

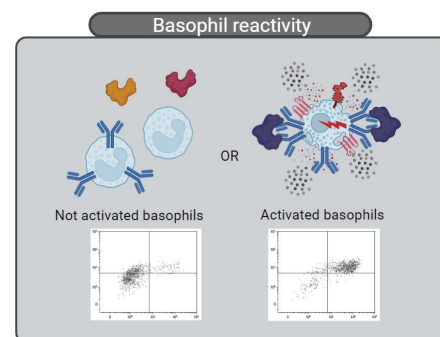


# Coming soon!

## Introducing the Next-Generation Basophil Activation Test (BAT)

### What is a Basophil Activation Test (BAT)?

BAT are functional assays measuring basophil activation in response to specific allergens or other stimuli



### Leveraging dry technology to overcome traditional BAT challenges



- **Fostering standardization:** Conjugated antibodies and allergens are pre-mixed at desired dose, and dried down together in a single tube. Negative and positive controls can be included.
- **Straightforward:** add your sample, vortex and acquire your data.
- **Easy to use:** no compensation, no centrifuge, compatible with any flow cytometer.
- **Customizable:** define your own combination of antibodies, allergens with specifications and format of your choice.

### A promising outlook for BAT

MEDIA ROOM | October 13, 2022

#### FARE Innovation Award Diagnostic Challenge Presented to Beckman Coulter Life Sciences

\$1,000,000 Awarded to Advance Innovation in Food Allergy Diagnosis

Our Next-Generation BAT technology led to the solution being selected as the most promising approach to create an accurate, accessible, user-friendly test for the study of food allergies by FARE, the US leading non-profit organization engaged in food allergy advocacy.

**Coming soon! Available from end Q1 2025**  
**Reach out to your BEC LS sales representative to learn more**

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



© 2024 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. All other trademarks are the property of their respective owners.

For Beckman Coulter's worldwide office locations and phone numbers, please visit Contact Us at [beckman.com](https://www.beckman.com)  
2024-GBL-EN-106664-V1

