

Dharmacon™ Trans-Lentiviral ORF Packaging Kit with Calcium Phosphate Transfection Reagent

Cat. #TLP5916, TLP5918, TLP5919, TLP5920

Product Description:

The Trans-Lentiviral Packaging System efficiently generates replication-incompetent, HIV-1-based lentivirus used to deliver and express your ORF clone of interest in either dividing or non-dividing mammalian cells. The multi-plasmid Trans-Lentiviral Packaging Mix is designed to express viral proteins at ratios optimal for lentiviral assembly and production. In combination with your choice of transfer vector, this packaging system produces active but replication-incompetent viral particles.

Quality Control:

All DNA is quantified by UV Spectrophotometry prior to shipment. Our Calcium Phosphate Transfection Reagent components are functionally tested using the Trans-Lentiviral Packaging System with HEK293T Packaging Cell Line and must meet minimum viral titer specifications. The HEK293T packaging cells are tested for viability and selected for high expressing virus titers; they are also tested for Mycoplasma contamination.

Shipping and Storage:

The Dharmacon Trans-Lentiviral ORF Packaging Kit includes the following components as listed in Table 1. The amount of Trans-Lentiviral Packaging Mix and Calcium Phosphate transfection reagent supplied is sufficient to perform 10-100 packaging events in 100 mm plates, depending on the kit size purchased.

Table 1. Components of the Trans-Lentiviral ORF Packaging Kit with Calcium Phosphate Transfection Reagent:

Kit Contents	TLP5916 (10 rxn)	TLP5918 (10 rxn + cells)	TLP5919 (50 rxn)	TLP5920 (100 rxn)	Shipping Conditions	Storage
Trans-Lentiviral Packaging Mix	285 µg	285 µg	285 µg × 5	285 µg × 10	Wet Ice	-20° C
HEK293T Packaging Cell Line*	not included	1.0 mL	not included	not included	Dry Ice	Liquid N ₂
CaCl ₂ Reagent	1.2 mL	1.2 mL	6 mL	6 mL × 2	Wet Ice	-20° C
2x HBSS Reagent	12 mL	12 mL	60 mL	60 mL × 2	Wet Ice	-20° C
Precision LentiORF RFP Control DNA	45 µg (0.45 µg/µL)	45 µg (0.45 µg/µL)	not included	not included	Wet Ice	-20° C

*Items can be ordered individually.

To allow any CO₂ that may have dissolved into the medium from the dry ice in shipping to dissipate, please store HEK293T cells at -80 °C for at least 48 hours before thawing. Warning: always wear protective eyewear when handling HEK293T cell cryovials stored in liquid nitrogen.

Important Safety Note:

Please follow the safety guidelines for use and production of vector-based lentivirus as set by your institution's biosafety committee. In general, the NIH Office of Biotechnology BSL2 or BSL2+ guidelines should be followed when handling any lentiviral particles. For guidance on containment for lentiviral vectors, please refer to the Recombinant DNA Advisory Committee (RAC) guidelines for research with lentiviral vectors: [Click here](#).

GE, imagination at work and GE monogram are trademarks of General Electric Company. Dharmacon is a trademark of GE Healthcare companies. All other trademarks are the property of General Electric Company or one of its subsidiaries. ©2017 General Electric Company—All rights reserved. First version published June 2014. GE Healthcare UK Limited, Amersham Place, Little Chalfont, Buckinghamshire, HP7 9NA, UK



Orders can be placed at:
gelifsciences.com/dharmacon

Customer Support: cs.dharmacon@ge.com
Technical Support: ts.dharmacon@ge.com or
1.800.235.9880; 303.604.9499 if you have any questions.