

Monolith Capillary Chips

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

Content

Name	Store at
Monolith Premium Capillary Chips (16 chips with 24 capillaries each)	Store dry and in a dark place at room temperature

Capillary chip handling and safety information

To be used with the Monolith Premium Capillary Chip Lid (CAT# MO-T022).
Keep chip container closed to ensure capillaries are clean and dust-free. Hold individual capillary chips on both ends, not in the center. Use gloves. Dispose of capillary chips according to applicable regulations and according to the sample contained in them.



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

Packaging disposal

Dispose of the plastic capillary chip container according to applicable regulations. The cardboard can be recycled with paper products.

Contact

For further information, visit nanotempertech.com/explore

