

# Andromeda Capillary Remover AN-051001

## General Information

NanoTemper Technologies offers dedicated capillaries for use with Andromeda instruments. Up to 48 single capillaries are used per run. The Andromeda Capillary Remover allows the quick and easy removal of all capillaries at once while avoiding direct contact of the user with capillaries and sample.

### Handling

- To remove capillaries after an experiment, open the Andromeda capillary tray gate and remove the magnetic lid from the capillary tray. Place the remover on the capillary tray with the NanoTemper logo facing upward. Slide the remover **slowly** towards the instrument body to collect the capillaries. When you have reached the end of the tray, tilt the open end of the remover upwards to collect the capillaries in the remover's reservoir.
- Clean the capillary remover regularly using water or ethanol.

**CAUTION** Danger of injury by broken glass. If sample has leaked from the capillaries during the experiment, capillaries may stick to the capillary tray. Remove capillaries slowly and with care to avoid breaking.

### General safety considerations



Be careful when using capillaries.

Broken capillaries are broken glass and can be harmful to skin and eyes.

For more information on capillary handling, please refer to the Andromeda instrument user manual or the capillary kit product insert.

### Help

For further information, please refer to our Andromeda User Manual or contact our NanoTemper application support through our Explorer Community:

<https://nanotempertech.force.com/explore/>