

# KILR<sup>®</sup> U118-MG Cell Pool

**Catalog Number:** 97-1018P032

**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:** U118-MG

**Cell Line Species:** Human

**Cell Line Source:** ATCC

**Cell Type:** Glioblastoma

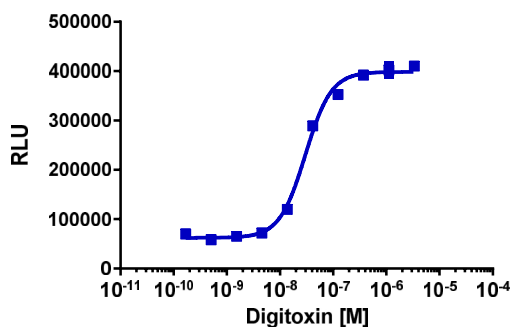
**Culture Mode:** Adherent

**Cytotoxicity Validation:** N/A

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

## Cytotoxicity Assay Performance

Cells were plated in a 96-well plate and incubated at 37°C and 5% CO<sub>2</sub> for indicated amount of time. Compound was 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>	20,000
<b>Control Compound:</b>	Digitoxin
<b>Cell Seeding Time (minutes):</b>	30
<b>Compound Incubation Time (minutes):</b>	960
<b>Compound Incubation Temperature (°C):</b>	37
<b>EC<sub>50</sub> for Compound:</b>	30.1
<b>Signal:Background Ratio:</b>	7.1
<b>Max % Lysis:</b>	78

**Recommended Culture Conditions:** Split ratio: 1:3 to 1:8; Medium renewal: 2-3 times per week.

### 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-105	92-3105G
Cell Plating	AssayComplete™ Cell Plating 39 Reagent	93-0563R39A
Cell Detachment	AssayComplete™ Cell Detachment Reagent	92-0009
Cell Thawing	AssayComplete™ Thawing Reagent T1	92-4101TR
Cell Freezing	AssayComplete™ Freezing Reagent F1	92-5101FR
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	500	92-0030

### Additional Ligand Information

**Control Compound:** Digitoxin

### 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.