



Infectious Diseases | Tuberculosis

2021 Global TB Clinical and Educational Summit

Poster
competition
Win an Apple®
iPad®

Poster submission information

All summit attendees are invited to take part in the poster competition. Posters must be submitted by a registered participant. Please submit a copy of your abstract/outline using the form at: www.qiagen.com/TB-Summit-2021 by Friday, September 17.

Abstracts

Abstracts or outlines may contain:

- Authors/Institution name
- Poster or project title
- Background/objectives
- Methods
- Results
- Conclusions

Posters

Final posters should be submitted (in PDF format) to events@qiagen.com by Tuesday, September 21, 2021 for upload onto the event platform.

- There are no restrictions regarding the number of words, font size, design and imagery.
- Participants may submit more than one abstract/poster.
- No oral presentations will take place, but we encourage you to be available in the networking lounge during session breaks to discuss your poster with attendees.
- Posters will be voted upon by summit attendees during the event and the winner will be announced at the end of the summit. Each registered attendee can vote only once.
- An Apple® iPad® will be awarded for the poster with the most peer votes.

Submission and deadlines

- The deadline for abstracts/outlines is Friday, September 17, 2021 via the form at www.qiagen.com/TB-Summit-2021
- The deadline to submit final poster PDFs is Tuesday, September 21, 2021. Send your poster PDF to events@qiagen.com. Posters will not be accepted without an abstract.

Notification

Submitting authors will be notified of acceptance.

Individuals or entities involved in the provision of health care services and/or items to patients, which purchase, lease, recommend, use, arrange for the purchase or lease of, or prescribe medical technologies, or who are classified as healthcare professionals or healthcare organizations under any applicable laws or industry codes of ethics, may not participate in the poster contest.

Presenters are responsible for confirming that participation is compliant with the internal policies of their employer.

Contact

Email: events@qiagen.com



Trademarks: QIAGEN®, Sample to Insight® (QIAGEN Group); Apple®, iPad® (Apple Inc.). Registered names, trademarks, etc. used in this document, even when not specifically marked as such, may still be protected by law.

PROM-18855-003 © 2021 QIAGEN, all rights reserved.

Ordering www.qiagen.com/shop | Technical Support support.qiagen.com | Website www.qiagen.com