



## Q&A ATCC® *Excellence in Research* Webinar “Microbiological Quality Control as Described in the Compendia”

### General Questions

1. Will we be able to download the presentation?  
This presentation will be available to watch on demand on the ATCC website, or [click here](#).
2. What percentage of glycerol is used for the ATCC® Minis stocks?  
ATCC uses 10% glycerol for our ATCC® Minis.
3. Does ATCC have ATCC® Minis that match my VITEK® 2 cards?  
Yes, ATCC plans on releasing strains later in the spring of 2015 in the ATCC® Mini format that are aligned with VITEK® 2 cards from bioMérieux.
4. The ATCC website says to store ATCC® Minis at -80°C. Can I store them at -20°C?  
Our real-time testing has so far demonstrated that ATCC® Minis are viable up to 2 months at -20°C. We plan on continuing this testing through 1 year to assess long-term viability at this temperature.
5. I currently purchase preceptrols. Why are the ATCC® minis more expensive?  
With ATCC® Minis, you are receiving 6 ready-to-use vials of a strain per order instead of one vial per order. This brings the unit price per vial to approximately \$38, which is comparatively less expensive than the \$60 per vial for each preceptrol. What’s more, because ATCC® Minis are provided in a ready-to-use glycerol stock format, you can do away with the time and expense associated with strain propagation and preservation, and go straight to your research.
6. Can a -70°C freezer be used to store ATCC® Minis?  
Yes, -70°C is an appropriate temperature to store ATCC® Minis.
7. What are the applications of ATCC® Minis?  
ATCC® Minis are ideal for quality control analyses. A number of our strains in the ATCC® Minis format are the same quality control strains recommended for use in the United States Pharmacopeia, European Pharmacopeia, and Japanese Pharmacopeia. Further, our newest products include strains that have applications in food testing analyses.
8. Can ATCC® Minis be used to make in-house control stocks?  
ATCC® Minis put an end to do-it-yourself banking with each six-pack of ready-to-use quality controls strains in glycerol stock. This new product format is the same frozen format that you

use in your own laboratory, so there is no need to make in-house control stocks. Just open, plate, and go!

9. What passage number are ATCC® Minis?

ATCC® Minis are at passage zero.

10. Is there data to support phenotypic or genotypic changes after 5 passages?

We recommend no more than 5 transfers or passages be made from an ATCC Genuine Culture®. Repeated passage or serial culture increases the risk of contamination and can result in undesirable genotypic and/or phenotypic changes. The recommendation of five passages or fewer from the ATCC reference culture is broadly accepted in the healthcare community and the pharmaceutical and biopharmaceutical industries. U.S. Pharmacopeia Chapter 51: Antimicrobial Effectiveness Testing. U.S. Pharmacopeia National Formulary, 2005 states: "The viable microorganisms used in the test must not be more than five passages removed from the original ATCC stock." ATCC agrees with this recommendation.

©2015 American Type Culture Collection. The ATCC trademark and trade name, and any other trademarks listed in this publication are trademarks owned by the American Type Culture Collection unless indicated otherwise. VITEK® 2 and bioMérieux are registered trademarks of bioMérieux, Inc. These products are for laboratory use only. Not for human or diagnostic use. ATCC products may not be resold, modified for resale, used to provide commercial services or to manufacture commercial products without prior ATCC written approval.