



Certificate of Purity

for

Refill Tip: 50 – 1000 µl
Cat. Number: 791002
Lot. Number: PR142652
Expiry Date: 2023 – 01
Manufacturing Date: 16.1.2018

This is to certify that the product described above has been manufactured in accordance with ISO 9001:2015, Certificate No.FI17/5101 and ISO 13485:2016, Certificate No.FI17/5103 quality management protocols and has passed all relevant Quality Assurance procedures related to this product.

The product is certified to be DNase, RNase and endotoxin free and is tested by an independent laboratory.

The expiry date of the product is 5 years from the manufacturing date and is stated on the product label of the packaging.

Endotoxins:

Sample preparation was carried out according to FDA guidelines for medical devices by extracting the pipette tips cut into pieces for 1 hour at 37°C with an optimized volume of pyrogen-free water. Endotoxins in the extractions were detected with Chromogenic kinetic method according to Ph.Eur. 2.6.14, Method D, which is based on development of colour after cleavage of a synthetic peptide-chromogen complex. Validated test result for endotoxin-free pipette tips is < 0.005 IU/ml (EU/ml).

RNase:

The tips were pipetted up and down in an optimized volume of pure water using maximal pipette tip capacity. RNase activity in the extraction was measured with fluorometric assay by detecting degradation of labelled RNase substrate. RNase-free pipette tips show no evidence of RNase activity in the assay with the detection level of $<3,125 \cdot 10^{-9}$ U/µl when RNase A was used as a standard.

DNase:

The tips were pipetted up and down in an optimized volume of pure water using maximal pipette tip capacity. DNase activity in the extraction was measured with fluorometric assay by detecting degradation of labelled DNase substrate. DNase-free pipette tips show no evidence of DNase activity in the assay with the detection level of $<6,25 \cdot 10^{-5}$ U/µl when DNase I was used as a standard.

23.2.2018



Seppo Riikonen

Director, Quality & Process Development