
Application form

We ask that you submit NON-CONFIDENTIAL INFORMATION ONLY.

All information you provide in this site will be governed by our Privacy Policy.

For questions or comments about the application process and form submission, please contact **Young.Investigator@qiagen.com**.

Instructions:

Please answer the following questions below and provide any supporting images that you'd like to have uploaded – 3 to 5 images would be optimal including one headshot.

Accepted file formats: .jpg and .tif of good quality and resolution, in landscape orientation only.

Name:

Affiliation:

Academic status/Position held:

Project title:

Email:

1. Tell us about your background and how you became interested in forensic science?

2. Can you provide a summary of the project you are working on?

3. Please describe your typical day in the lab.

4. What do you find most interesting about your project? Have you seen any surprising results?

5. What are the benefits of your project?

6. What are the major challenges faced while working on your project and how do you overcome them?

7. Which QIAGEN products do you use and what do you like about the products?

8. Outside of forensic science, what are your hobbies?



Please sign below indicating that you have read and agree to the use of your data and photo for QIAGEN's Young Investigator Blog. Without this signature we are unable to process your submission.

VIDEO/PHOTOGRAPH RELEASE FORM

I hereby grant QIAGEN and its affiliates the irrevocable right and permission to use photographs and/or video recordings of me arising from my participation in the QIAGEN's Young Investigator Blog on QIAGEN's websites and in publications, promotional flyers, educational materials, derivative works, or for any other similar purpose without compensation to me. I agree that all such portraits, pictures, photographs, video and audio recordings, and any reproductions thereof, and all plates, negatives, recording tape and digital files, are and shall remain the sole property of QIAGEN.

I understand and agree that such photographs and/or video recordings of me may be placed on the Internet. I also understand and agree that I may be identified by name and/or title in printed, Internet or broadcast information that might accompany the photographs and/or video recordings of me. I understand and agree that I will not have an opportunity to review the final products before they are published or distributed.

I hereby release, acquit and forever discharge QIAGEN, and its current and former trustees, agents, officers, employees, and affiliates, from any and all claims, demands, rights, promises, damages and liabilities arising out of or in connection with the use or distribution of said photographs and/or video recordings, including but not limited to any claims for invasion of privacy, appropriation of likeness or defamation.

I acknowledge that my participation in the photo and video portions of the QIAGEN's Young Investigator Blog is completely voluntary, as is my consent to this release. I understand that no action will be taken against me should I decline to participate in the photo/video portions of these events.

This release is binding on me and my heirs, assigns and personal representatives.

Agreed and accepted:

Signature of Participant

Date

Printed Name of Participant