

Application acknowledgement

Upon successful application to this course, you will receive an automatically generated email acknowledging your application. If you do not receive this email within two business days of applying, please email sartorius.academy@sartorius.com to confirm receipt of your application.

Application confirmation

Should your application be confirmed, you will receive an email by September 10, 2021 to confirm your application and to inform you on the payment and registration process.

Payment and Registration

Please note that we accept payments only from your company or institute via PO and not via online payment gateways. The payment is due within 10 days of application confirmation.

On confirmation of application, the applicant is responsible to complete registration by sending the PO and ensure payment is completed within 10 days of application confirmation.

If no payment has been received within 10 business days of application confirmation, Sartorius reserves the right to cancel registration.

Registration is subject to our General Terms and Conditions of Business.

Once registered, confirmed course participants will receive an email with detailed information regarding course logistics, platform (MS Teams), and a short pre-course questionnaire. If you do not receive this email before Sep 15, please email sartorius.academy@sartorius.com

Sartorius will not be responsible for inability of participant to join the course due to problems with internet connectivity and/ or the learning platform (MS Teams).

Any additional questions prior to payment and registration should be directed to sartorius.academy@sartorius.com

Change of Scope of Services and Cancellations

Cancellation by learner

Cancellations from learners are not accepted. However, registered participants always have the option to appoint a substitute to attend.

SARTORIUS

Substitutions may be made up to two business days prior to the course start date. Notify us by sending an e-mail to sartorius.academy@sartorius.com including the name and email contact for the replacement from the same company or institute.

Cancellation by us

We reserve the right to modify course, material, or coaches, while preserving the overall nature of the course or to limit course registration. In certain circumstances, we may even need to cancel the course. Registrants will be notified by Sartorius if a course is cancelled. We will offer in these situations an alternative date/ course to attend or full refund of registration fees paid. We will not be responsible for any other costs incurred. If we need to cancel the course due to acts of God, or other adventitious events, the registrants will have no claim.

Copyrights

The information is provided for educational and informational purposes only and should not be construed as legal advice. It is intended for the sole purpose of individuals participating in this course. Any republication, retransmission, reproduction, downloading, storing or distribution of all or part of any materials found here is expressly prohibited unless obtained with the written consent of Sartorius.