

SpheroRuler

PROTOCOL



distributed in the US/Canada by:

EAGLE BIOSCIENCES, INC.

20A NW Blvd, Suite 112 Nashua, NH 03063

Phone: 617-419-2019 FAX: 617-419-1110

www.EagleBio.com info@eaglebio.com



EAGLE
BIOSCIENCES



Check videos
of protocol, examples of
results and much more on:
[idylle-labs.com/spheroruler-
by-fluoref](http://idylle-labs.com/spheroruler-by-fluoref)

idylle-labs.com
30, rue de Campo-Formio
Paris - France
+ 33 1 84 25 51 44
contact@idylle-labs.com



SPHR-P.16.11.22

idylle

1. Material you need

- The SpheroRuler suspension (7.10⁸ particles/mL in PBS 1X, pH 7.4, 0.02% azide)
- A blinking buffer

2. Storage

- Keep the SpheroRuler suspension in the fridge (2-8°C) for 7 months.
- Keep it in the dark in its opaque bag.

❗ **Do not freeze.**

3. SpheroRuler mounting

- Thaw the SpheroRuler suspension at room temperature.
- Sonicate for 1-2 minutes and vortex the SpheroRuler 5 minutes before using it – for homogeneous dispersion and to avoid aggregates.

💡 **Check that it is well resuspended: you should obtain an opalescent suspension of SpheroRuler after the sonication and vortex step.**

- Add 5 µL of the resuspended SpheroRuler in a 22mm/500µL Willco dish.
- The SpheroRuler concentration can be adjusted according to your preferred bead density.

- Add the blinking buffer.
- Put the coverslip on the cavity filled with the blinking buffer without making any bubble.
- Apply the sealant.
- Wait for 10-15 minutes to let the SpheroRuler beads settle down.
- Proceed with image acquisition.

4. Image acquisition

- Find a field of view with isolated SpheroRuler beads.
- Do a gentle z-scan to focus on the larger diameter.
- Record reference images by standard epifluorescence and brightfield microscopy.
- Select a ROI and perform dSTORM acquisition.
- Apply dSTORM image processing to reconstruct the SpheroRuler beads images.

💡 **The SpheroRuler beads are coated with a high density of fluorophores. We recommend using a multi-emitter fitting approach, or any equivalent, to guarantee an efficient localization of individual blinking events.**

