

# Andromeda – Capillary Chips Kits

## General Information

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

### Content

- 16 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 only
- 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
AN-042001	Andromeda Capillary Chips (16 chips of 24 capillaries each)

### 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/>