



**THE ONLY RESOURCE YOU NEED  
FOR ANTIBODY COCKTAILS IN BULK.**

YOU DESIGN IT. WE MAKE LOTS OF IT. SO IT'S THERE WHEN YOU NEED IT.



**Contract  
Manufacturing  
Services**

**EVERY**  
*event matters.*



# Introducing one resource you need if you want to optimize your lab's productivity.

It's your singular resource for bulk orders of made-to-specification antibody cocktails\*—from experts in flow cytometry.

CMS options include:

- Antibodies from any source (ours, yours, or our competitors)
- Dried or liquid formats
- Exclusive DURA Innovations
- Flexible delivery terms
- QC testing on request (per your validation method and specified acceptance criteria)

## Example of Typical Workflow vs

10 Specimens with single-color panels from receipt to first results

> 25 Reagent vials



Re-Titration  
Re-Validation

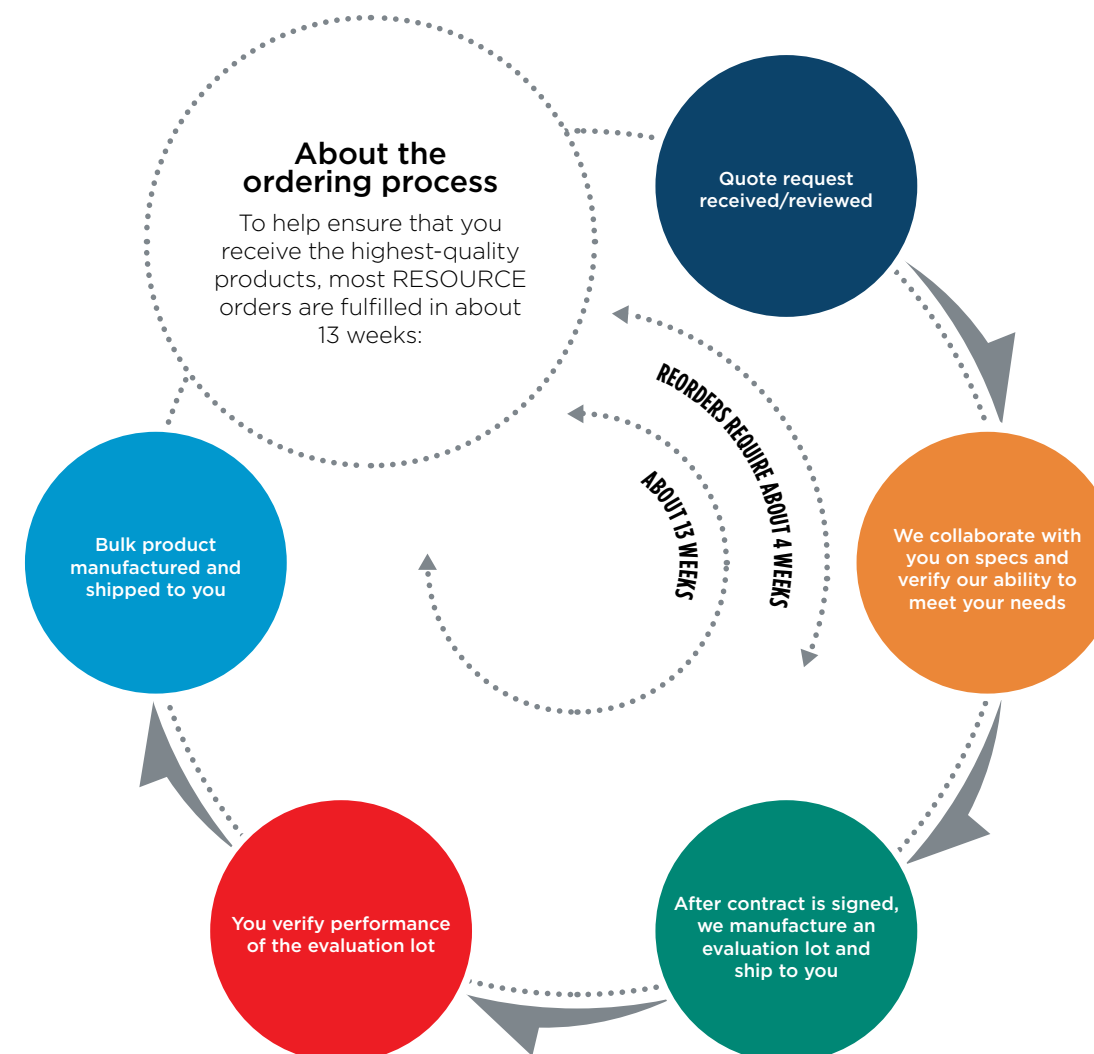


= 4 Hours



= 0 Hours

You can see an instant savings by eliminating labor costs for creating large quantities of cocktails. 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.](#)



## Why choose Beckman Coulter Life Sciences as a trusted RESOURCE for bulk antibody cocktails?

- Your cocktails will be manufactured following GMP\*\* in ISO 9001-compliant facilities.
- Your order will include a complimentary evaluation lot for you to validate and approve.
- You can optimize QC by minimizing the potential for errors/variability.
- You can save time by simplifying inventory management and streamlining your workflows.

## Frequently asked questions

### Can any lab use this contract service?

Yes. It's available to clinical flow labs, contract research organizations, biopharmaceutical companies that perform clinical trials, or any facility that performs the same test multiple times each day.\*

### Can I use this service if Beckman Coulter Life Sciences doesn't offer the antibodies I need?

Yes. You simply provide us with the antibodies and instructions. (Note: volumes required for some formulated reagents from other vendors may preclude use in dry-down formats.)

### Can you manufacture ASR- and/or IVD-labelled products for us?

No. We will take your proprietary cocktail and manufacture it to your specifications following good manufacturing practices. As part of this strictly confidential process, you'll receive your products labelled only with:

- A brief description
- Unique part and lot numbers
- Volume
- Any applicable warning

### Will we get a sample of our product to evaluate before committing to manufacturing it in bulk?

Yes. After you give us your panel formulation, we'll provide you with 25 tests that you can evaluate for up to 4 weeks.

### Can you assist us with titrations?

No. Products contracted under this service are manufactured to your specifications. Your laboratory should validate the products as necessary to meet your specific needs.

### Can we use this service even if we don't have a flow cytometer from Beckman Coulter Life Sciences?

Absolutely.

### Can we establish a delivery schedule that meets our needs?

Yes. After you've approved and received your first order, we'll work with you to create a suitable delivery schedule for recurring orders of the same product.

\*These products have no regulatory status. Customer is solely responsible for validating them  
\*\* GMP as defined under India Drugs & Cosmetic Act 1940.



Contract  
Manufacturing  
Services

## What is DURA Innovations?



## DURA Innovations

Dry Unitized Reagent Assays

### 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

## Be more productive by being resourceful.

Visit [www.beckman.com/coulter-flow-cytometry/reagents/custom-reagent-services](http://www.beckman.com/coulter-flow-cytometry/reagents/custom-reagent-services) today to request a quote.

*\*This service is currently available only in the United States. If your lab is outside the U.S., consult with your local Beckman Coulter Life Sciences representative about your interest in making this service available in your country.*



Products under DURA Innovations are intended for both RUO and IVD. Kindly refer to the product specifications for more information.

© 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-2598SB04.17