

Cat Nr: PR-P002

Prometheus® Capillary Sealing Applicators

Applicators for the sealing of capillaries for high temperature or long-term experiments for use in Prometheus®

CONTENT AND STORAGE

Prometheus Capillary Sealing Applicators are shipped at room temperature. Each kit contains two reusable applicators.

Expiry date: none

IMPORTANT INFORMATION BEFORE STARTING

This kit is designed to seal capillaries for experiments using Prometheus instruments. For standard thermal unfolding experiments (measurement times < 3 h, temperatures ≤ 95 °C), sealing of capillaries is not required. Longer measurements or higher temperatures require sealing of capillaries to prevent air-bubble formation and/or extensive evaporation.

Two sealing paste applicators are included with the Prometheus Capillary Sealing Kit. Sealing paste (Cat. No. PR-P001) must be purchased separately to seal capillaries. Sealing will not influence your measurements in any way since the sealing paste is chemically inert, silicone-free, and water-insoluble.

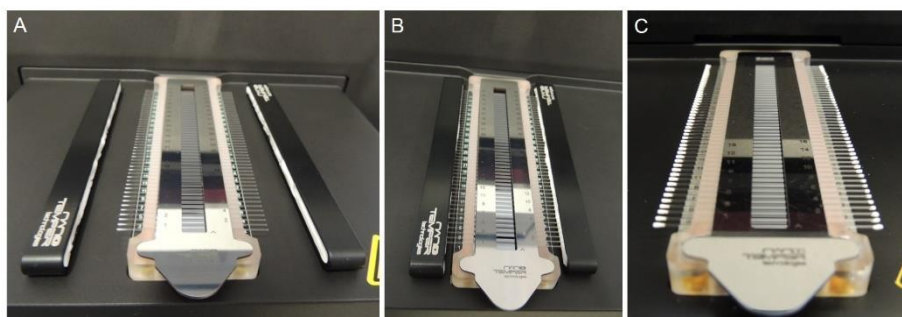
For safety instructions for handling of the capillary sealing paste, please consult the respective information sheet provided with the paste and the Safety Data Sheets (SDS). SDS are available from NanoTemper Technologies upon request.

CAPILLARY SEALING PROCEDURE

1. Fill capillaries with sample and place them on the capillary tray. When working with individual capillaries on the Prometheus NT.48 or Panta, make sure capillaries are aligned and fixed with the magnetic lid (see the Prometheus Series Manual for more information on sample loading).
2. Always wear protective gloves when handling the sealing paste. Apply sealing paste to the cavity of both sealing paste applicators.



3. (A) Place the applicators next to the capillary tray with the NanoTemper Technologies logos facing upward and the sealing paste toward the capillaries. (B) Slide the applicators toward the tray until all capillaries dip into the sealing paste. (C) Withdraw applicators.



4. Let the first layer dry for five minutes, then repeat. Always apply two layers of sealing paste. One loading of paste to the applicators is sufficient for the complete sealing process, as the sealing paste on the applicators will not dry within the application time.
5. Let the sealing paste dry another five minutes and remove the applicators from the Prometheus drawer before starting the experiment.
6. Execute the experiment using your Prometheus software.
7. When the sealing paste remaining in the applicators has dried (1-2 h), remove it and discard. The applicators can be reused.

NOTE

When the capillary sealing paste tube is opened after prolonged storage, the first few drops to emerge may be a colorless liquid. Discard this liquid and use only the white sealing paste for capillary sealing.

Contact

TECHNICAL SUPPORT

Please get in touch with us for specific questions concerning the product performance.

NanoTemper Technologies GmbH
Flößergasse 4
81369 München
Germany
Tel.: +49 (0)89 4522895 0
info@nanotempertech.com

PURCHASER NOTIFICATION

NanoTemper Technologies grants the buyer the non-transferable right to use the purchased product for research conducted by the buyer. The buyer cannot sell or otherwise transfer this product or its components for commercial purposes.

Limited warranty:

NanoTemper Technologies will replace any product that does not meet the specifications. This warranty limits NanoTemper Technologies' liability only to the cost of the product when within the expiration date. No warranty is applicable unless all product components are stored in accordance with instructions. NanoTemper Technologies assumes no responsibility or liability for any indirect or consequential loss or damage whatsoever. The above limited warranty is sole and exclusive.

For information contact NanoTemper Technologies GmbH
Flößergasse 4
81369 München
Germany
Tel.: +49 (0)89 4522895 0
info@nanotempertech.com

For ordering NanoTemper products:
Phone: +49 (0)89 4522895 0
Fax: +49 (0)89 4522895 60
E-mail: order@nanotempertech.com

NanoTemper®, Prometheus® and nanoDSF® are
registered trademarks.

V07_2021-03-24