



INSTRUCTIONS FOR USE

ATCC® MICROQUANT™ LOW CFU KIT

DESCRIPTION

ATCC® MicroQuant™ strains are precisely quantitated controls in a single-use format that rehydrates immediately, requires minimal handling, and is convenient to store. The low CFU kits comprise five vials of an ISO 17034 reference material and five vials of rehydration buffer. Once rehydrated, each cryopreserved preparation provides reference cultures with 100 to 1000 CFU per 1.0 mL (10 to 100 CFU per 100 µL).

INTENDED USE

This kit is intended for research use only (RUO). It was developed to support USP General Chapters.

PRECAUTIONS

- Aseptic procedures are recommended while handling this kit.
- The product should be handled under the proper biosafety level conditions. ATCC determines the biosafety level of a material based on our risk assessment as guided by the current edition of *Biosafety in Microbiological and Biomedical Laboratories (BMBL)*, U.S. Department of Health and Human Services. It is your responsibility to understand the hazards associated with the material per your organization's policies and procedures as well as any other applicable regulations as enforced by your local or national agencies.
- Biohazard disposing procedures must be followed while discarding kit contents.
- The rehydration buffer contains the animal by-product sodium chloride peptone. The buffer was formulated to meet EU, JP, and USP guidelines (concentration of animal origin products <1% w/v).
- The rehydration buffer for spore-forming fungal organisms (ATCC® 16404-LQ-PACK™, ATCC® 10106-LQ-PACK™, and ATCC® 11709-LQ-PACK™) is formulated differently than the other low CFU kits and does not contain animal by-product.

RECOMMENDED STORAGE CONDITIONS

ATCC® MicroQuant™ Low CFU kits can be stored at 2-8°C for 6-12 months. Verify the expiry for individual products on the respective certificate of analysis.

LIMITED LABEL LICENSE FOR MICROQUANT™

Use of ATCC's MicroQuant™ products are subject to the ATCC Material Transfer Agreement. Notwithstanding, ATCC hereby grants you the further right to use this single-use consumable MicroQuant™ product as a reagent for a single quality control assay in your laboratory only, whether as a commercial service or internally. ATCC's MicroQuant™ products contain trade secrets and intellectual property of ATCC, and you may not reverse engineer, replicate, alter, or tamper with the products or authorize any third party to do any of the foregoing. ATCC's MicroQuant™ products shall not be used for any other use restricted by the ATCC Material Transfer Agreement, including any animal or human therapeutic use, any human or animal consumption, any diagnostic use, or use to manufacture any products.

For information on obtaining additional rights, please contact ATCC Licensing at licensing@atcc.org.



INSTRUCTIONS FOR USE

1. Remove the desired number of cryopreserved reference culture vials (gray crimp) and rehydration buffer vials (green crimp) from 2–8°C storage and place into a biosafety cabinet. One vial of rehydration buffer should be used for each vial of reference culture.
2. Uncap each vial. Add 1.0 mL of rehydration buffer to the reference culture vial. Recap the vial and allow the pellet to dissolve for 30 seconds.
3. Vortex on the high setting for 30 seconds to ensure the suspension is well mixed. The suspension is now useable and contains 100 to 1000 CFU per 1.0 mL (10 to 100 CFU per 100 µL). The suspension can be used immediately or stored at 2–8°C for up to 8 hours. Note: Before plating, gently pipette the suspension up and down five times to ensure thorough mixing. If the suspension has been left standing for more than five minutes, or if it was stored at 2–8°C after rehydration, vortex the suspension again to ensure uniformity prior to use.
4. Uncap the vial and pipette 100 µL of the resuspended reference culture directly on a non-selective media agar plate. If following a specific pharmacopeial chapter, use the appropriate media specified. Note: ATCC quantitation results were obtained using 100 mm diameter Tryptic Soy Agar plates with 1.5% Tryptone, 0.5% Soytone, 0.5% Sodium Chloride, and 1.5% Agar (plates were procured from Teknova).
5. Use a sterile plate spreader (not provided) to spread the sample evenly over the surface of the media plate. Incubate the plates in an incubator at the appropriate temperature (see the product page or product sheet for details). If following a specific pharmacopeial chapter, use the incubation conditions specified.
6. Following incubation, count the number of visible colonies on the plate using a standard colony counting procedure.

To learn more about MicroQuant™, visit us online at www.atcc.org/MicroQuant. For assistance, please contact ATCC customer support at tech@atcc.org or your local distributor.

 10801 University Boulevard
Manassas, Virginia 20110-2209

 703.365.2700

 703.365.2701

 sales@atcc.org

 www.atcc.org

LCFU-112025-v05

©2025 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. Teknova is a registered trademark of Alpha Teknova, 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.