

Monolith NT.LabelFree Capillary Kit

Specially-designed and precisely engineered, Monolith NT.LabelFree Capillaries are developed exclusively for use with the Monolith NT.LabelFree 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, Monolith NT.LabelFree Capillaries ensure optimal performance, maximum reproducibility and the highest quality results.

Content

Name	Store at
Monolith NT.LabelFree Capillaries	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/explore

