## **Finntip Filter**

Verification of protection against DNA aerosols and acid vapours

- For preventing of aerosol contamination of pipette and sample
- For protection of inner parts of the pipette
- Optimized for use with Finnpipettes
- Finntip Filter Micro up to Finntip Filter 1000 µl are guaranteed free of human DNA, DNase and RNase
- Ideal for PCR work
- Irradiated tip racks to guarantee sterility



Finntip Filter 200 µl sterile

These test proved that Finntip Filter is an effective barrier against DNA aerosols and acid vapours. This feature is especially important for DNA and microbiological applications where even a small amount of carryover can disturb the test and cause false results.

North America: USA/Canada +1 866 984 3766
Europe: Austria +43 1 801 40 0, Belgium +32 2 482 30 30,
Finland +358 9 329 100, France +33 2 2803 2000, Germany
national toll free 08001-536 376, Germany international +49 6184
90 6940, Italy +39 02 95059 1, Netherlands +31 76 571 4440,
Russia/CIS +7 095 225 11 15, Spain/Portugal +34 93 223 3154,
Switzerland +41 44 454 12 12, UK/Ireland +44 870 609 9203
Asia: China +86 21 6865 4588 or +86 10 5850 3588, India +91
22 5542 9494, Japan +81 45 453 9220, Other Asian countries
+852 2885 4613

**Countries not listed**: +49 6184 90 6940 or +33 2 2803 2000 **www.thermofisher.com** 

	Aim	Method	Result
DNA test	To study if Finntip Filter is an effective barrier against DNA aerosols.	DNA-solutions with concentrations of 20, 50 and 100 $\mu$ g/ $\mu$ l were pipetted using Finntip Filter tips. To detect if DNA aerosol had gone through the filter PCR was used.	No DNA was detected at the other side of the filter. This means that Finntip Filters prevented DNA aerosol coming through filter into the interior of the pipette.
Acid test	To study if Finntip Filter is an effective barrier against acid aerosols.	Different trifluoroacetic acid (TFA) solutions (10, 20, 30, 35, 50, 75 and 100 %) were pipetted using Finntip Filter tips with their maximum volumes. To detect if acid vapour had gone through the filter, pH paper was used.	Finntip Filter prevented acid vapours coming through the filter. For example FT Filter micro and FT Filter 100 prevented a 75 % and FT Filter 1000 prevented a 30 % TFA solution coming through.