

KILR[®] NCI-H292 Cell Pool

Catalog Number: 97-1008P024

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: NCI-H292

Cell Line Species: Human

Cell Line Source: ATCC

Cell Type: NSCLC (Squamous)

Culture Mode: Adherent

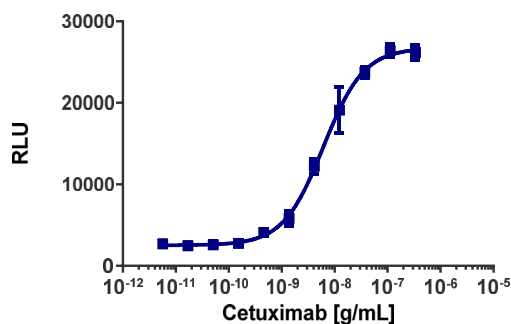
ADCC Validation: Cetuximab

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 PBMCs (frozen) |
| Effector Species: | Human |
| Effector Cell Number/Well: | 200,000 |
| Effector to Target Ratio: | 40:1 |
| Control Antibody: | Cetuximab |
| 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: | 6.1 |
| Signal:Background Ratio: | 9.5 |
| Max % Lysis: | 20 |

Recommended Culture Conditions: Split ratio: 1:3 to 1:8; Medium renewal: every 2-3 days.

Note: cells in confluent monolayer may require additional time to detach with trypsin (e.g. up to 10' at 37°C is normal).

Passage Stability

This cell line has been confirmed to be stable through 15 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-101 | 92-3101G |
| Cell Plating | AssayComplete™ Cell Plating 39 Reagent | 93-0563R39A |
| Cell Detachment | AssayComplete™ Cell Detachment Reagent | 92-0009 |
| 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 | 300 | 92-0030 |

Additional Ligand Information

Control Compound: Cetuximab

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.