

Prometheus - Capillary Chips

General Information

NanoTemper Technologies offers special Capillary Chips with high quality and verified functionality. Our Capillary Chips are suited for use in Prometheus NT.Plex and Prometheus Panta. It is highly recommended to use original Prometheus Capillary Chips for qualitative and reproducible measurements. The Prometheus warranty is only valid if the instrument is used with original Prometheus Capillary Chips.

Content

- 8 Capillary Chips (24 capillaries each)

Handling and Storage

- Store at room temperature
- Store in a dark and dry place
- Do not touch the Capillary Chips in the capillary center. Hold them from the ends
- Use gloves
- Capillary Chips are for single use
- Used Capillary Chips must be disposed of in accordance with federal, state and local environmental control regulations
- **Important:** Always close Capillary Chip box to keep Capillary Chips dust-free

Security Advice



Be careful when using Capillary Chips.

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

Capillary Type

Cat #	Capillary Type
PR-AC002	Standard
PR-AC006	High Sensitivity

Help

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

Call: +49-(0)89-4522895-0

Email: info@nanotempertech.com