



MANAGE REGULATORY NOTIFICATION

- Any notification awaiting a response will be listed under My Open Notifications. Click on **View** for more details for each notification.

- From here you can open and print the notification letter as well as any other attachments.

Subject: FAN Project TESTING Customer Notification >> IMMEDIATE ACTION REQUIRED

Beckman Coulter is initiating a product notice FA-700-7177-A Launch for External website.

Please take the below listed actions:

- Download and view the attached notice (PDF).
- Download and view the Replacement Order Form (if applicable).
- Please acknowledge by selecting Respond.

IT IS MANDATORY THAT YOU COMPLETE THE RESPONSE WITHIN 10 DAYS.

Attachments

[130-UAT-FAN-intFANint-001.xlsx](#)

- Select the site(s) you are responding for in the box to the left of **Site Number** and click on **Respond**.

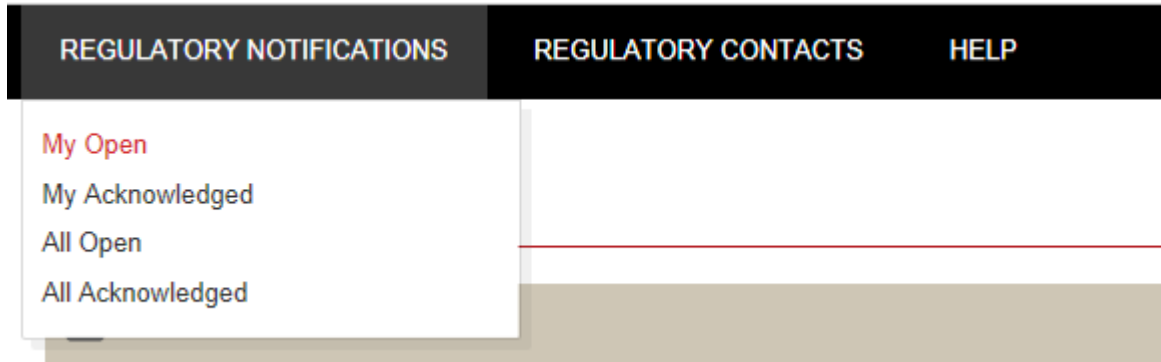
Upon closure of this field action by Beckman Coulter the regulatory notification will be removed from your view

Site Number	Address Line1	Address Line2	City	Postal Code	State/ Province	Country
<input type="checkbox"/> > 300	FERRY HILL RD	MEDICAL CENTER	MONTGOMERY	36109-3725	Alabama	United States
<input type="checkbox"/> > 300	LOOP RD E	MEDICAL CENTER 679	TUSCALOOSA	35404-5099	Alabama	United States
<input type="checkbox"/> > 300	HOSPITAL RD	MEDICAL CENTER 680	TUSKEGEE	36083-5001	Alabama	United States

- Select either one or both of the options. Click on Save.

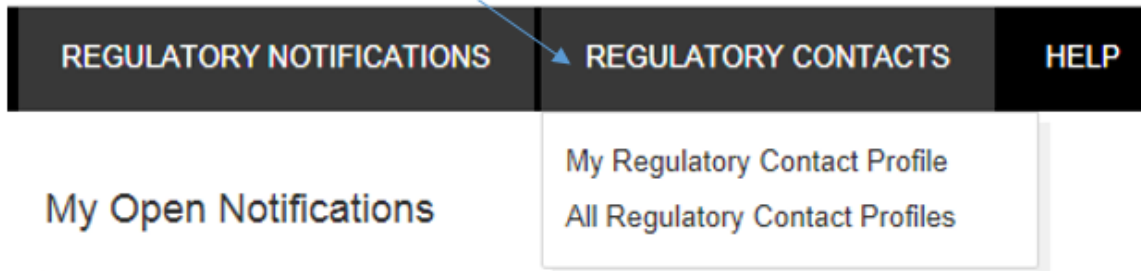


- 5. You will then be taken back to the **My Open Notifications** page. As long as there are open notifications you may keep responding for the sites you are responsible for. When all are answered for, no further notifications will be listed. You may also view your acknowledged or other acknowledged notifications for your site(s). Hover with the mouse cursor over **Regulatory Notifications** and click on the one you wish to view.



MANAGE REGULATORY CONTACTS

- 6. You can also maintain Regulatory Contact information.



- 7. Under My Regulatory Contact Profile you may update your Contact Details, Opt-in Products and Opt-in Locations. Click on **Add** to add products and/or locations you want to be notified of; click on the trash can to remove products and/or locations.

Regulatory notifications opt-in

To allow Beckman Coulter to provide this contact with regulatory notifications, please opt-in the contact for a product AND location below.

Opted-In Products [Add](#)

Product	Actions
All Products	

Opted-In Locations [Add](#)

Country	State	City	Actions
United States	AL	BIRMINGHAM	



8. If you have made changes to the **Contact Details**, click on **Save** to keep your changes, or **Cancel** to revert back to the previously stored information. If you don't click on Save, you will lose any of the changes made.

My Regulatory Contact Profile

Cancel **Save**

▼ Contact Details

Customer Name U S VETERANS ADMINISTRATION

*** Contact First Name** Mr.

9. You can also view All Regulatory Contacts currently registered with your account, as well as update their info and opt-ins, and create new contacts.

Contact First Name	Contact Last Name	Contact Status	Job Title	Contact Type	Customer Name	
YAP		Active	Lead Technologist	Regulatory Contact	U S	View
ADA		Active	Chemistry Super...	Regulatory Contact	U S	View

Clicking on the plus (+) sign next **Filters** allows you to search by **Country** and **State/Province**.

Filters

Country State/Province City

[Search](#)

10. Under Help you will find this Tutorial as well as information on how to contact us.

REGULATORY NOTIFICATIONS **REGULATORY CONTACTS** **HELP**

[Tutorial](#)
[Contact Us](#)