

KILR[®] SKBR3 Cell Pool

Catalog Number: 97-1002P018

Lot Number:

See Vial

Contents: 1 x 10⁶ 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: SKBR3

Cell Line Species: Human

Cell Line Source: ATCC

Cell Type: Breast adenocarcinoma

Culture Mode: Adherent

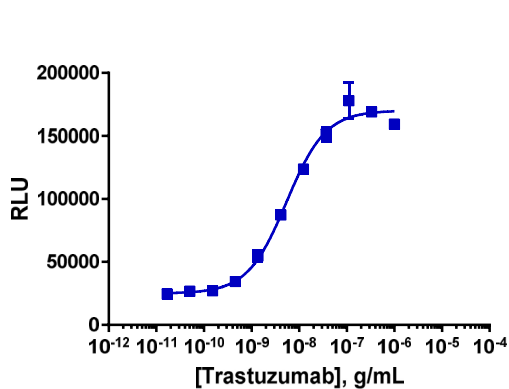
ADCC Validation: Trastuzumab

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% CO₂ 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% CO₂ 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.



Target Cell Number/Well:	5,000
Effector Type:	primary PBMC's (frozen)
Effector Species:	Human
Effector Cell Number/Well:	125,000
Effector to Target Ratio:	25:1
Control Antibody:	Trastuzumab
Cell Seeding Time (minutes):	30
Antibody Incubation Time (minutes):	30
Antibody Incubation Temperature (°C):	37
Assay Incubation Time (minutes):	180
Assay Incubation Temperature (°C):	37
Incubation with KILR detection reagent (hours):	1
EC₅₀ for Antibody:	5.5
Signal:Background Ratio:	5.7
Max % Lysis:	20

Recommended Culture Conditions: Split ratio: 1:2; Medium renewal: 2-3 times per week. Seeding density: 3.0×10^4 viable cells/cm². Note: These cells recover slowly from cryopreservation and tend to pile on each other. Do not trypsinize cells until they have reached 70-80% confluence. In the first week of growth, it is important to retain the floating cells by gentle centrifugation and adding the cells back to the same flask containing the attached cells when replacing the medium.

Passage Stability

This cell line has been confirmed to be stable through 30 passages with no significant drop in assay window or change in EC₅₀.

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 [®] 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	250	92-0030

Additional Ligand Information

Control Compound: Trastuzumab

Vendor: Eurofins DiscoverX[®] (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.