

Research Use Only. Not for use in diagnostic procedures.

Opal™ 4-Color anti-Rabbit Manual IHC Kit 50 slides

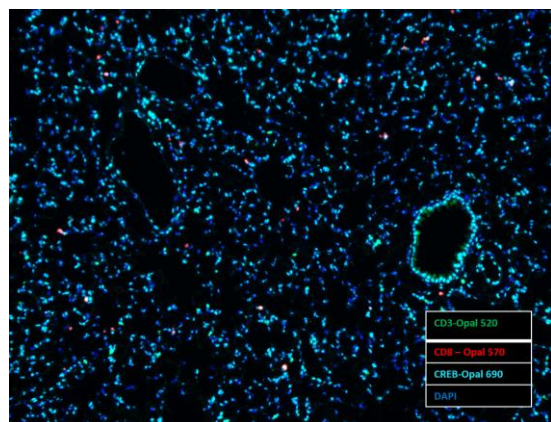
Product Information

| | |
|------------------------|--|
| Storage | Store dry Opal reagent at -20 °C. Upon reconstituting in DMSO, store at 2–8°C. Store remaining kit components at 2–8°C |
| Stability | This product is stable for a minimum of 3 months when stored at 2–8 °C |
| Application | The Opal 4-Color anti-Rabbit Manual IHC Kit is intended for multiplex fluorescent IHC. |
| Safety Note | DMSO is classified as hazardous and combustible. Some reagents in this kit contain Proclin® 300 that is classified as corrosive to metals and skin, a skin and eye irritant, and hazardous to the aquatic environment. DAPI is considered corrosive to the skin and an irritant to the eye. All other reagents are classified as nonhazardous. 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. |
| Quality Control | We certify that QC results of these reagents meet our quality release criteria. |
| Slide Number | When using this kit's recommended Opal dilution of 1:100, it will enable a 4-color assay on 50 slides. |

What is the Opal Method?

The Opal workflow allows simultaneous detection of multiple biomarkers in tissue. This Opal protocol was written specifically for a 4 or 7-Color IHC in formalin-fixed paraffin-embedded (FFPE) tissue. The approach involves detection with Opal fluorophores, followed by microwave treatment (MWT) for: removal of primary and secondary antibodies¹; removal of any non-specific staining; and reduction of tissue auto-fluorescence. The Opal signal is largely unaffected by MWT and antibody removal. After MWT, another round of staining can be performed for additional target detection without risk of antibody cross reactivity.

Opal allows staining of multiple IHC targets using unlabeled primary antibodies raised in the same species². Combining Opal with multispectral imaging and analysis enables simultaneous, quantitative results for up to 6 biomarkers in fluorescence, even with co-localized markers, plus nuclear counterstain (DAPI). **Fluorescent multispectral imaging (usually with the Mantra™ or Vectra® systems) is required for successful analysis of more than 3 Opal fluorophores at once.**



PerkinElmer provides assistance with assay development and offers [multiplex Opal IHC and IF services](http://www.perkinelmer.com/Opal). Visit: www.perkinelmer.com/Opal.

Material Provided

| | Format* | Catalog # | Kit Components |
|---|-----------|-------------|--|
| Opal 4-Color anti-Rabbit Manual IHC Kit | 50 slides | NEL840001KT | <ul style="list-style-type: none"> • 1X Plus Amplification Diluent (1 X 50 mL) • Opal 520 Fluorophore • Opal 570 Fluorophore • Opal 690 Fluorophore • DMSO (1 X 500 µL) • Spectral DAPI solution (1 X 1.5 mL) • Opal Polymer anti-Rabbit HRP (1 X 10mL) • Opal Polymer anti-Rabbit HRP Diluent (1x 40mL) • Blocking/Ab Diluent (1 X 100 mL) • 10X AR6 buffer (2 X 250ml) |

*The format of the kit is based on ~150 µL per slide of Opal Working Solution.

Reagents and Materials

Required Materials

- Baths and solvents for deparaffinization and rehydration of FFPE tissue. Xylene is recommended for deparaffinization. Histological grade ethanol is required for rehydration.
- Standard microwave oven with carousel, rated at 1,000 W or higher with 10 or more power settings
- Standard staining dishes
- Opal slide processing jars (Perkin Elmer catalogue number STJAR4) or equivalent
- Slide incubation/humidity tray
- Hydrophobic barrier pen
- Glass coverslips (No. 1.5)
- Control tissues
- Charged slides

Required Reagents

- TBST wash buffer
- Neutral buffered formalin (NBF)
- Peroxidase-free water.
Note: This specification may be met by commercial “cell culture grade” water or ultra-pure (i.e. Milli-Q™) water.
- Antibody diluent & blocking reagent
 - Antibody Diluent / Block (PerkinElmer catalog number ARD1001EA) is recommended
 - Mouse tissue often requires more robust blocking methods, such as goat or horse serum.
 - Other options should be validated independently
- Primary antibodies for targets of interest
- Mounting medium for fluorescence (i.e. ProLong® Diamond Antifade Mountant (Thermofisher)).
- AR9 Buffer (PerkinElmer catalog number AR900250ML) may be required for certain antigens requiring a higher pH antigen retrieval buffer

Solutions to prepare

TBST Wash Buffer

25 mM TRIS-HCl, pH 7.5
150 mM NaCl
0.05% Tween®20 (v/v)

AR6 Buffer Working Solution:

Dilute 10X AR6 buffer at 1:10 with peroxidase-free water.

Antibody Diluent

Antibody Diluent / Block from PerkinElmer is supplied as a ready-to-use solution.

Primary Antibody Working Solution

Dilute primary antibody in PerkinElmer Antibody Diluent / Block at optimal concentration for Opal detection as determined below.

Secondary Antibody Working Solution

Opal Polymer anti-Rabbit HRP is supplied as a concentrate. Working dilution recommendation for manual assays is 1:5 in the supplied Opal Polymer anti-Rabbit HRP Diluent.

Opal Reagent Stock Solution

Dissolve the Opal Reagents in 75 μ L of DMSO. Carefully dispense DMSO along the sides of the vial several times to dissolve any Opal Reagent that might coat the sides of the vial. (Minimize bubbles while mixing.)

Opal Fluorophore Working Solution

Before each procedure, dilute Opal Fluorophore in 1X Amplification Diluent to make Opal Fluorophore Working Solution. Recommend to start diluting the Opal fluorophore at 1:100. Optimize your assay according on Opal Assay Development Guide. Generally, 100-300 μ L of Opal Working Solution is required per slide. Discard any unused portion of Opal Working Solution.

DAPI working solution

Add two drops of DAPI solution into 1ml of TBST. Approximately 150 μ L of DAPI Working Solution is required per slide. Discard any unused portion of DAPI Working Solution.

Recommendations

- Use xylene for removal of paraffin from FFPE tissue sections. Do not let slides dry out between steps.
- Before the first use, spin down the Opal Fluorophore tubes with a standard microcentrifuge to make sure that all of the solution is at the bottom of the tubes.
- A humidified chamber is recommended for all incubation steps.
- Drain off as much of the incubation solutions as possible before addition of the next solution, to prevent reagent dilution and uneven staining. Blot area around, but not on, tissue section using absorbent paper.
- Be sure to use enough volume of each reagent to completely cover sections or cells.
- Optimize your monoplex staining according to the Opal Assay Development Guide, with all MWT steps incorporated.
- Spectral DAPI in this kit is formulated for optimal separation from other fluorophores. Exposure time may be somewhat longer than other DAPI formulations.
- Before attempting multiplexed staining, assay conditions for each analyte should be optimized singly with Opal detection (refer to Opal Assay Development Guide for more details).
- Microwave treatment (MWT) as outlined in this protocol performs antigen retrieval, quenches endogenous peroxidases, and removes antibodies from earlier staining procedures.
- This protocol was developed with specified reagents. Other options should be independently validated.
- If the antigen requires higher pH retrieval, it is recommended to purchase PerkinElmer's AR9 Buffer.

Opal Optimization Strategies

Microwave Optimization

Microwave treatment (MWT) is used in the Opal method to quench endogenous peroxidase activity, for antigen retrieval, and to remove antibodies after a target has been detected. Timing for each step in the procedure may have to be modified for the microwave oven that you are using. Slides are placed vertically in an Opal Slide Processing Jar which is then filled to the top (~140 mL) with AR6 or AR9 buffer and covered loosely. One jar is placed in the microwave at a time, near the edge of the carousel to ensure even distribution of energy. The microwave procedure consists of two steps:

1. 100% power until the boiling point is reached. The time for this step may have to be increased or decreased depending upon the performance of the microwave in your lab. This will usually take 45-90 seconds.
2. 20% power for 15 minutes.

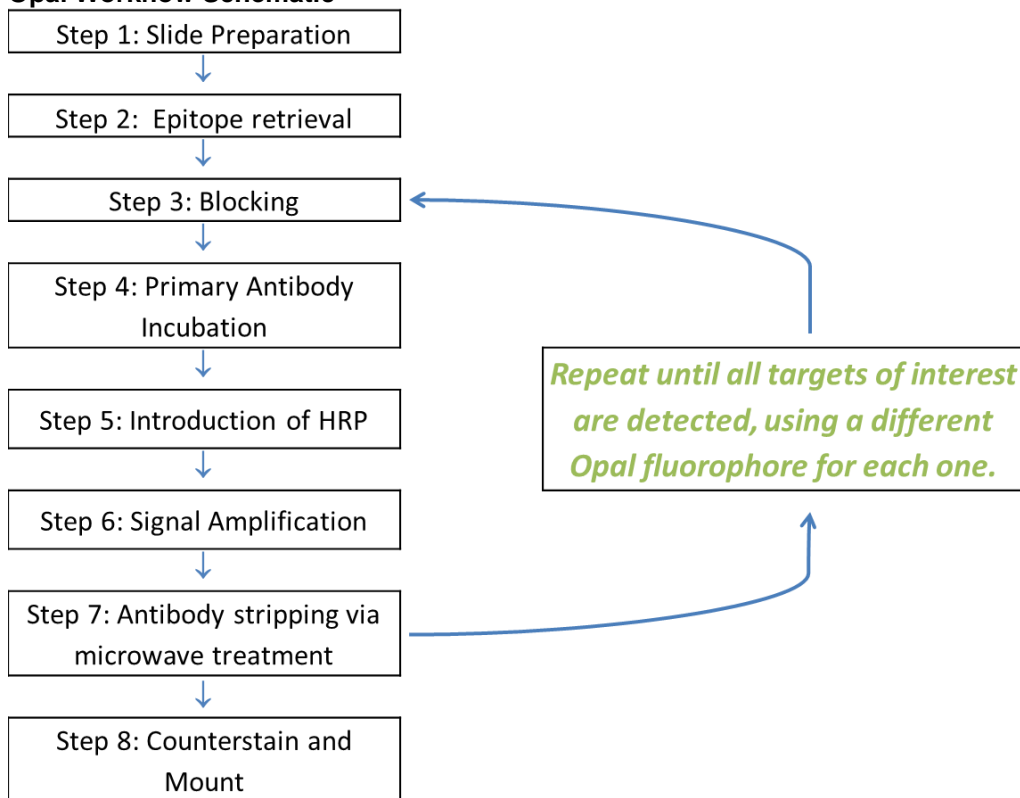
Do not operate the microwave unattended and keep the oven chamber clean and clear of debris.

Opal Multiplexed IHC

Single analyte Opal IHC methods may be combined for multiplexed detection within a single tissue section. After signal amplification, MWT is performed to strip away detection antibodies. The Opal fluorophore is largely unaffected by MWT because it is covalently bound. Then the process is repeated using another Opal fluorophore.

Importantly, a different Opal fluorophore should be used for each target.

Opal Workflow Schematic



Step by Step Opal IHC Protocol (Single Analyte) Refer to Opal Assay Development Guide for more detailed instruction:

Single analyte Opal IHC assays should be optimized before combining for use in multiplexed detection. Concentration for each primary antibody should be optimized with the selected Opal fluorophore to yield intensity between 10 to 30 counts. Optimized single fluorophore images (without DAPI counterstain) will subsequently be used for spectral library development.

The following protocol details the workflow for a single analyte, and can subsequently be employed for multiplexed IHC. In multiplexed IHC, the order of target/fluorophore detection may be a point of optimization, and must be independently validated.

Step 1: Slide Preparation

Prepare tissues or cells for detection with Opal using standard fixation and embedding techniques. We recommend running an isotype control slide with all steps replacing the primary antibody with corresponding isotype control for each experiment. For each slide, baked in the oven at 65°C for 1 hour; dewax with xylene (3 x 10 min) and rehydrate through a graded series of ethanol solutions: (100% 1 x 10 min; 95% 1 x 10 min; and rinse in 70%). After rehydration, briefly rinse slides in distilled water and fix in 10% neutral buffered formalin for 20min. Longer times of fixation in NBF may be needed for certain tissues such as skin.

Rinse slides in distilled water and then in the appropriate AR buffer.

Step 2: Microwave treatment

Place slides in an Opal Slide Processing Jar and fill it completely with the appropriate AR buffer. Loosely cover the jar with lid, place it in microwave for 45 sec at 100% power; (may require optimization as described). Microwave for an additional 15 min at 20% power. Allow slides to cool down at room temperature before proceeding (15 – 30 min). Importantly, do not let slides dry out. Rinse slides in distilled water followed by TBST.

Step 3: Blocking

Use a hydrophobic barrier pen to completely surround the tissue section on the slide. Cover tissue sections with blocking buffer and incubate slides in a humidified chamber for 10 min at room temperature.

- *Note: This protocol was developed with PerkinElmer Antibody Diluent / Block for blocking. Other options should be independently validated.*

Step 4: Primary Antibody Incubation

Drain off the blocking buffer and apply Primary Antibody Working Solution. Incubate according to the manufacturer's instructions regarding incubation time and temperature requirements or conditions optimized within your lab. Use enough volume to completely cover the tissue section (generally 100-300 µL per slide).

Rinse slides in TBST. Wash the slides 3 x 2 min in TBST at room temperature, preferably with agitation.

Step 5: Introduction of Secondary-HRP

Incubate slides in the working solution Opal Polymer anti-Rabbit HRP for 10 min at room temperature. Use adequate reagent volume to cover the tissue section, generally 100-300 µL per slide.

- *Note: Opal Polymer anti-Rabbit HRP working concentration is recommended at 1:5 with the provided Opal Polymer Diluent. Other concentrations need to be validated.*

Rinse slides in TBST. Wash the slides 3 x 2 min in TBST at room temperature preferably with agitation.

Step 6: Opal Signal Generation

Drain off excess wash buffer and pipette 100-300 µL of Opal Fluorophore Working Solution onto each slide. Incubate the slides at room temperature for 10 mins.

Rinse slides in TBST. Wash the slides 3 x 2 min in TBST at room temperature preferably with agitation.

Rinse slides in the appropriate AR buffer.

Step 7: Microwave treatment

Place slides in an Opal Slide Processing Jar and fill it completely with the appropriate AR buffer. Loosely cover the jar with lid, place it in microwave for 45 sec at 100% power; (may require optimization as described). Microwave for an additional 15 mins at 20% power. Allow slides to cool down at room temperature before proceeding (15 – 30 min). Importantly, do not let slides dry out. Rinse slides in distilled water followed by TBST.

This microwave step strips the primary-secondary-HRP complex allowing introduction of the next primary antibody. For detection of the next target, restart the protocol at Step 3: Blocking.

If all targets have been detected, continue to Step 8.

Step 8: Counterstain and Mount

Apply DAPI Working Solution for 5 min at room temperature in a humidity chamber. Wash the slides for 2 min in TBST buffer and then for 2 min in water. Coverslip slides with mounting medium (i.e. ProLong® Diamond Antifade Mountant (ThermoFisher)). *(Note: do not counterstain monoplex slides to be used for spectral library development.)*

Imaging and Analysis

Visualization of 4- or 7-color Opal slides can be performed using Mantra or Vectra Quantitative Pathology Imaging Systems. The systems use multispectral imaging for quantitative unmixing of many fluorophores and tissue autofluorescence, enabling advanced analysis including *in situ* cellular phenotyping. For more information, please see: <http://www.perkinelmer.com/quantitative-pathology>.

Important Notes:

1. All standard Vectra or Mantra epi-fluorescent cubes should be used for imaging Opal slides: DAPI, FITC, CY3, Texas Red, and CY5.
2. If the Opal fluorophores are not found in the Stain Store Manager in your version of inForm, please visit <http://www.perkinelmer.com/resources/software-downloads.xhtml> to download the latest inForm update that contains the novel Opal stains. If you believe that you have the latest version of inForm and you still cannot find the stains in the store manager, please contact your Field Application Specialist.

References

- ¹ Toth, Zsuzsanna E., and Eva Mezey. "Simultaneous visualization of multiple antigens with tyramide signal amplification using antibodies from the same species." *Journal of Histochemistry & Cytochemistry* 55.6 (2007): 545-554
- ² Stack, E.C., Wang, C., Roman, K., and Hoyt, C.C. "Multiplexed immunohistochemistry, imaging, and quantitation: a review, with an assessment of Tyramide signal amplification, multispectral imaging and multiplex analysis." *Methods*: (2014) doi:10.1016/j.ymeth.2014.08.016.

Troubleshooting

Technical Support Resources

- **Email:** global.techsupport@perkinelmer.com
- **Telephone**
 - **USA toll-free** **800-762-4000**
 - **Worldwide** **+1 203-925-4602**
 - **Fax** **+1 203-944-4904**
 - **Local contact numbers:** <http://www.perkinelmer.com/corporate/locations>

IHC Troubleshooting

| PROBLEM | REMEDY |
|-----------------|---|
| Low Signal | <ul style="list-style-type: none"> • Titer primary and Opal dye to determine optimum concentration for Opal detection. • Increase primary antibody and Opal Working Solution incubation time. (Suggested range is 3-10 minutes.) • Use additional rounds of microwave treatment to unmask the target. |
| Excess Signal | <ul style="list-style-type: none"> • Decrease concentration of primary antibody. • Decrease concentration of Opal dyes • Decrease Opal Working Solution incubation time. (Suggested range is 3-10 minutes.) |
| High Background | <ul style="list-style-type: none"> • Confirm that microwave treatment step has fully quenched endogenous peroxidases. • Titer primary and/or Opal dyes to determine optimum concentration for Opal detection. • Freshly prepare buffers. • Evaluate laboratory water source for contamination with HRP. • Check for endogenous biotin (if using streptavidin conjugates) • Increase number and/or length of washes. |

Opal Fluorophore Excitation and Emission Maxima

| Fluorophore | Opal Multicolor IHC Kits | | Excitation | Emission | Cap color |
|---------------|--------------------------|---------|------------|----------|-----------|
| | 4-color | 7-color | | | |
| Spectral DAPI | ✓ | ✓ | 358nm | 461 nm | Blue |
| Opal520 | ✓ | ✓ | 494 nm | 525nm | Green |
| Opal540 | | ✓ | 523nm | 536nm | Yellow |
| Opal570 | ✓ | ✓ | 550 nm | 570 nm | Red |
| Opal620 | | ✓ | 588 nm | 616 nm | Amber |
| Opal650 | | ✓ | 627 nm | 650 nm | Orange |
| Opal690 | ✓ | ✓ | 676nm | 694 nm | Clear |

Related Products

Opal Multiplex IHC Detection Kits

| | SIZES | PRODUCT NUMBER |
|--|-----------|----------------|
| Opal 4-Color Automation IHC Kit* | 50 slides | NEL820001KT |
| Opal 7-Color Automation IHC Kit* | 50 slides | NEL821001KT |
| Opal 4-Color Manual IHC Kit | 50 slides | NEL810001KT |
| Opal 7-Color Manual IHC Kit | 50 slides | NEL811001KT |
| Opal 4-Color anti-Rabbit Automation IHC Kit* 50 Slides | | NEL830001KT |
| Opal 4-Color anti-Rabbit Manual IHC Kit | 50 Slides | NEL840001KT |
| Opal 4 Lymphocyte Kit | 50 slides | OP4LY2001KT |
| Opal 7 Immunology Discovery Kit | 50 slides | OP7DS2001KT |
| Opal 7 Tumor Infiltrating Lymphocyte Kit | 50 slides | OP7TL3001KT |
| Opal 7 Solid Tumor Immunology Kit | 50 slides | OP7TL4001KT |

*Optimized for Leica Biosystems BOND RX System

Opal Reagent Packs

| | PRODUCT NUMBER |
|-----------------------|----------------|
| Opal 520 Reagent Pack | FP1487001KT |
| Opal 540 Reagent Pack | FP1494001KT |
| Opal 570 Reagent Pack | FP1488001KT |
| Opal 620 Reagent Pack | FP1495001KT |
| Opal 650 Reagent Pack | FP1496001KT |
| Opal 690 Reagent Pack | FP1497001KT |

Ancillary

| | PRODUCT NUMBER |
|--|----------------|
| 1X Plus Automation Amplification Diluent 1 X 50 mL | FP1609 |
| 1X Plus Amplification Diluent 1 x 50 mL | FP1498 |
| Opal Polymer anti-Rabbit HRP Kit | ARR1001KT |
| AR6 buffer (10X) 4 x 250 mL | AR6001KT |
| AR6 buffer (10X) 250 mL | AR600250ML |
| AR9 buffer (10X) 4 x 250 mL | AR9001KT |
| AR9 buffer (10X) 250 mL | AR900250ML |
| Antibody Diluent / Block 100 mL | ARD1001A |
| Opal Polymer HRP Ms + Rb 50 mL | ARH1001A |

For the latest product listings, please go to www.perkinelmer.com/opal.

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.

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.