

## KILR<sup>®</sup> SKOV3 Cell Pool

**Catalog Number:** 97-1007P023

**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:** SKOV3

**Cell Line Species:** Human

**Cell Line Source:** ATCC

**Cell Type:** Ovarian adenocarcinoma

**Culture Mode:** Adherent

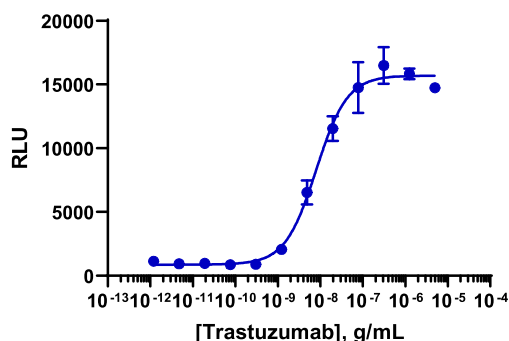
**ADCC Validation:** Trastuzumab

**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>	5,000
<b>Effector Type:</b>	KILR CD16 effectors
<b>Effector Species:</b>	Human
<b>Effector Cell Number/Well:</b>	50,000
<b>Effector to Target Ratio:</b>	10:1
<b>Control Antibody:</b>	Trastuzumab
<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>	240
<b>Assay Incubation Temperature (°C):</b>	37
<b>Incubation with KILR detection reagent (hours):</b>	16
<b>EC<sub>50</sub> for Antibody:</b>	8
<b>Signal:Background Ratio:</b>	17
<b>Max % Lysis:</b>	27

**Recommended Culture Conditions:** Split ratio: 1:2 to 1:6; Medium renewal: 2-3 times per week.

**Note:** these cells are relatively slow-growing; typically split once 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-102	92-3102G
Cell Plating	AssayComplete™ Cell Plating 39 Reagent	93-0563R39A
Cell Detachment	AssayComplete™ Cell Detachment Reagent	92-0009
Cell Thawing	AssayComplete™ Thawing Reagent T8	92-4108TR
Cell Freezing	AssayComplete™ Freezing Reagent F6	92-5106FR
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:** Trastuzumab

**Vendor:** Eurofins DiscoverX<sup>®</sup> (Catalog No. 92-1359)

### 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.