

Certificate Of Analysis

Background

KILR Bioassay cells are engineered to express an enhanced Prolabel (ePL) tagged housekeeping gene; some cell models may also overexpress an untagged version of an antigen of interest. Incubation of the KILR Bioassay cells with an antibody that recognizes the antigen of interest, and the desired effector cells results in immune cell-mediated death of the KILR Bioassay cells, releasing the ePL -tagged protein into the media. Addition of exogenous enzyme acceptor (EA) to the medium results in complementation of the β -galactosidase enzyme fragments, EA and ePL. The resulting functional enzyme will hydrolyze its substrate to generate a chemiluminescent signal .

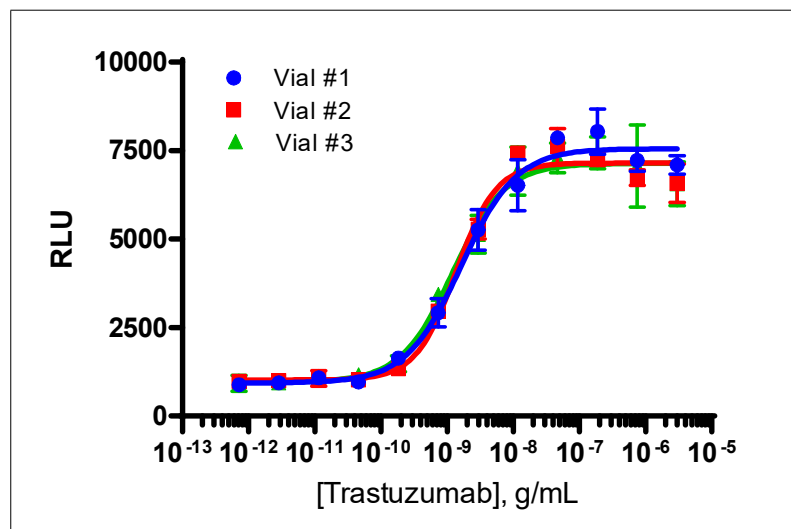
Product Name	KILR NCI-N87 Bioassay Cells
Cryovial Label	KILR NCI-N87 Bioassay Cells
Bioassay Catalog #	97-1004Y021
Bioassay Manufactured Lot #	23B0205
Passage # @ Freezing	3

Assay Information	
Target Species	Human
Cell	NCI-N87
KILR Cell Type	Gastric cancer
CP Reagent	AssayComplete™ Cell Plating 39 Reagent (DiscoverX, 93-0563R39A)
Control Antibody	Trastuzumab
Ligand Diluent	AssayComplete™ Cell Plating 39 Reagent
Detection Kit	KILR® Detection Kit (DiscoverX, 97-0001M)
ADCC Target Cell Number	5,000
ADCC Effector Type	KILR CD16
ADCC Effector Species	Human
ADCC Effector Cell Number	50,000
ADCC Effector to Target Ratio	10:1
ADCC Seeding Time	N/A
ADCC Compound Inc Time	30
ADCC Compound Inc Temp (°C)	37C
ADCC Assay Inc Time	240
ADCC Assay Inc Temp (°C)	37C
KILR Detection Reag. Inc Time	1

Certificate Of Analysis

 Product Name **KILR NCI-N87 Bioassay Cells**

Cell Density Information	
Cell Number (millions)	1
Fill Volume per Vial (mL)	0.1
Cell Viability	
Viability at Initial Thaw (%)	96
Recovery After 72 Hours (%)	152
Mycoplasma and Sterility	
Mycoplasma Test	Passed
Sterility Test	Passed
Functional Performance (3 manufactured vials)	
S:B Ratio	Vial 1 8.2
	Vial 2 6.9
	Vial 3 7.7
EC ₅₀ Check	Vial 1 1.6 x 10 ⁻⁹
	Vial 2 1.4 x 10 ⁻⁹
	Vial 3 1.2 x 10 ⁻⁹



Certificate Of Analysis

Shipping and Storage Information	
Shipping Conditions	Dry Ice
Storage Conditions	Short term (<24 hours): -80°C; Long term (>24 hours): Vapor phase of liquid nitrogen.
Manufacturing Date	February 2023
Expiration Date	February 2026

Shelf life of over 3 years has been established for DiscoverX cell lines and Assay-Ready Cells in general, when stored in the vapor phase of liquid nitrogen.

Documented by / Date: _____

Approved by / Date: _____