

AssayComplete™ Cell Culture Kit 117

Materials Provided

Catalog Number:	92-3117G	
Components:	AssayComplete Cell Culture Reagent 117	500 mL
	Component B	57 mL
	Component G	5.7 mL
Lot Number:	See product label	

Description

AssayComplete Cell Culture Kit 117 is validated to provide an optimal culture environment for propagating DiscoverX KILR® CD16 effector cells.

Product Information

Storage Conditions:	Store at -20°C. Thaw contents at room temperature and mix well by gently inverting the bottle prior to use. Once thawed and mixed, store the final product at 4°C for up to 4 weeks. Avoid multiple freeze/thaw cycles.
Shelf Life:	See product label for expiration date.
Shipping Conditions:	Frozen on dry ice (-70°C).
Instructions on Use:	Refer to the Instructions for Use below and the effector cell user manual for usage.

Notice to Purchaser

AssayComplete Cell Culture Kit 117 has been optimized for use with DiscoverX KILR CD16 effector cells. Occasionally, the reagent may be yellow or pink in color, indicating a slight change in the pH. It has been determined that using discolored media does not adversely impact growth of cultured cell lines or assay performance. Refer to the FAQs section for a detailed explanation of this phenomenon.

For guaranteed performance, use only DiscoverX products optimized with the specific Cell Culture Kit in your assay.

For research use only. Not intended for use in diagnostic or therapeutic procedures.

Quality Control Data

AssayComplete Cell Culture Kit 117 is tested for performance with DiscoverX KILR CD16 effector cells. In addition, each lot is tested for the absence of bacterial, fungal and mycoplasma contaminants.

This product includes high quality serum that has been tested for optimal performance. Testing includes sterility, endotoxin (≤ 5 EU/ml), pH (≥ 6.9 to ≤ 7.8), osmolality (≥ 280 to ≤ 340), total protein (≥ 3.0 to ≤ 5.0), hemoglobin (≤ 10 mg/dl) and absence of mycoplasma, viruses, and bacteriophages.

Instructions for Use

The following procedures are for preparing the AssayComplete Cell Culture Kit.



Refer to the KILR CD16 Effector Cells Certificate of Analysis prior to preparing the culture media.

1. AssayComplete Cell Culture Kit must arrive in a frozen state on dry ice.
2. Thaw individual components of AssayComplete Cell Culture Kit in a 37°C water bath.



For optimal performance avoid longer incubation time. Do not let kit components remain at 37°C after initial thawing.

3. Mix each component separately by gently inverting bottles prior to use under the tissue culture hood.

4. Using aseptic techniques, prepare complete medium by adding the entire contents of Component B and Component G to the 500 mL of AssayComplete™ Cell Culture Reagent. Place a checkmark next to each component.
5. Note down the name of the cells and the date added in the allotted spaces found on the AssayComplete Cell Culture Reagent bottle label.
6. For optimal performance, prepare complete medium using components provided in the kit. In rare instances, precipitates might be observed. Continue with the assay as this does not impact assay performance.
7. Tighten the cap thoroughly to avoid spillage and gently invert the medium bottle several times to mix.
8. Store AssayComplete Cell Culture Kit at 4°C for up to 4 weeks to ensure maximal assay performance.



For cell culture maintenance procedures, please refer to KILR® CD16 Effector Cells user manual.

FAQs

[After the reagent was thawed, the media was discolored, is this normal?](#)

Variations in color of the media reagent can occur. These variations result from CO₂ vapor from the dry ice in the shipping package altering the concentrations of dissolved CO₂ in the media reagents. This can result in slight changes in pH that will induce a yellowing of the phenol red indicator dye present in the media.

We recommend incubating an aliquot of the discolored media in a 5% CO₂ tissue culture incubator until it regains its color, to re-balance the media to its proper pH prior to use in the assay.

For additional information or Technical Support, see contact information below.