

## Tycho™ NT.6 Capillaries

Specially-designed and precisely engineered, Tycho NT.6 Capillaries are developed exclusively for use with the Tycho NT.6 system. These single-use capillaries enable researchers to analyze samples using small, microliter volumes. Manufactured in a state-of-the-art facility following stringent protocols and rigorous testing, Tycho NT.6 Capillaries ensure optimal performance, maximum reproducibility and the highest quality results.

### Content

Cat#	Name	Content	Store at
TY-C001	Tycho NT.6 Capillaries	200 Count	Store dry and in a dark place at room temperature

### Capillary handling and safety information

Keep vial closed to ensure capillaries are clean and dust-free. Hold individual capillaries at one end, not in the center. Use gloves. Dispose of capillaries according to applicable regulations and according to the sample contained in them.



**WARNING** Broken glass can cut skin. Handle capillaries carefully to avoid breaking. Do not touch broken capillaries.

### Packaging disposal

All packaging, including the cardboard insert with foil, can be recycled with paper products.

### Contact

For further information, visit [nanotempertech.com](http://nanotempertech.com) or email us at [support@nanotempertech.com](mailto:support@nanotempertech.com).

