

Certificate of Purity

ATP-free pipette tips

Sartorius purity certified Optifit tips and Safetyspace™ filter tips have been tested by a luminometric assay for absence of adenosine triphosphate (ATP). The test result proves the products are suitable and ready for use in monitoring ATP, testing hygiene and cleanliness and other applications where ATP contamination from plasticware is detrimental to results.

ATP:

An optimized volume of pure and ATP-free water was pipetted up and down using maximal pipette tip capacity. Presence of ATP in the extraction was detected with a luminometric assay by measuring the emitted light from a chemiluminescence reaction. ATP free tips were tested to contain $<3.0 \times 10^{-12}$ mg/ul of ATP.

These results are based on a one-time analysis and are for guidance to the probable level of ATP in Sartorius tips

Signed:  Date: 15. June 2020

Full Name: Seppo Riikonen
Title: Quality manager, Pipettes and Dispensing