



Frequently Asked Questions

Q What is the purpose of the Lab StartUp Program?

A The Lab StartUp program has been developed to help you maximize your new lab's performance while optimizing spend. [Register now](#) to enjoy incentives and specials including grant matching, discounts, promotional items, resources, and tips on select instruments and consumables. As well as access to our expansive scientific and technical expertise.

Q Where do I sign up for the Lab StartUp Program?

A Complete the [Registration form](#) on the Registration page to sign up.

Q Who is eligible for the Lab StartUp Program?

A To be eligible for the Lab StartUp Program, you must have done within the last 12 months or will be planning to do in the next 12 months one of the following:

- Opened or are opening a new lab within your institution
- Opened or are opening a new lab at different institution
- Starting a new company
- Moving your current lab into a new space
- Received your first research grant

Upon registering for the program using the [Registration form](#), you will be contacted shortly thereafter confirming eligibility.

Q I am not sure if I am eligible for the Lab StartUp Program?

A Complete the [Registration form](#) and the PerkinElmer Lab StartUp Program team will review your application and contact you shortly thereafter with your eligibility.

Q What products does the Lab StartUp Program support:

A The Lab StartUp Program involves a wide range of PerkinElmer instruments, accessories and consumables. Specific incentives within the program are for a Limited Time, we recommend registering soon to maximize access incentives and specials including discounts, promotional items, resources, and tips on select instruments and consumables.

Q How will I know if I have meet eligibility for the Lab StartUp Program?

A You will be contacted by a member of the PerkinElmer program team shortly after registration with eligibility

Q How will I know what specials and/or promotions I am eligible for in the Lab StartUp Program?

A The Lab StartUp Program involves a wide range of PerkinElmer instruments, accessories and consumables. Specific incentives within the program are for a Limited Time, for specific products, and/or specific geographies. A PerkinElmer program team contact will guide you to maximize your benefits. We recommend [registering](#) soon to maximize access incentives and specials including discounts, promotional items, resources, and tips on select instruments and consumables.

Q Is the Lab StartUp Program available in my country?

A The New Lab StartUp program is available in most countries where permissible by law. Upon registering for the program using the [Registration form](#), you will be contacted shortly thereafter on eligibility and any conditions.

Q Can I purchase multiple products using the Lab StartUp Program?

A Yes, you can purchase multiple products

Q Can I combine the Lab StartUp Program specials with other promotions?

A New Lab Start up Program and Incentives and other promotions cannot be combined

Q What is the duration of the Lab StartUp Program

A Currently there is no end date to the PerkinElmer New Lab StartUp program. Specific incentives within the program are Limited Time promotions and can end at any time. PerkinElmer reserves the right to end or change the program and promotion within the program at any time without notification. We recommend [registering](#) soon to maximize access incentives and specials including discounts, promotional items, resources, and tips on select instruments and consumables.

Q Is there cost or commitment on my part for registering for the Lab StartUp Program

A The New Lab StartUp Program is complimentary; nor are there any commitments on your part. The New Lab StartUp program has been developed to help you maximize your new lab's performance while optimizing spend.