

Prometheus NT.48 Capillary Remover

General Information

NanoTemper Technologies offers dedicated capillaries for use with Prometheus series instruments. Up to 48 single capillaries are used per run. The Prometheus NT.48 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 Prometheus NT.48 capillary tray gate and remove the magnetic lid from the capillary tray. Place the remover on the capillary tray with the NanoTemper Technologies 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 Capillary Chips.
Broken capillaries are harmful to skin and eyes.

For more information on capillary handling, please refer to the Prometheus Instruments User Manual or the capillary kit product inserts.

Help:

For further information, please refer to the Prometheus Instruments User Manual or contact NanoTemper Technologies application support:

Call: +49-(0)89-4522895-0
Email: info@nanotempertech.com

V02_2018-12-10