

## KILR<sup>®</sup> WIL2-S Cell Pool

**Catalog Number:** 97-1013P027

**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  $\beta$ -galactosidase enzyme fragments, EA and ePL. The resulting functional enzyme will hydrolyze its substrate to generate a chemiluminescent signal.

### Product Information

**Cell Background:** WIL2-S

**Cell Line Species:** Human

**Cell Line Source:** ATCC

**Cell Type:** B lymphoblast (hereditary spherocytosis)

**Culture Mode:** Suspension

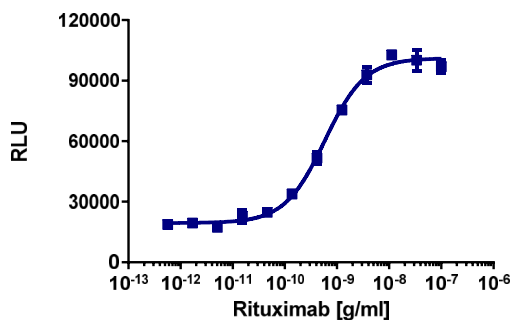
**ADCC Validation:** Rituximab

**CDC Validation:** N/A

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

### ADCC Assay Performance

Cells were plated in a 96-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>	10,000
<b>Effector Type:</b>	primary PBMCs (frozen)
<b>Effector Species:</b>	Human
<b>Effector Cell Number/Well:</b>	250,000
<b>Effector to Target Ratio:</b>	25:1
<b>Control Antibody:</b>	Rituximab
<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>	180
<b>Assay Incubation Temperature (°C):</b>	37
<b>Incubation with KILR detection reagent (hours):</b>	1
<b>EC<sub>50</sub> for Antibody:</b>	0.58
<b>Signal:Background Ratio:</b>	6
<b>Max % Lysis:</b>	20

**Recommended Culture Conditions:** Start cultures at  $2 \times 10^5$  cells/mL; maintain between  $1 \times 10^5$  and  $1 \times 10^6$  cells/mL. Add fresh medium every 2 to 3 days, Cells perform best if cultured without antibiotics for 48h prior to use in ADCC assay.

### Passage Stability

This cell line has been confirmed to be stable through 15 passages with no significant drop in assay window or change in  $EC_{50}$ .

### 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:** Rituximab

**Vendor:** InvivoGen (Catalog No. hcd20-mab13)

### 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.