

Cat# PR-030001

Molecular Weight Reference Kit

For Prometheus Panta with Static Light Scattering Optics

CONTENTS 10 vials of PEG-40 solution (50 µL each)

STORAGE AND HANDLING

- Store at -20°C
- Store in a dark and dry place
- Wear gloves
- Vials are single use only

ADDITIONAL MATERIAL REQUIRED

- Prometheus Capillaries (PR-C002 or PR-C006)
- Capillary Chips (PR-AC002 or PR-AC006)

IMPORTANT INFORMATION BEFORE STARTING

Static light scattering (SLS) measures the average scattering intensity of a sample solution and can be used to calculate the average sample molecular weight. This reference sample ensures that measurements are accurate and should be run alongside with every molecular weight measurement.

PROCEDURE

1. Pull one vial from the freezer and allow it to fully thaw at room temperature.
2. Spin the vial for a few seconds to ensure that the sample is mixed well before use.
3. Open the single-use vial and fill 3 capillaries with the PEG-40 Solution.
4. Place the capillaries on the tray with your samples to be analyzed. Ensure that the same capillary type is used for all samples in the run.
5. Open the PR.Panta Control Software and assign the positions of the capillaries containing PEG-40 capillaries as the reference samples.
6. Discard of the vial once the measurement is complete.

Note: For more information on how to preform molecular weight analysis in PR.Panta Control, please refer to the software FAQ.

Help:

For information, contact our NanoTemper application support:

Call: +49-(0)89-4522895-0

Email: info@nanotempertech.com