

Cat Nr: PR-P001

Prometheus® Capillary Sealing Paste

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

CONTENT AND STORAGE

Prometheus Capillary Sealing Paste is shipped at room temperature. Each kit contains material sufficient for approximately 50 sealing procedures.

5 x 15 g Capillary Sealing Paste

Expiry date: see kit cover.

Capillary sealing paste can be used for at least three months after opening the tube.

SAFETY INSTRUCTIONS

Prometheus Capillary Sealing Paste is an oil-based liquid rubber. Keep away from skin and eyes. Do not eat. Easily flammable. Wear protective gloves when handling the sealing paste.

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.

Prometheus Capillary Sealing Applicators (PR-P002, sold separately) are required for capillary sealing.

The sealing paste contained in this kit will not influence your measurements in any way since it is chemically inert, silicone-free, and water-insoluble.

SAFETY INFORMATION

CAPILLARY SEALING PASTE

Hazard statements



H225	Highly flammable liquid and vapour.
H304	May be fatal if swallowed and enters airways.
H315	Causes skin irritation.
H319	Causes serious eye irritation.
H336	May cause drowsiness or dizziness.
H361f	Suspected of damaging fertility.
H373	May cause damage to organs through prolonged or repeated exposure.
H412	Harmful to aquatic life with long lasting effects.



Precautionary statements



P210	Keep away from heat, hot surfaces, sparks, open flames and other ignition sources. No smoking.
P235	Keep cool.
P260	Do not breathe dust/ fume/ gas/ mist/ vapours/ spray.
P264	Wash hands thoroughly after handling.
P271	Use only in well-ventilated area.
P273+P501	Avoid release to the environment. Dispose of content/container to an appropriate recycling and disposal facility.
P280	Wear protective gloves/ protective clothing/ eye protection/ face protection.
P301+P310	IF SWALLOWED: Immediately call a POISON CENTER/ doctor.
P331	Do NOT induce vomiting.

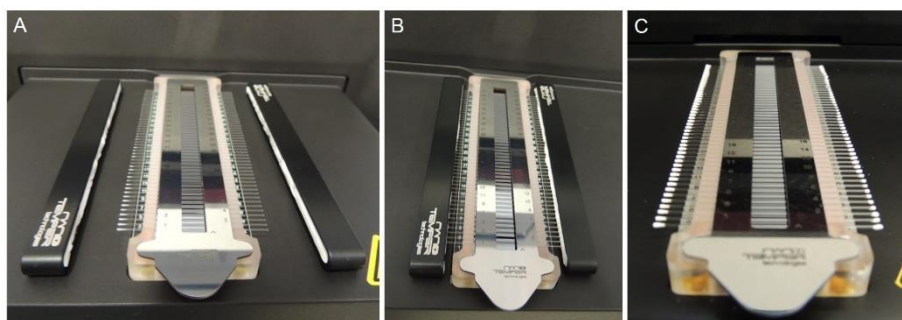
For more information, please consult the respective Safety Data Sheets (SDS). SDS are available from NanoTemper Technologies upon request.

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.

V06_2021-03-24