

## KILR<sup>®</sup> MM-1R Cell Pool

**Catalog Number:** 97-1045P052

**Lot Number:**

See Vial

**Contents:** 1 x 10<sup>6</sup> cells per vial in 1 mL

### Background

KILR cell lines are engineered to express an enhanced Prolabel (ePL) tagged housekeeping gene and may sometimes overexpress an untagged version of a receptor. Once the cells have been lysed the ePL-tagged protein is released into the media. Addition of enzyme acceptor (EA) will cause the complementation of the β-galactosidase enzyme fragments, EA and ePL. The resulting functional enzyme will hydrolyze its substrate to generate a chemiluminescent signal.

### Product Information

**Cell Background:** MM-1R

**Cell Line Species:** Human

**Cell Line Source:** ATCC

**Cell Type:** B lymphoblast

**Culture Mode:** mixed, suspension and lightly attached cells

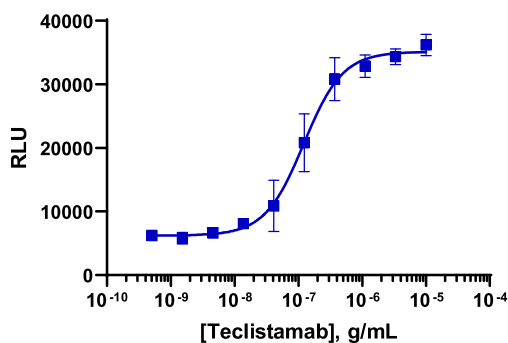
**ADCC Validation:** Teclistamab (BCMA x CD3 mAb)

**CDC Validation:** N/A

**Storage:** Short term (<24 h): Store at -80°C; Long term (>24 h): Store in vapor phase of liquid nitrogen.

### T Cell Redirection Assay Performance

Cells were plated in a 384-well plate and incubated at 37°C and 5% CO<sub>2</sub> for the indicated amount of time. Antibody was added and opsonized for the indicated time (below). Effector cells were added and the plate was incubated at 37°C/5% CO<sub>2</sub> using the assay conditions described below. Target cell death was detected using the KILR detection reagent according to the recommended protocol. Additional reagents needed are noted on this document.



<b>Target Cell Number/Well:</b>	2,500
<b>Effector Type:</b>	pan T cells
<b>Effector Species:</b>	Human
<b>Effector Cell Number/Well:</b>	25,000
<b>Effector to Target Ratio:</b>	10:1
<b>Control Antibody:</b>	Teclistamab
<b>Cell Seeding Time (minutes):</b>	30
<b>Antibody Incubation Time (minutes):</b>	30
<b>Antibody Incubation Temperature (°C):</b>	37
<b>Assay Incubation Time (minutes):</b>	2,880
<b>Assay Incubation Temperature (°C):</b>	37
<b>Incubation with KILR detection reagent (hours):</b>	1
<b>EC<sub>50</sub> for Antibody:</b>	121.5
<b>Signal:Background Ratio:</b>	5.7
<b>Max % Lysis:</b>	N/A

MM.1R cells are semi-adherent. For passaging, transfer all culture volume to a fresh tube (10 mL or 50 mL conical tube) then rinse the flask with PBS and lift adherent population using 0.05% - 0.1% Trypsin-EDTA solution. Use the cell suspension which is set aside to harvest detached cells. Dilute in fresh medium to density of  $> 0.5 \times 10^6$  cells/mL. These cells will stop proliferating at a low density ( $< 0.5 \times 10^6$  cells/mL).

### Passage Stability

This cell line has been confirmed to be stable through 15 passages with no significant drop in assay window or change in EC<sub>50</sub>.

### Mycoplasma Testing

This lot was tested and found to be free of mycoplasma contamination. Data available upon request.

### Required Materials

The following additional materials are required but not provided:

Product Use*	Product Description	Catalog Number
Detection	KILR <sup>®</sup> Detection Kit	97-0001M
Cell Culture	AssayComplete™ Cell Culture Kit-101	92-3101G
Cell Plating	AssayComplete™ Cell Plating 0 Reagent	93-0563R0A
Cell Detachment	AssayComplete™ Cell Detachment Reagent	92-0009
Cell Thawing	AssayComplete™ Thawing Reagent T6	92-4106TR
Cell Freezing	AssayComplete™ Freezing Reagent F5	92-5105FR
Ligand Dilution	AssayComplete™ Protein Dilution Buffer	92-0023

\*Please inquire about our cell line-specific AssayComplete Starter Packs to get you started with your cell culture needs.

### Required Antibiotics

Antibiotic Name	Concentration (µg/mL)	Catalog Number
AssayComplete™ Puromycin	Not Applicable	Not Applicable
AssayComplete™ Hygromycin B	Not Applicable	Not Applicable
AssayComplete™ G418	250	92-0030

### Additional Ligand Information

**Control Compound:** Teclistamab (BCMA x CD3 mAb)

**Vendor:** IchorBio (Catalog No. ICH5209) or equivalent

### Limited Use License Agreement

These products may be covered by issued US and/or foreign patents, patent application and subject to Limited Use Label License.

Please visit <https://www.discoverx.com/legal-policies/> for limited-use label license agreement and trademark information for this product.

© 2024 Eurofins DiscoverX, a Eurofins Discovery Company. All Rights Reserved.