

LUCID PLATES

DRY-REAGENT SOLUTION ADAPTED FOR
HIGH-THROUGHPUT LABS

LUCID
DURA Innovations
Now in Microplate
Format



In combination with DURA Innovations technology, Beckman Coulter has expanded its **LUCID - Custom Panel Design & Cocktail Services** offering to include microplates. Now, even high-throughput flow cytometry laboratories can benefit from the advantages that dry-reagent cocktails provide—without needing to adapt their workflow to the tube formats.

These ready-to-use cocktails are tailored to your specific research needs. They are easy to order, easy to use, and free your lab staff from unnecessary, time-consuming tasks.

The result for your lab is a more **streamlined, efficient and productive workflow**.

Specifications

- Two 96-well plates per kit: 2 pouches of 1 microplate each and 1 pouch with Compensation Kit
- Shallow- and deep-well options of microplate format
- 2 to 15 conjugates per panel
- Same reagent panel across all wells
- Minimum order quantity: 2 kits (4 microplates)
- 1 plate of *Evaluation* or *Pre-lot* provided for validation of every new panel ordered
- *DURA Innovations technology*—dry, unitized reagents that can be stored at room temperature
- Barcoded for easy traceability

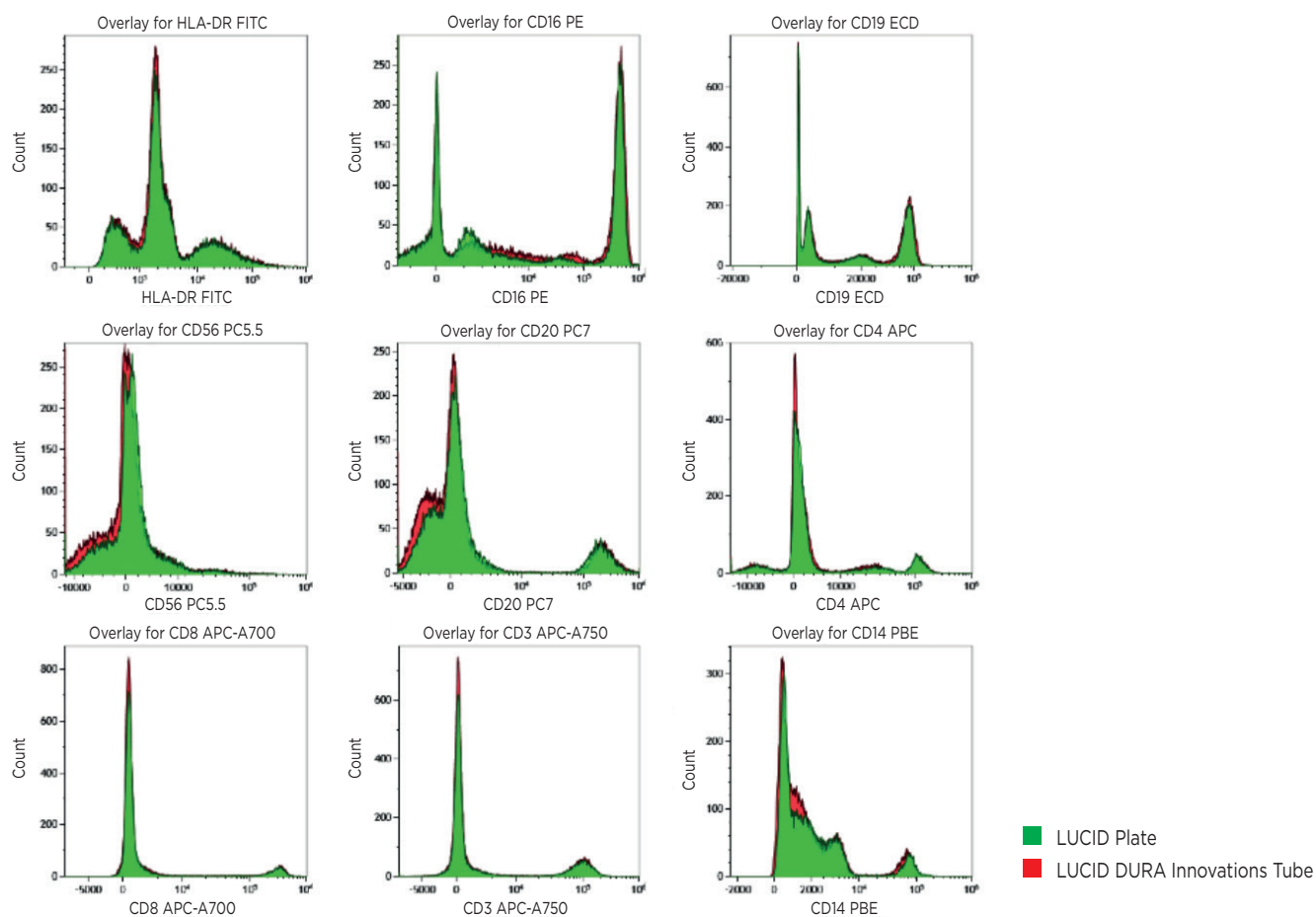


Figure 1: Shallow-well Microplate (330 µL)



Figure 2: Deep-well Microplate (1000 µL)

Function Equivalent to dried customized panel in Tubes



Recruitment Percentage

	HLA-DR FITC	CD16 PE	CD19 ECD	CD56 PC5.5	CD20 PC7	CD4 APC	CD8 AA700	CD3 AA750	CD14 PBE	CD45 KrO
LUCID Plate	20.59	3.09	13.27	2.91	10.9	16.53	9.61	17.81	6.41	93.89
LUCID DURA INNOVATIONS Tube	20.94	3.05	15.36	2.95	12.86	16.16	9.74	18.62	6.2	87.03



Customization need not be synonymous with expensive and time-consuming.

Find out how much we can do for you.

Visit beckman.com/coulter-flow-cytometry/custom-design-services to request a quote.

For Research Use Only. Not for use in diagnostic procedures.



© 2019 Beckman Coulter, Inc. All rights reserved. Beckman Coulter, the stylized logo, and the Beckman Coulter product and service marks mentioned herein are trademarks or registered trademarks of Beckman Coulter, Inc. in the United States and other countries.

For Beckman Coulter's worldwide office locations and phone numbers, please visit "Contact Us" at beckman.com

FLOW-5452FLY05.19