

## Certificate Of Analysis

### Product Description

KILR® Retroparticles for Suspension Cells (G418) contains MMLV particles, that can transduce a wide variety of suspension cells with a housekeeping gene tagged with ProLabel® (ePL), a β-gal reporter fragment. This results in a high level of expression of the fusion protein inside the target cells. Target cell death in a cytotoxicity assay results in the release of the ePL-tagged protein into the medium. Addition of KILR detection reagent, containing the complementing β-gal reporter fragment, EA, results in complementation of the two enzyme fragments (EA and ePL). The resulting functional enzyme hydrolyzes a substrate to generate a chemiluminescent signal.

| Product Information |  |
|---------------------|--|
| Product Name        | KILR® Retroparticles for Suspension Cells (G418) |
| Cryovial Label      | KILR Retroparticles (G418) - Suspension          |
| Lot #               | 23C1408  |
| Vial Contents       | 1.0 ml   |

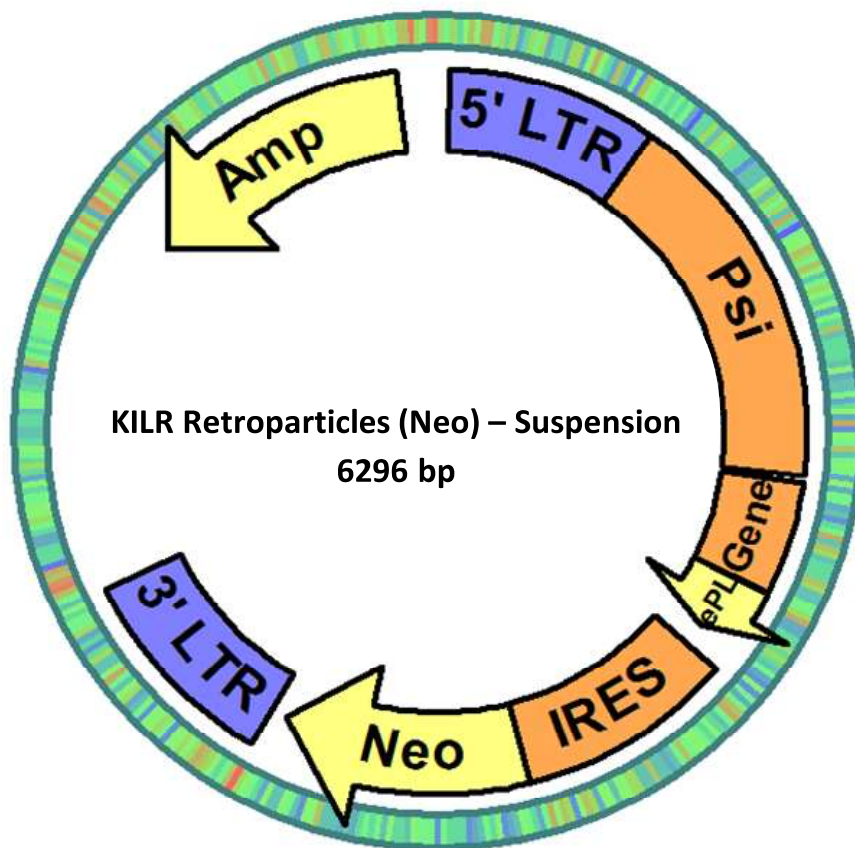
| Shipping and Storage Information |                                     |
|----------------------------------|-------------------------------------|
| Storage Conditions               | Store at -80°C. Do not freeze/thaw. |
| Shipping Conditions              | Dry ice (-80°C)                     |
| Expiration Date                  | 6/15/2025                           |

**IMPORTANT SAFETY NOTE:** Replication-defective retroviral particles, such as provided in this product are not known to cause any diseases in humans or animals. However, retroviral particles can transduce, express protein and/or integrate into human cells. Accordingly, this material is in Risk Group 2 and should be handled under BSL2 controls as defined by the US Public Health Service. Please refer to the CDC Biosafety Manual: <http://www.cdc.gov/biosafety/publications/bmbl5/index.htm> for details.

| Use and Handling    |  |
|---------------------|--|
| Biosafety Level     | 2 (Biosafety classification is based on US Public Health Service guidelines)                 |
| Product User Manual | KILR™ Retroparticles For Cytotoxicity Assays   |
| Single Use          | For one time use only. Repeated freeze/ thaw will result in loss of activity.                |
| Recommended Use     | Transduction of suspension cells to generate KILR cell lines, for use in cytotoxicity assays |
| Acceptable Use      | Research Use Only. Not for use in Humans.  |

## Certificate Of Analysis

| Retroparticles Vector Information |   |  |
|-----------------------------------|---|--|
| Vector                            | pMLV backbone   | Vector identification was confirmed by sequencing  |
| Viral Elements                    | 5' and 3' LTRs  |  |
| Viral Replication Status          | Replication incompetent retroviral particles - helper virus free.<br>Retrovirus can only infect dividing cells. |  |
| Antibiotic Resistance             | Neomycin  | Expression driven by 5' retroviral LTR promoter  |
| KILR Reporter                     | Housekeeping Gene tagged with ePL   | Hybrid murine moloney leukemia virus/murine sarcoma virus (MMLV/MSV) retroviral LTR promoter |
| Viral Pseudotype                  | VSV-G envelope  | Suitable for infecting all mammalian cell types  |



## Certificate Of Analysis

### Quality Control Data

Titer: The titer of viral particles was determined by plaque formation at limiting dilution on adherent cells after 7 days under selection with the appropriate antibiotic concentration. Titer (plaque forming units/mL, abbreviated pfu/mL) was calculated by multiplying the number of colonies per well by the dilution factor divided by the volume (in mL) used in the experiment. Additional details of QC tests available upon request.

| Analytical QC Tests |                              |
|---------------------|------------------------------|
| Viral Titer         | 7.5 X 10 <sup>7</sup> cfu/ml |
| Mycoplasma          | Passed                       |
| Sterility           | Passed                       |

Functional: KILR™ ePL-reporter protein expression was functionally assessed in indicated cell line(s) to confirm transduction of the KILR Retroparticles. Following transduction and antibiotic selection for 7 days the target cells were treated as described for "Total Lysis Control" wells by addition of lysis buffer and KILR Detection Reagents (DiscoverX, Cat. # 97-0001).

| Functional Test |                   |                   |           |                   |
|-----------------|-------------------|-------------------|-----------|-------------------|
| Cell Line       | Average RLU (-EA) | Average RLU (+EA) | S:B Ratio | Days in Selection |
| Jurkat          | 31220             | 1294730           | 41        | 7                 |

### Signatures

Signature: \_\_\_\_\_

Documented by Rene Hoffman  
Scientist I

Date: 4/7/2023

Signature:  \_\_\_\_\_

Approved by Chao-Tsung Yang, PhD  
Principal Scientist

Date: 04/06/2023

Certificate Of Analysis