

Andromeda – Capillary Kits

General Information

NanoTemper Technologies offers special Capillaries with high quality and verified functionality. Our Capillaries are developed for usage in Andromeda. It is highly recommended to use original Andromeda Capillaries for qualitative and reproducible measurements. The Andromeda warranty is only valid if the instrument is used with Andromeda Capillaries.

Content

- 1 vial of Capillaries (200 count)

Handling and Storage

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

Security Advice



Be careful when using Capillaries.
Broken capillaries are broken glass and can be harmful to skin and eyes.

Capillary Type

Cat #	Capillary Type
AN-041001	Andromeda Capillaries (200 count)

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/>