

Certificate of Quality Assurance

Bacterial Filtration Efficiency

Sartorius Safetyspace™ Filter tips protect your work from contamination. The products have been tested by an accredited independent laboratory for bacterial filtration efficiency (BFE).

The average BFE for Safetyspace™ Filter tips was >99,9%.

This shows that in addition to the protective Safetyspace™ gap the pure polyethylene pipette tip filter protects from cross-contamination.

Bacterial Filtration Efficiency:

Two product variants (Order codes: 790301F and 791001F) and total of three production batches were tested using an increased challenge level in a test setup adapted from ASTM F2101: "Standard Test Method for Evaluating the Bacterial Filtration Efficiency (BFE) of Medical Face Mask Materials, Using a Biological Aerosol of Staphylococcus aureus". A suspension of S. aureus, ATCC #6538, was delivered to the pipette tips from the smaller orifice at a challenge level of $>10^4$ colony forming units. The challenge was aerosolized using a nebulizer and delivered at a flow rate of 0,5 liters per minute. The adapted method was chosen as there is no standard test method to determine the filtration efficiency of pipette tip filters.

These results are based on one-time analysis and are for guidance to the BFE of Sartorius Safetyspace™ Filter tips.

Signed:



Date: 14. May 2020

Full Name:

Seppo Riikonen

Title:

Quality manager, Pipettes and Dispensing