Chem Biol*. S2451-9456(20)30236-1. doi: 10.1016/j.chembiol.2020.06.013.
5. Tachie-Menson, T. et al. (2020). Characterisation of the biochemical and cellular roles of native and pathogenic amelogenesis imperfecta mutants of FAM83H. *Cell Signal*. doi: 10.1016/j.cellsig.2020.109632.
6. Grabowska, K. et al. (2020). Alphaherpesvirus gB Homologs Are Targeted to Extracellular Vesicles, but They Differentially Affect MHC Class II Molecules. *Viruses*. **12**(4). pii: E429. doi: 10.3390/v12040429.
7. Wu, K.Z.L. et al. (2019). Pathogenic FAM83G palmoplantar keratoderma mutations inhibit the PAWS1:CK1 α association and attenuate Wnt signalling. *Wellcome Open Res*. **4**:133. doi: 10.12688/wellcomeopenres.15403.1.
8. Wenta, T. et al. (2019). HtrA4 Protease Promotes Chemotherapeutic-Dependent Cancer Cell Death. *Cells*. **8**(10):1112. doi: 10.3390/cells8101112.
9. Terada, Y. et al. (2019). Human Pluripotent Stem Cell-Derived Tumor Model Uncovers the Embryonic Stem Cell Signature as a Key Driver in Atypical Teratoid/Rhabdoid Tumor. *Cell Rep*. **26**(10):2608-2621.e6. doi: 10.1016/j.celrep.2019.02.009.
10. Vargas, G. et al. (2019). ERR α promotes breast cancer cell dissemination to bone by increasing RANK expression in primary breast tumors. *Oncogene*. **38**(7):950-964. doi: 10.1038/s41388-018-0579-3.
11. Hafner-Bratkovič, I. et al. (2018). NLRP3 lacking the leucine-rich repeat domain can be fully activated via the canonical inflammasome pathway. *Nat Commun*. **9**(1):5182. doi: 10.1038/s41467-018-07573-4.
12. Macartney, T. J. et al. (2017). An Affinity-directed Protein Missile (AdPROM) System for Targeted Destruction of Endogenous Proteins. *Bio-protocol*. **7**(22): e2614. doi: 10.21769/BioProtoc.2614.
13. Sušjan P, Roškar S, Hafner-Bratkovič I. (2017). The mechanism of NLRP3 inflammasome initiation: Trimerization but not dimerization of the NLRP3 pyrin domain induces robust activation of IL-1 β . *Biochem Biophys Res Commun*. doi: 10.1016/j.bbrc.2017.01.008.
14. Yang, W. S. et al. (2016). Peroxidation of polyunsaturated fatty acids by lipoxygenases drives ferroptosis. *Proc Natl Acad Sci U S A*. doi:10.1073/pnas.1603244113.
15. Gallant, J. N. et al. (2015). EGFR kinase domain duplication (EGFR-KDD) is a novel oncogenic driver in lung cancer that is clinically responsive to afatinib. *Cancer Discov*. **5**:1155-1163.

16. Zhang, T. et al. (2015). Homoharringtonine binds to and increases myosin-9 in myeloid leukemia. *Br J Pharmacol*. doi:10.1111/bph.13359.
17. Amagai, Y. et al. (2015). A point mutation in the extracellular domain of KIT promotes tumorigenesis of mast cells via ligand-independent auto-dimerization. *Sci Rep*. **5**:9775.
18. Abram, M. E. et al. (2015). A cell-based strategy to assess intrinsic inhibition efficiencies of HIV-1 reverse transcriptase inhibitors. *Antimicrob Agents Chemother*. **59**:838-848.
19. Gallaher, Z. R. et al. (2014). Neural proliferation in the dorsal root ganglia of the adult rat following capsaicin-induced neuronal death. *J Comp Neurol*. **522**:3295-3307.
20. Handorf, A. et al. (2014). Endogenously Produced Indian Hedgehog Regulates TGF β -Driven Chondrogenesis of Human Bone Marrow Stromal/Stem Cells. *Stem Cells Dev*. doi:10.1089/scd.2014.0266.
21. Kobayashi, J. et al. (2014). Directed differentiation of patient-specific induced pluripotent stem cells identifies the transcriptional repression and epigenetic modification of NKX2-5, HAND1, and NOTCH1 in hypoplastic left heart syndrome. *PLoS One*. **9**:e102796.
22. Higuchi, A. et al. (2014). Preparation of induced pluripotent stem cells on dishes grafted on oligopeptide under feeder-free conditions. *J Taiwan Inst Chem E*. **45**:295-301.
23. Amagai, Y. et al. (2013). Stem cell factor contributes to tumorigenesis of mast cells via an autocrine/paracrine mechanism. *J. Leukoc. Biol*. **93**:245-250.
24. Okamoto, K. et al. (2012). Dengue virus strain DEN2 16681 utilizes a specific glycochain of syndecan-2 proteoglycan as a receptor. *J. Gen. Virol*. **93**:761-770.
25. Oida, T. et al. (2010). Overexpression of TGF- β ₁ gene induces cell surface localized glucose-regulated protein 78-associated latency-associated peptide/TGF- β . *J. Immunol*. **185**:3529-3535.

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: tech@cellbiolabs.com
www.cellbiolabs.com

©2013-2021: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.