

For Laboratory Use. A research method for research purposes only.

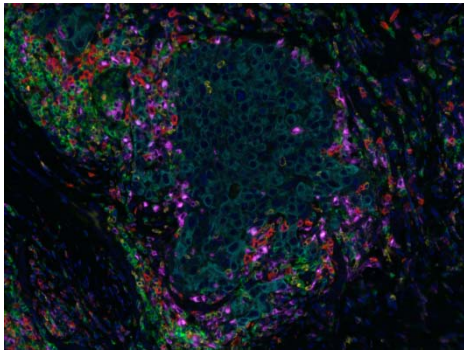
Opal™ 4-Color Automation IHC Kit 50 slides Opal™ 7-Color Automation IHC Kit 50 slides

Optimized for Leica Biosystems BOND RX System

Product Information

Storage	Store kits in the dark at 4°C.
Stability	See kit label on outside of box for expiration date
Application	The Opal 4 or 7-Color kits are intended for multiplex fluorescent IHC.
User Manual	Download the latest version of the User Manual at www.perkinelmer.com/opal
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.

What is the Opal Method?



Human breast cancer tissue was stained with CD4, CD8, CD20, FoxP3, CD68 and PanCK using Opal 7-color Automation Kit on BOND RX automated IHC & ISH stainer. Image was captured using Vectra 3.0 automated quantitative pathology imaging system and image analysis was performed using inForm software.

The Opal workflow allows simultaneous detection of multiple biomarkers in tissue. This Opal protocol was written specifically for 4 or 7-Color IHC on formalin fixed paraffin embedded (FFPE) tissue section*. The approach involves detection with Opal reactive fluorophores, followed by heat-induced stripping 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 stripping and antibody removal. After stripping step, 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 4 Opal fluorophores at once.**

***Please contact us if you would like to work with other types of samples. 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 Automation IHC Kit	50 slides	NEL800001KT	<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) • Blocking/Ab Diluent (1 X 100 mL) • Opal Polymer HRP Ms + Rb (1 x 50 mL)
Opal 7-Color Automation IHC Kit	50 slides	NEL801001KT	<ul style="list-style-type: none"> • 1X Plus Amplification Diluent (2 X 50 mL) • Opal 520 Fluorophore • Opal 540 Fluorophore • Opal 570 Fluorophore • Opal 620 Fluorophore • Opal 650 Fluorophore • Opal 690 Fluorophore • DMSO (1 X 500 µL) • Spectral DAPI solution (1 X 1.5 mL) • Blocking/Ab Diluent (1 X 100 mL) • Opal Polymer HRP Ms + Rb (2 x 50 mL)

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

Troubleshooting

Technical Support Resources

- Email: 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>

IHC Troubleshooting

PROBLEM	REMEDY
Low Signal	<ul style="list-style-type: none"> • Titer primary and/or Opal dyes to determine optimum concentration for Opal detection. • Increase primary antibody and/or Opal Working Solution incubation time. • Change staining order of multiplex and move this specific target detection towards the end of staining procedure..
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"> • Add extra step to block endogenous peroxidases. • Titer primary and/or Opal dyes to determine optimum concentration for Opal detection. • Make fresh buffers. • Evaluate laboratory water source for contamination. • 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	NEL800001KT
Opal 7-Color Automation IHC Kit*	50 slides	NEL801001KT
Opal 4-Color IHC Kit	50 slides	NEL794001KT
Opal 5-color fIHC Kit	50 slides	NEL795001KT
Opal 6-color fIHC Kit	50 slides	NEL796001KT
Opal 7-color fIHC Kit	50 slides	NEL797001KT
Opal 4 Lymphocyte Kit	50 slides	OP4LY1001KT
Opal 7 Immunology Discovery Kit	50 slides	OP7DS1001KT
Opal 7 Tumor Infiltrating Lymphocyte Kit	50 slides	OP7TL1001KT
Opal 7 Solid Tumor Immunology Kit	50 slides	OP7TL2001KT

*Optimized for Leica Biosystems BOND RX System

Opal Reagents

	SIZES	PRODUCT NUMBER
Opal520 Reagent, powder	50 slides	FP1487A
Opal540 Reagent, powder	50 slides	FP1494A
Opal570 Reagent, powder	50 slides	FP1488A
Opal620 Reagent, powder	50 slides	FP1495A
Opal650 Reagent, powder	50 slides	FP1496A
Opal690 Reagent, powder	50 slides	FP1497A
Spectra DAPI	50 slides	FP1490A

Ancillary

	PRODUCT NUMBER
1X Plus Amplification Diluent 1 X 50 mL	FP1498A
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.

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.