

# Human cancer subset TRC lentiviral gene family —glycerol

Cat. #RHS5143

**Vector** pLKO.1  
**Species** *Homo sapiens*  
**Source** Custom Manual Curation

## Product description

The Expression Arrest™ TRC Library is the result of a collaborative research effort based at the Broad Institute of MIT and Harvard, and includes six MIT and Harvard associated research institutions and five international life sciences organizations. The goal of TRC is to create lentiviral shRNA libraries targeting 15,000 human and 15,000 mouse annotated genes with multiple constructs per gene. We have partnered with the TRC to make these shRNA libraries available to researchers worldwide.

## Custom manual curation

The Cancer Gene Sets are a compilation of many published cancer related gene lists centered around the National Cancer Institutes Center for Bioinformatic's (NCICB) Cancer Gene Data Curation Project. These published and curated lists include findings from multiple bioinformatic studies.

## Quality control

All cultures are checked for growth prior to shipment.

## Shipping and storage

Collections are in 96-well plate format as bacterial cultures of *E. coli* (DH5α) in 2x LB broth (low salt)\* medium with 8% glycerol, 100 µg/mL carbenicillin. Collections are shipped on dry ice and should be stored at -80 °C. \*1x LB medium can be used instead of 2x LB broth medium.

To allow any CO<sub>2</sub> that may have dissolved into the media from the dry ice in shipping to dissipate, please store constructs at -80 °C for at least 48 hours before thawing.

## 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.

## 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](http://horizondiscovery.com/contact-us) **or** [dharmacon.horizondiscovery.com/service-and-support](http://dharmacon.horizondiscovery.com/service-and-support)

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