



Poster Submission

Submit a Poster

- Your abstract must be submitted by **August 31st, 2019**, to be included in symposium materials. Late posters will not be listed in symposium materials.
- Please complete the attached poster submission form, and email to symposium@genscript.com.
- An approval email will be sent to you within one week of your submission.

Benefits

- Receive complementary registration
- Gain extra exposure of your research work

Criteria

- Scientific and relevant to symposium theme
- Clearly presented data

On-site Information

- All posters should be portrait orientation, with maximum dimensions of 36 inches wide (3 feet) by 48 inches tall (4 feet).
- Poster locations are available at a first-come, first-serve basis.
- Presenters must be available to present during poster viewing time.
- Posters must be set up by 11am, and taken down by 6pm.
- You are responsible for printing, transportation and logistics of your poster.
- 4 posters will be selected to present 5 minute lightning talks.

Contact Information

Name of Person Submitting Poster Form:

Company/Organization:

Phone:

Email:

Poster

Poster Title

By

(Include all co-authors and their affiliations.)

Presented By

(List the person who will present the poster at the symposium.)

Abstract (attach additional sheets if necessary)

DISCLAIMER: GenScript reserves the right to publish your poster title and abstract, in whole or in part, in any Symposium marketing materials or products, including but not limited to conference brochures, websites, program books, speaker presentation links, videos, social media posts and Symposium satellite events.

By submitting content, you hereby represent and warrant that you have the required authority to submit said content and that it does not infringe upon the patent, trademark, copyright, or other intellectual property rights of a third party.

INQUIRIES: Please direct poster submission inquiries to symposium@genscript.com.