

DharmaconTM Trans-Lentiviral ORF Packaging kit with calcium phosphate transfection reagent

Cat. #TLP5916, TLP5918, TLP5919, TLP5924

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 cells and must meet minimum viral titer specifications. The HEK293T cells are tested for viability and selected for high expressing virus titers; they are also tested for Mycoplasma contamination.

Shipping and storage

The Trans-Lentiviral shRNA 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 shRNA 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](#).

If you have any questions, contact

t +44 (0) 1223 976 000 (UK) **or** +1 800 235 9880 (USA); +1 303 604 9499 (USA)

f + 44 (0)1223 655 581

w horizondiscovery.com/contact-us **or** dharmacon.horizondiscovery.com/service-and-support

Horizon Discovery, 8100 Cambridge Research Park, Waterbeach, Cambridge, CB25 9TL, United Kingdom

All trademarks are the property of Horizon Discovery Company unless otherwise specified. ©2018 Horizon Discovery Group Company—All rights reserved. First published November 2016. UK Registered Head Office: Building 8100, Cambridge Research Park, Cambridge, CB25 9TL, United Kingdom.