

Research Use Only. Not for use in diagnostic procedures.

## Opal™ 520 Reagent Pack (FP1487001KT)

## Material Provided

- Opal 520 Reagent, 1 vial
- Dimethyl sulfoxide, 1 x 100 uL

Please note that Opal 520 Reagent is provided dry. DMSO has been included to reconstitute this dye. Please follow the protocol provided below.

## Product Information

<b>Protocol</b>	<ol style="list-style-type: none"><li>1. Dissolve the Opal 520 Reagent in 75 uL of DMSO</li><li>2. Carefully dispense DMSO along the sides of the vial several times to dissolve any Opal Reagent that might coat the sides of the vial. <i>Minimize bubbles while mixing.</i></li></ol>
<b>Storage</b>	Store dry reagent in the dark at -20 °C. Upon reconstituting in DMSO, store in the dark at 4°C.
<b>Stability</b>	See label on outside of box for expiration date
<b>Safety Note</b>	DMSO is classified as hazardous and combustible. It is strongly recommended to wear disposable gloves and safety glasses while working with the items in this kit. Thorough washing of hands after handling is also recommended.
<b>Quality Control</b>	We certify that QC results of these reagents meet our quality release criteria.

### Opal Fluorophore Working Solution

Before each procedure, dilute Opal Fluorophore in 1X Amplification Diluent to make Opal Fluorophore Working Solution. The following dilution ratios are recommended:

- 1:100 for manual kits
- 1:150 for automated kits

Generally, 100-300 µL of Opal Working Solution is required per slide. Discard any unused portion of Opal Working Solution.

## Troubleshooting

### Technical Support Resources

- **Email:** [global.techsupport@perkinelmer.com](mailto:global.techsupport@perkinelmer.com)
- **Telephone**
  - **USA toll-free** **800-762-4000**
  - **EU toll-free** **00800 33 29 0000**
  - **China toll-free** **800 820 5046**
  - **Local contact numbers:** <http://www.perkinelmer.com/corporate/locations>

TSA and its use are protected under U.S. Patents 5,688,966, 5,863,748, 5,767,287, 6,372,937, 6,399,299, 6,593,100, 6,617,125 and patents pending, and foreign equivalents thereof.

For Research use only. This product is distributed and sold for research purposes only by the end-user in the research market, and, to that extent, by purchasing this product the end-user is granted a limited license to use this product for research use only. This product is not intended for diagnostic or therapeutic use and no license or right is granted for use of this product for diagnostic or therapeutic purposes. Purchase does not include or carry any right or license to use, develop or otherwise exploit this product commercially. Any commercial use, development or exploitation of this product without the express prior written authorization of PerkinElmer is strictly prohibited and may constitute infringement of the intellectual property rights of PerkinElmer under the aforementioned patents.

TSA is a registered trademark of PerkinElmer. Opal is a trademark of PerkinElmer. Other trademarks are property of their respective owners