

Description

The PSMA Lentiviruses are replication incompetent, HIV-based, VSV-G pseudotyped lentiviral particles that are ready to transduce nearly all types of mammalian cells, including primary and non-dividing cells. The particles contain a human PSMA (NM_004476.3) driven by a CMV promoter and a puromycin selection marker (Figure 1).

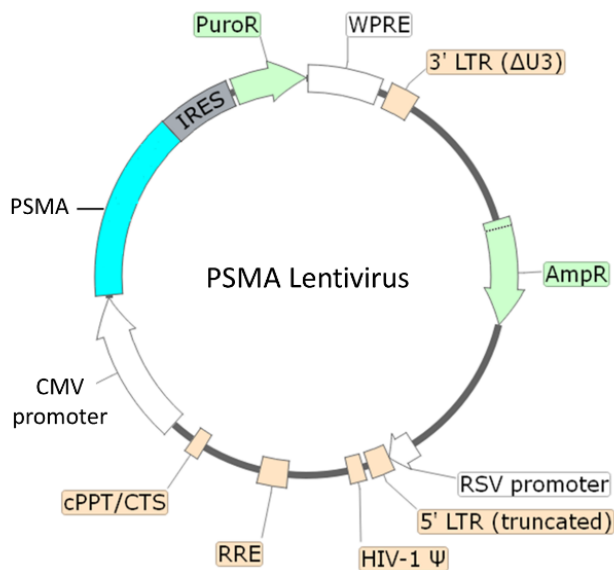


Figure 1: Schematic of the lenti-vector used to generate the PSMA Lentivirus.

Background

PSMA (prostate-specific membrane antigen, also known as Folate hydrolase 1, FOLH1), is highly expressed in prostate cancer cells and is used as a diagnostic and prognostic indicator of prostate cancer. The enzyme has folate hydrolase and peptidase activity. It plays a role in prostate cancer progression through the PI3K-Akt and MAPK-ERK1/2 pathways. PSMA/FOLH1 is used in the clinic as a target for PET (positron emission tomography) imaging of prostate cancer whereas radioactive analogs are used for the treatment of prostate cancer. It is also expressed in other tumor types and in a few normal tissues. PSMA/FOLH1 is the target of CAR-T cells and bi-specific antibodies currently under development.

Application(s)

Generate stable cell line expressing human PSMA with puromycin selection

Formulation

The lentivirus particles were produced from HEK293T cells. They are supplied in cell culture medium containing 90% DMEM + 10% FBS.

Titer

Two vials (500 μ l x 2) of lentivirus at a titer $\geq 10^7$ TU/ml. The titer will vary with each lot; the exact value is provided with each shipment.

Storage



Lentiviruses are shipped with dry ice. For long-term storage, it is recommended to store the lentiviruses at -80°C . Avoid repeated freeze/thaw cycles. Titers can drop significantly with each freeze/thaw cycle.

Biosafety



The lentiviruses are produced with SIN (self-inactivation) lentivector which ensures self-inactivation of the lentiviral construct after transduction and integration into the genomic DNA of the target cells. None of the HIV genes (*gag*, *pol*, *rev*) will be expressed in the transduced cells, as they are expressed from packaging plasmids lacking the packing signal and are not present in the lentivirus particle. Although the pseudotyped lentiviruses are replication-incompetent, they require the use of a Biosafety Level 2 facility. BPS Bioscience recommends following all local federal, state, and institutional regulations and using all appropriate safety precautions.

Notes

To generate a PSMA stable cell line, remove the growth medium 48 hours after transduction and replace it with fresh growth medium containing the appropriate amount of puromycin (as pre-determined from a killing curve) for antibiotics selection of transduced cells. Visit: <https://bpsbioscience.com/cell-line-faq> for guidelines on performing a kill curve.

Figures and Validation Data

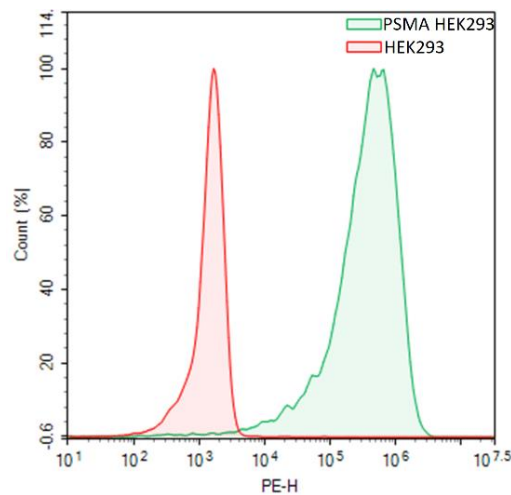


Figure 2. Transduction of HEK293 using PSMA Lentivirus.

Approximately 50,000 HEK293 cells were transduced with 500,000 TU of PSMA lentiviruses. After 48 hours of transduction, the cells were stained using PE anti-PSMA (FOLH1) Antibody (Biolegend #342504) and analyzed by flow cytometry.

Sequence

Human PSMA (NM_004476.3)

```
MWNLLHETDSAVATARRPRWLCAGALVLAGGFFLLGFLFGWFIKSSNEATNITPKHNMKAFLDELKAENIKKFLYNFTQIPHLAG
TEQNFQLAKQIQSQWKEFGLDSELAHYDVLVLLSYPNKTHPNYISIINEDGNEIFNTSLFEP PPPGYENVSDIVPPFSAFSPQGMPEG
DLVYVNYARTEDFFKLERDMKINCSGKIVIARYGKVFVRGNKVKNAQLAGAKGVILYSDPADYFAPGVKSYPDGWNLPGGGVQRG
NILNLNGAGDPLTPGYANEYAYRRGIAEAVGLPSIPVHPIGYDAQKLEKMGGSAPPDSSWRGSLKVPYINVPGFTGNFSTQK
VKMHIHSTNEVTRIYNVIGTLRGAVEPDRYVILGGHRDSWVFGIDPQSGAAVVHEIVRSFGTLKKEGWRPRRTILFASWDAEEF
GLLGSTEWAEENSRLQERGVAYINADSSIEGNYTLRVDCTPLMYSLVHNLTKELKSPDEGFEGKSLYESWTKKSPSEFGMPRIS
KLGSGNDFEVFFQRLGIASGRARYTKNWETNKFSGYPLYHSVYETVELVEKFDPMFKYHLTVAQVRGGMVFELANSIVLPFDCR
DYAVVLRKYADKIYSISMKHPQEMKTYSVSFDLSFAVKNFTEIASKFSERLQDFDKSNPIVLRMMNDQLMFLERAFIDPLGLPDR
PFYRHVIYAPSSHKNKYAGESFPGIYDALFDIESKVDPSKAWGEVKRQIYVA AFTVQAAAETLSEVA
```

Troubleshooting Guide

Visit bpsbioscience.com/lentivirus-faq for detailed troubleshooting instructions. For all further questions, please email support@bpsbioscience.com.

Related Products

<i>Products</i>	<i>Catalog #</i>	<i>Size</i>
Trop2 Lentivirus	78710	500 µl x 2
GPC3 Lentivirus	78711	500 µl x 2
Nectin-4 Lentivirus	78712	500 µl x 2
BCMA Lentivirus	78714	500 µl x 2
FcRL5 Lentivirus	78715	500 µl x 2
GPRC5D Lentiviruses	78716	500 µl x 2
Claudin-9 Lentivirus	78721	500 µl x 2
Claudin-3 Lentivirus	78722	500 µl x 2
Claudin-4 Lentivirus	78723	500 µl x 2
LYPD1 Lentivirus	78724	500 µl x 2