



# CUSTOM DESIGN SERVICES

YOU CAN DO MORE BECAUSE WE DO MORE FOR YOU.



**Custom Panel Design  
& Cocktail Services**

With a spectrum of multicolor options for you to choose from, our panel design experts can create a custom cocktail to clearly deliver the information you seek.

**EVERY**  
*event matters.*

 **BECKMAN  
COULTER**  
*Life Sciences*

# Do more research. Do more data analysis. Do more testing. Do more of what you're passionate about—and less manual cocktailing.

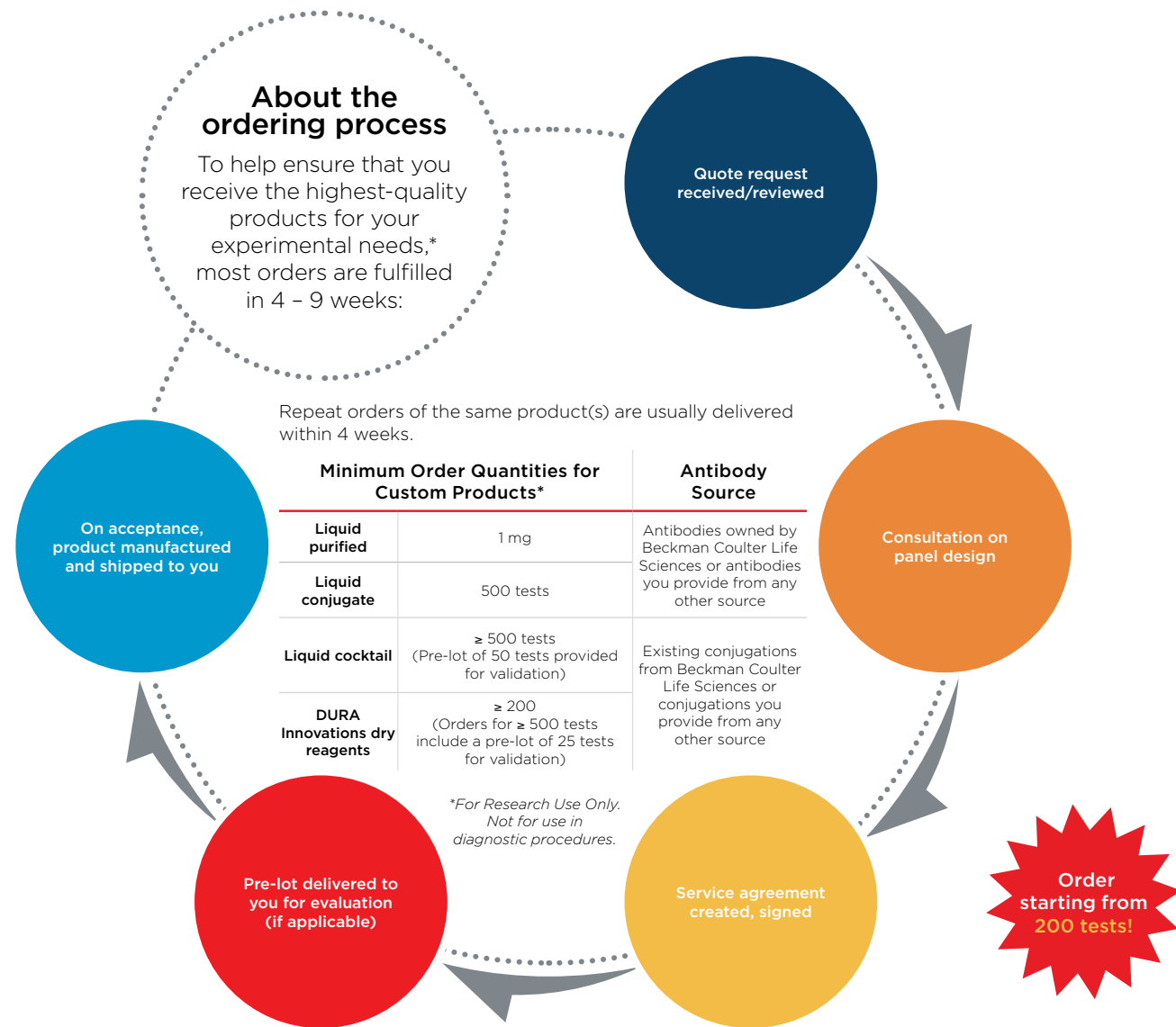
Whatever you want to do more of, we can help. Because we can do more for you.

We can be a time- and cost-saving, single-source provider of off-the-shelf and customized products and services to support all your flow cytometry needs.

With our flexible, collaborative approach, small lot size requirements, and waste-reducing DURA Innovations-based dry reagents, we can help you do more, save more and achieve more.



Save time and reduce the potential for human errors from manual cocktailing.



## If we don't currently offer the product you need, our panel design experts can create an optimized panel that clearly meets your specific criteria.

Whether you need a different conjugate of one of our 630+ ASR-labeled human antibodies—or a custom conjugation of one of your own—we can provide affordable options for all your experimental design requirements:

- Single-color conjugations
- Optimized multi-conjugate combinations
- Bulk purified antibodies (you specify concentration and buffer)
- Exclusive DURA Innovations dry (not lyophilized) custom reagents
  - Stable at room temperature (18°C - 23°C)
  - Warranted for 1 year
  - Include lot-matched compensation kits



## Frequently asked questions

### How long does it take to get a custom order?

It depends on availability of raw materials, but we can usually deliver single-color conjugates in 4 - 6 weeks, and more complex multicolor reagents in 9 weeks to 12 months. With advance notice, repeat orders can be delivered in 4 weeks.

*If antibodies required for an order are unavailable in either a purified state or as ascites fluid, delivery may take longer.*

### Will we get a pre-lot of our custom product to evaluate?

Yes. With a minimum order of 500 tests, you'll get a pre-lot of 25 tests for evaluation.

### Will we receive a compensation kit with our order?

Yes. We include a compensation kit with every 25 tests,\* regardless of order quantity.

### Can my custom product be labeled as CE/ASR?

No. Our CDS products are designed and manufactured in facilities whose quality systems have been certified as compliant with the ISO 9001 and/or ISO 13485 quality system standards. However, unlike our ASR and CE-IVD products, our CDS products are intended for research use only, and are not intended to be used in diagnostic procedures.

### Can I obtain sodium azide-free formulations?

Yes. Purified antibodies can be formulated without azide for use in conjugation and functional studies, as they can be stored frozen to minimize risk of microbial contamination. Or conjugates can be formulated at high concentrations with low amounts of sodium azide.

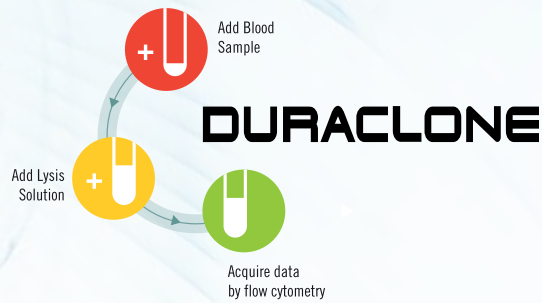
### Is there a performance difference between liquid reagents and dry DURA Innovations reagents?

No. The proprietary DURA Innovations process yields a dry, stable product that delivers the same performance as a liquid product.

**Expedite and simplify  
your sample preparation.**



**Custom Panel Design  
& Cocktail Services**



*For DURA Innovations products, you can also save money by reducing reagent waste. Visit [Contact](#) your local **Beckman Coulter** sales rep to understand the wastage for your specific lab to create a detailed cost-savings report specifically for your lab*

**What is DURA Innovations?**



### **DRY**

- Drying creates a uniform reagent layer at the bottom of the tube.
- DURA dry reagents require no refrigeration.
- You can produce a single batch and store it at room temperature for the study duration.

### **UNITIZED**

- Custom panels are supplied in a single-use cocktail per tube/plate, ready for addition of sample.

### **REAGENT**

- Stable reagent layer at the bottom of the tube helps ensure optimum tube-to-tube and batch-to-batch consistency.
- Cocktails can include calibration beads for absolute cell counting (optional).

### **ASSAYS**

- Conjugates are suitable for:
  - Flow cytometry
  - Microscopy
  - Immunoassays

**Customization need not be synonymous with expensive and time-consuming. Find out how much we can do for you. Visit [www.beckman.com/coulter-flow-cytometry/custom-design-services](http://www.beckman.com/coulter-flow-cytometry/custom-design-services) to request a quote.**

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



© 2017 Beckman Coulter Life Sciences. 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](http://beckman.com)  
FLOW-2597SB04.17