

## KILR<sup>®</sup> THP-1 Cell Pool

**Catalog Number:** 97-1011P014

**Lot Number:**

See Vial

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

### Background

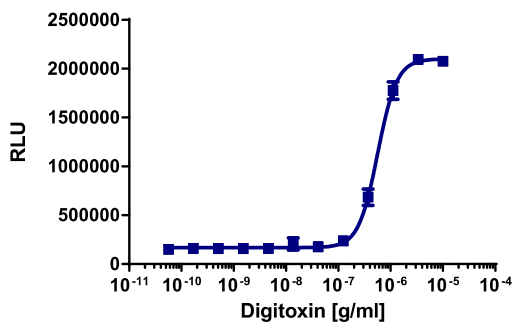
KILR cell lines are engineered to express an enhanced Prolabel (ePL) tagged housekeeping gene and may sometimes overexpress an untagged version of a receptor. Once the cells have been lysed the ePL-tagged protein is released into the media. Addition of enzyme acceptor (EA) will cause the complementation 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:** THP-1  
**Cell Line Species:** human  
**Cell Line Source:** ATCC  
**Cell Type:** Monocyte (AML)  
**Culture Mode:** Suspension  
**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>	10,000
<b>Control Compound:</b>	Digitoxin
<b>Cell Seeding Time (minutes):</b>	30
<b>Compound Incubation Time (minutes):</b>	360
<b>Compound Incubation Temperature (°C):</b>	37
<b>EC<sub>50</sub> for Compound:</b>	595
<b>Signal:Background Ratio:</b>	14.1
<b>Max % Lysis:</b>	60%

**Recommended Culture Conditions:** Start cultures at 2-4 x 10<sup>5</sup> viable cells/mL in complete medium containing 0.05mM β-mercaptoethanol. Subculture when cell concentration reaches 8x10<sup>5</sup> cells/mL. Do not allow the cell concentration to exceed 1 x 10<sup>6</sup> cells/mL. Add fresh medium every 2-3 days.

### 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 39 Reagent	93-0563R39A
Cell Detachment	Not Applicable	Not Applicable
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	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.