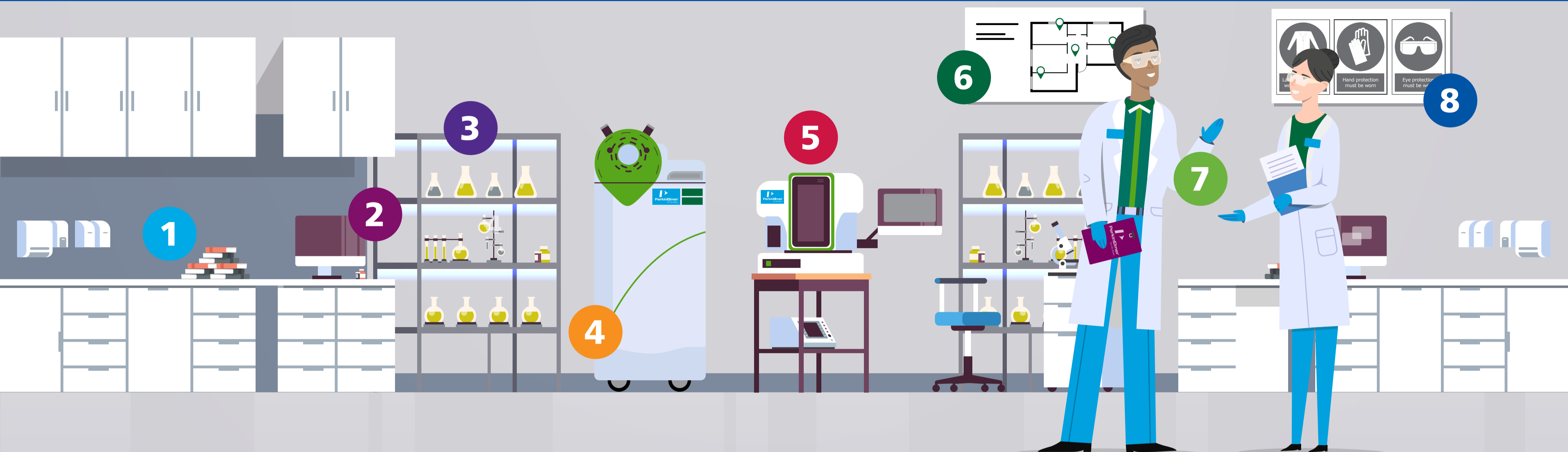


# LABStartUp Program Guide



1

## Training and education

We offer education and training that helps you unlock the full potential of your scientists and your instruments.

[LEARN MORE](#)

2

## Automation and data management

We deliver automated processes and analytics between the laboratory and operations.

[LEARN MORE](#)

3

## Reagents and consumables

Our Lab StartUp program provides numerous benefits, including up to 30% off your first consumables order.

[LEARN MORE](#)

4

## Maintenance

Preventative maintenance plus our multivendor approach that helps incorporate any vendor's instrumentation into your lab.

[LEARN MORE](#)

5

## Instrumentation

You benefit from significant financial assistance, including grant-matching, discounts, and special financing options – plus free advice and quotes.

[LEARN MORE](#)

6

## Ergonomics and setup

Our asset location and optimization services are delivered through our skilled OneSource Laboratory Services team.

[LEARN MORE](#)

7

## Customized support

OneSource support and maintenance for your lab startup with customized solutions – so you can focus on your core scientific activities.

[LEARN MORE](#)

8

## Safety

Ensure that post-COVID return-to-work assessments are performed and standard operating procedures are in place.

[LEARN MORE](#)

[APPLY TO LAB STARTUP PROGRAM HERE](#)