

Product Information

MISSION® TRC3 Human Whole Genome Lentiviral ORF Pool

Catalog Number **ORFPOOLWG**

Storage Temperature **-70 °C**

TECHNICAL BULLETIN

Product Description

An open reading frame (ORF) is a continuous segment of DNA beginning with an initiation codon, methionine ATG, and ending with one of the three termination codons, TAA, TAG, or TGA, that is coded into a polypeptide chain or a protein. An ORF contains the coding sequence of a gene (CDS) and lacks both the 5' and 3' UTRs. These ORFs are inserted into expression vectors (containing an artificial stop codon) and their expression may be modulated (overexpressed) to understand the corresponding gene/protein function. ORFs also play an important role in RNAi and CRISPR rescue experiments, where gene expression is restored (gain-of-function) for gene validation.

Sigma-Aldrich offers the largest collection of fully sequenced and pre-cloned human open reading frames (ORFs) to-date. Our ready-to-use MISSION® TRC3 Human LentiORFs allow for stable integration, enrichment of cells, and long-term gene expression in difficult-to-transfect cell lines utilizing our best-in-class lentiviral manufacturing. This collection provides researchers with the unique tools to gain insights into gene function through modulating gene and protein expression. The Complete Mission TRC3 Human LentiORF collection consists of over 50,000 pre-cloned ORFs, spanning over 14,000 human genes.

Whether looking to perform gene overexpression studies or validate gene knockdown/knockout results from RNAi/CRISPR experiments, LentiORFs are the ideal shortcuts to protein expression and tools for gene analysis. Researchers need to validate "hits" from a screen to eliminate false positives. This means confirming whether the observed phenotype is due the target gene of interest. Re-expressing the gene of interest through ORFs is the best and most accepted way to validate a gene target.

These libraries were introduced in the 2011 *Nature Methods* paper: "A public genome-scale lentiviral expression library of human ORFs." (DOI: 10.1038/nmeth.1638). Supplementary Table 4, containing a summarized view of the contents of the library, is available in two formats ([pdf](#), [xls](#)). For protocols detailing the use of this library, see the Protocols section found below.

The MISSION TRC3 Human Lentiviral ORF Pool consists of plasmid-based, puromycin resistant ORF constructs expressing over 14,000 human genes. On average, there are 1.2 ORFs for each gene in the ORFeome pool. The ORF plasmids are further processed into lentiviral particles to facilitate stable gene expression in both dividing and quiescent cells.

Gene Overexpression Studies

While RNA interference technology allows scientists to "turn off" or silence certain genes, ORFs can be used to "turn on" or "overexpress" particular genes. Flipping the switches on genes one at a time can help reveal the functions of individual genes, such as those that play a role in cancer. The lentiviral expression system has been used in several studies where gene screening, expression analysis, and genetic rescue are important. Overexpression of genes and proteins is widely used in functional genomics, proteomics, and system biology studies.

Some of the studies include:

1. [COT drives resistance to RAF inhibition through MAP kinase pathway reactivation](#)
2. [KRAS and YAP1 converge to regulate EMT and tumor survival](#)

The MISSION TRC3 Human Whole Genome

Lentiviral ORF Pool is a lentiviral pool produced using a proprietary process. Assay for representation of each individual ORF ensures robust library coverage. The pool is provided in ready-to-use lentiviral format at titers of at least 5×10^8 TU/ml via p24 assay. See Table 1 for approximate number of ORF constructs.

Table 1.

Clone and Gene Coverage by Human Whole Genome Pooled ORF Library

| Product | Approximate Total Clone Coverage | Approximate Number of Genes |
|-----------|----------------------------------|-----------------------------|
| ORFPOOLWG | 17,000+ | 14,000+ |

Quality Control

Each MISSION TRC3 Human Whole Genome Lentiviral ORF Pool has been tested for representation of each individual ORF through an internal QC process. The pool has also been tested via p24 assay and has demonstrated a titer of at least 5×10^8 TU/ml.

Components

The MISSION Human Whole Genome Lentiviral ORF Pool includes eight aliquots - enough virus to perform multiple screens. Each aliquot is provided at 5×10^8 TU/ml or higher in a volume of 25 μ l. Please refer to the Certificate of Analysis for the exact titer.

Additionally, each ORF Pool kit will include a USB Flash Drive containing the MISSION Human ORF pool platemap file and the complete list of ORF clone information.

Note: Primers for amplification and sequencing of MISSION ORF are not included, but sequence information is provided. See Table 2 for details.

Included Materials

Positive Target Overexpression Controls and Transduction Efficiency Optimization

- ORF Lenti Human Puro BFP Control Virus High Titer, Catalog Number ORFBFPH

This positive control can be used for measuring transduction efficiency and optimizing ORF delivery in cell lines that do not survive any puromycin exposure. The BFP Control contains a gene encoding BFP driven by the EF1a promoter. This control provides fast visual confirmation of successful transduction.

- ORF Lenti Human Puro TurboGFP Control High Titer, Catalog Number ORFGFPH

This positive control can be used for measuring transduction efficiency and optimizing ORF delivery in cell lines that do not survive any puromycin exposure. The TurboGFP Control contains a gene encoding GFP driven by the EF1a promoter. This control provides fast visual confirmation of successful transduction.

Negative Target Overexpression Controls

- ORF Lenti Human Puro TurboGFP Control High Titer, Catalog Number ORFGFPH
- ORF Lenti Human Puro BFP Control High Titer, Catalog Number ORFBFPH

Either the BFP or GFP control can also be used as a negative control in cell lines that don't already express either fluorophore. The control virus serves as a negative control of lentivirus treatment for the same lentiviral backbone. It allows for validation of the specificity of target expression in target cells, and observation of potential cellular effects due to the transduction process.

Materials And Reagents Required But Not Provided For Transduction of MISSION TRC3 Human Whole Genome Lentiviral ORF Pool into Target Cells

- Hexadimethrine bromide (chemical equivalent of Polybrene), Catalog Number H9268
- Tissue Culture Plates and Related Tissue Culture Supplies
- Puromycin dihydrochloride, Catalog Number P9620

For Purification of Genomic DNA

- GenElute™ Mammalian Genomic DNA Miniprep Kit 70 preps, Catalog Number G1N70

For PCR Amplification of ORF Inserts

- JumpStart™ Taq ReadyMix™, Catalog Number P2893
- Magnesium chloride solution, Catalog Number M8787
- 1% 1× TAE Agarose gel
- DirectLoad™ PCR 100 bp Low Ladder, Catalog Number D3687
- Thermal Cycler

For Purification of Amplified ORF Inserts

- GenElute PCR Clean-Up Kit, Catalog Number NA1020

Precautions And Disclaimers

For R&D use only. Not for drug, household, or other uses. Please consult the Safety Data Sheet for information regarding hazards and safe handling practices.

Though the lentiviral transduction particles produced are replication incompetent, it is recommended they be treated as Risk Group Level 2 (RGL-2) organisms for laboratory handling. Follow all published RGL-2 guidelines for laboratory handling and waste decontamination.

Storage and Stability

All components are stable for at least 6 months after receipt when stored at -70°C . Avoid repeated freeze/thaw cycles, which will severely reduce transduction efficiency.

Procedures

Please refer to the Pooled shRNA Technical Bulletin for more information:

<http://www.sigmaaldrich.com/content/dam/sigma-aldrich/docs/Sigma/Bulletin/shph01bul.pdf>

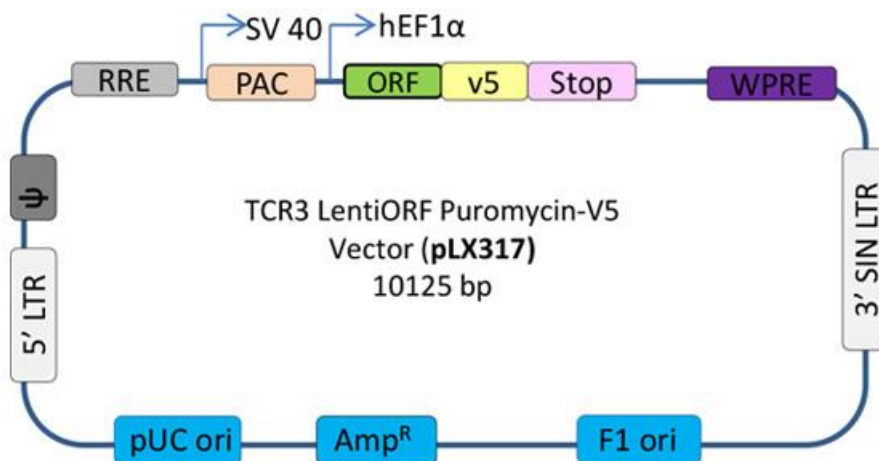
Table 2.

Suggested Materials (**Not Included**) – Primers and Positive Control

| Description | Sequence | Minimum Volume | Approximate Concentration |
|----------------------------------|-----------------------|-------------------|---------------------------|
| Amplification Primer 1 (Forward) | ctgtacaaagtgggtgatc | 750 μl | 20 μM |
| Amplification Primer 2 (Reverse) | caactcctcataaagagacag | 750 μl | 20 μM |
| Sequencing Primer | ttgtacaaagtgggtgatc | 750 μl | 5 μM |

Figure 1.

Mission TRC3 LentiORF Expression Vector Map



The puromycin based LentiORF expression vector systems allow for additional peptide sequences or tags to modify properties of the native protein. Affinity tags can be used to purify the protein further for further structural and characterization studies. Epitope tags (such as V5 found in the blasticidin and puromycin vectors) allow for simple detection or purification via antibodies. The V5 epitope tag is also useful for Western blotting, immunofluorescence, and immunoprecipitation experiments. The addition of reporters and fluorescent tags also allows for gene expression monitoring and cellular localization.

| Vector Label | Function |
|---|---|
| attR1 | Gateway recombination site 1 |
| ORF | Open Reading Frame Insert |
| attR2 | Gateway recombination site 2 |
| pUC ori | High copy replication and maintenance in <i>E. coli</i> |
| Spect | Spectinomycin resistance gene |
| RRE (Rev Response Element) | Increased full-length viral packaging |
| SV40 Promoter | Simian virus 40 (SV40) promoter |
| PAC | Puromycin resistance gene |
| PGK Promoter | Human phosphoglycerate kinase (PGK) promoter |
| Bsd | Blasticidin resistance gene |
| hEF1α Promoter | RNA Polymerase II |
| hCMV Promoter | RNA Polymerase II |
| V5 | V5 Epitope Tag |
| WPRE (Woodchuck Posttranscriptional Regulatory Element) | Enhance transcription of transgenes |
| 5' LTR | 5' long terminal repeat |
| 3' SIN LTR | *Safety* 3' Self inactivating long terminal repeat |
| F1 ori | Phage-derived ORI that allows for the replication and packaging of ssDNA into phage particles |
| AmpR | Ampicillin bacterial selection marker |
| Ψ | Psi viral genome packaging sequence |
| *Barcode (not shown) | Pooled Screening application barcode-Puromycin only |

MISSION® Functional Genomics Licenses

This product is for non-clinical research use only. Use of this product for Commercial Purposes requires a license from Sigma-Aldrich® Corporation or an affiliate thereof. The purchase of this product conveys to the buyer the nontransferable right to use the purchased amount of the product and components of the product in research conducted by the buyer (whether the buyer is an academic or for-profit entity). The buyer cannot sell or otherwise transfer (a) this product (b) its components or (c) materials made using this product or its components to a third party, or otherwise use this product or its components or materials made using this product or its components for Commercial Purposes. "Commercial Purposes" means any activity by a party for consideration, but excludes not-for-profit core facilities providing services within their own research institutions at cost.

One or more of the following licenses may be relevant for MISSION® functional genomic products:

Lentiviral Licenses

This product is licensed under U.S. Pat. Nos. 6,235,522; 6,312,682; 6,541,219; 6,669,936; 6,924,123; 7,056,699; and 7,198,784 and corresponding patents and applications, including foreign equivalents, from Oxford BioMedica (UK) Ltd., Oxford, UK, and is provided for use in academic and commercial *in vitro* and *in vivo* research for elucidating gene function, and for validating potential gene products and pathways for drug discovery and development, but excludes any use of LentiVector® technology for: creating transgenic birds for the purpose of producing useful or valuable proteins in the eggs of such transgenic birds, the delivery of gene therapies, and for commercial production of therapeutic, diagnostic or other commercial products not intended for research use where such products do not consist of or incorporate a lentiviral vector. Information about licenses for commercial uses excluded under this license is available from Oxford BioMedica (UK), Ltd., Medawar Center, Oxford Science Park, Oxford OX4 4GA UK enquiries@oxfordbiomedica.co.uk or BioMedica Inc., 11622 El Camino Real #100, San Diego CA 92130-2049 USA. LentiVector is a registered U.S. and European Community trademark of Oxford BioMedica plc.

This product (based upon the lentikat system) is sub-licensed from Invitrogen Corporation under U.S. Patent Nos. 5,994,136 and 6,165,782 and corresponding patents and applications, including foreign equivalents, for internal research purposes only. Use of this technology for gene therapy applications or bioprocessing other than for nonhuman research use requires a license from Cell Genesys, Inc. Please contact Cell Genesys, Inc. at 342 Lakeside Drive, Foster City, California 94404 USA. Use of this technology to make or sell products or offer services for consideration in the research market requires a license from Invitrogen Corporation, 1600 Faraday Ave., Carlsbad, CA 92008 USA.

Any Sigma-Aldrich® lentiviral backbone-containing products containing a specific genetic component (WPRE) are licensed from the Salk Institute for Biological Studies and covered under the following patents: U.S. Patent Nos. 6,136,597; 6,284,469; 6,312,912; and 6,287,814; and foreign equivalents.

Licensee has a license to sell the Product containing WPRE, under the terms described below. Any use of WPRE outside of Licensee's Product or the Product's intended use, requires a license as detailed below. Before using the Product containing WPRE, please read the following license agreement. If you do not agree to be bound by its terms, contact Licensee within 10 days for authorization to return the unused Product containing WPRE and to receive a full credit. Licensee grants you a non-exclusive license to use the enclosed Product containing WPRE in its entirety for its intended research use. The Product containing WPRE is being transferred to you in furtherance of, and reliance on, such license. Any use of WPRE outside of Licensee's Product or the Product's intended use including for Commercial Purposes, requires a license from the Salk Institute for Biological Studies. Commercial Purposes means any activity by a party for consideration, but excludes not-for-profit core facilities providing services within their own research institutions at cost. This license agreement is effective until terminated. You may terminate it at any time by destroying all Products containing WPRE in your control. It will also terminate automatically if you fail to comply with the terms and conditions of the license agreement. You shall, upon termination of the license agreement, destroy all Products containing WPRE in your control, and so notify Licensee in writing. This License shall be governed in its interpretation and enforcement by the laws of the State of California.

Contact for WPRE Licensing:

The Salk Institute for Biological Studies
10010 North Torrey Pines Road
La Jolla, CA 92037
Attn.: Office of Technology Management
Phone: (858) 453-4100 extension 1703
Fax: (858) 546-8093

This product and its use are the subject of one or more of the following patents 6,682,907; 7,981,671; 8,093,042; 8,367,068; 8,450,087; 8,460,678; 8,512,993; 8,512,994; and 9,238,824; and/or other pending foreign patent applications licensed to Sigma-Aldrich. The purchase of this product conveys to the buyer the non-transferable right to use the purchased amount of the product and components of the product in research conducted by the buyer (whether the buyer is an academic or for-profit entity). The buyer cannot disclose information, sell or otherwise transfer this product, its components or materials made using this product or its components to a third party or otherwise use this product or its components or materials made using this product or its components for any commercial purposes. If the buyer is not willing to accept the limitations of this limited use statement, Sigma-Aldrich is willing to accept return of the product with a full refund. For information on purchasing a license to the DNA-Flap technology for purposes other than research, contact the Transfer of Technology Office, Institut Pasteur, 28 rue du Docteur Roux, 75 724 Paris Cedex 15 (www.pasteur.fr).

DirectLoad, GenElute, JumpStart, and ReadyMix are trademarks and MISSION is a registered trademark of Sigma-Aldrich Co. LLC.

JC,RR,PHC,MAM 12/19-1