

# User guide

## Testing service portal for AerosolSense Sampler

Accessing portal, logging orders, shipping sample cartridges, and reviewing results

01

### ACCESSING PORTAL

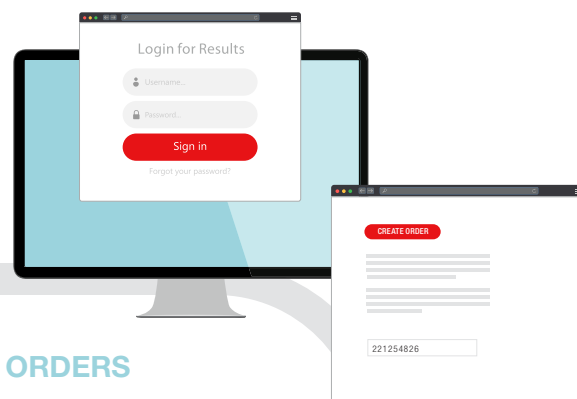
- **If you are a system admin:** when testing service is purchased, you will receive an email to login to your GoMeyra admin account
- Once you log in, add the users (under the “user” tab) that you want to have access to the portal



02

### LOGGING ORDERS

- Log in to your GoMeyra portal
- Click “Instruments” on the left toolbar, then select the instrument you’re entering an order for
- Click “Create Order”. Submit an order for each sample cartridge ID shipped to the lab
- Enter the barcode serial number from your sample cartridge
- If you would like to track your shipment, enter the shipping label barcode



03

### SHIPPING CARTRIDGE

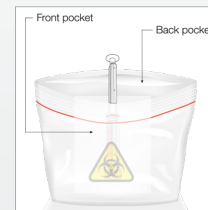
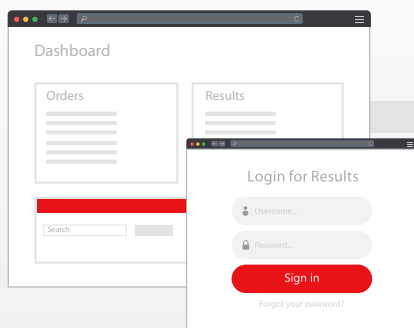
- Place the sample cartridge in the pocket of the biohazard bag containing the absorbent pad. Cap must be on cartridge
- Put the bagged and sealed sample cartridge into the shipping box provided by testing service, and place shipping label on box
- Drop off at the nearest shipping location



04

### REVIEWING RESULTS

- Log in to your GoMeyra portal
- Results will be posted at top of dashboard within 24 hours
- Export your results



For further support, visit [www.thermofisher.com/apssupport](http://www.thermofisher.com/apssupport)