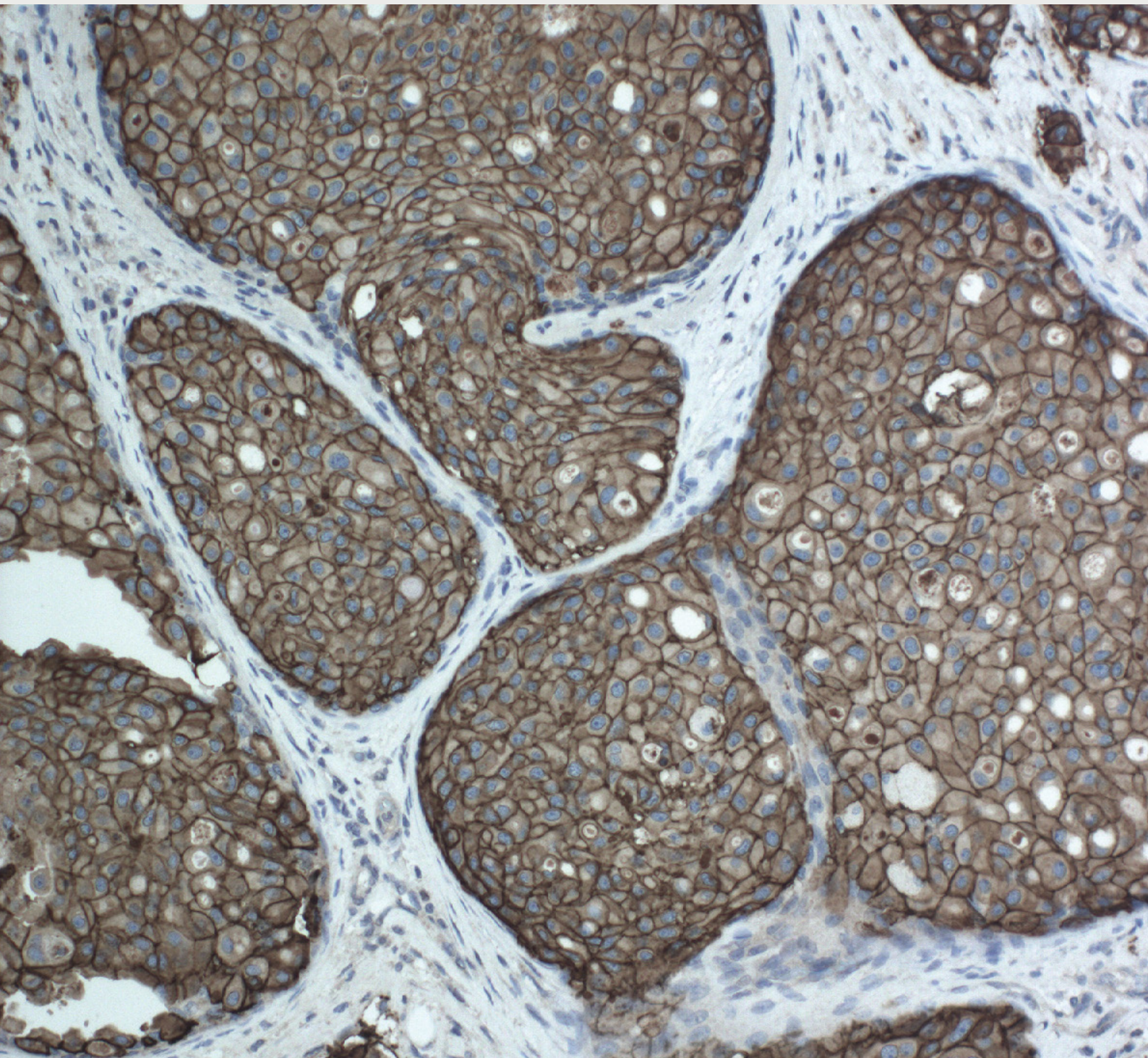




# Immunohistochemistry

Complete Solutions for IHC Analysis



[enzolifesciences.com](http://enzolifesciences.com)



# ELIMINATE DATA AMBIGUITY

**High Sensitivity. Low Background. True Results.**

For decades, immunohistochemistry (IHC) has been an invaluable tool for the detection, localization, and quantification of antigens in preserved tissue for research and diagnostic purposes. Enzo's complete set of tools improve the quality and efficiency of your IHC process from start to finish. Avoid ambiguity with high-sensitivity and low background reagents that are compatible with manual and automated platforms. Combined with the market's most extensive palette of unique chromogens, Enzo provides the flexibility needed for simple or complex protocols.

*"The POLYVIEW® PLUS HRP-DAB (Anti-Rabbit) kit worked very well; there was no background and the protocol is faster than many other IHC kits available commercially. I would surely recommend it. I was really happy with the POLYVIEW® PLUS HRP-DAB (Anti-Mouse) kit too."*

– Columbia University

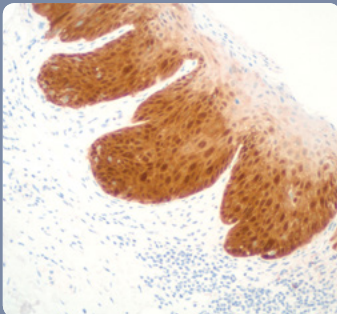
*"We tested Enzo's POLYVIEW® IHC detection reagents and HIGHDEF® chromogens on our immunohistochemical autostainers against similar available detection reagents and chromogens. We found that Enzo's detection reagents and chromogens matched or exceeded what we currently used, allowing for increased sensitivity and specificity when used against both human and murine tissue."*

– The Beatson Institute

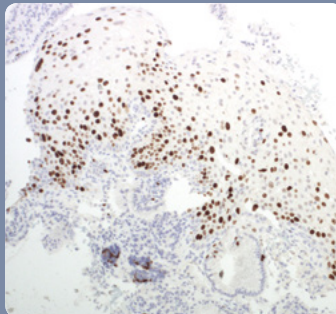
## ENZO'S IHC PLATFORM

- Eliminates false positive results
- Consistent, reliable, and sharp results
- Compatible with both manual and automated platforms
- Cost-effective and time-saving

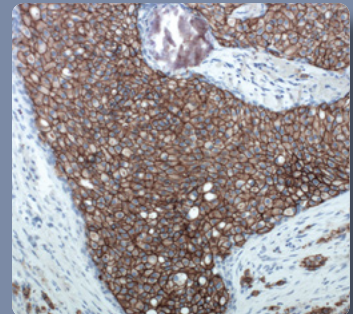
## IHC ANTIBODIES FOR CRITICAL MARKERS



**p16 Antibody**  
ENZ-ABS377



**Ki-67 Antibody**  
ENZ-ABS678



**HER2/Neu Antibody**  
ENZ-ABS381

# PROVIDING THE WHOLE ASSAY: RETRIEVAL TO VISUALIZATION

**Antigen  
Retrieval**

**Diluted  
Primary  
Antibody**

**Detection**

**Chromogens**

**Hematoxylin Counterstain**

**Mounting**

Category	Product Name	Product #
<b>Antigen Retrieval Reagents</b>	Antigen Retrieval Reagent, pH 6 (10X)	ENZ-ACC112
	Antigen Retrieval Reagent (Citrate, pH 7.0)	ADI-950-271
	Antigen Retrieval Reagent (Citrate, pH 8.0)	ADI-950-272
	Antigen Retrieval Reagent (EDTA, pH 8.0)	ADI-950-273
	Antigen Retrieval Reagent, pH 9 (10X)	ENZ-ACC113
	Antigen Retrieval Reagent (Tris, pH 10.0)	ADI-950-275
	IHC Enzyme Antigen Retrieval Reagent	ADI-950-280
<b>Diluted Primary Antibody</b>	Antibody Blocking Buffer	ENZ-ACC108
	IHC Diluent (Primary Ab)	ADI-950-244
	CD138 Monoclonal Antibody	ENZ-ABS683
	CD3 (Human) Monoclonal Antibody (TR66)	ALX-804-822
	Ki-67 Monoclonal Antibody (Ready-to-Use)	ENZ-ABS682
	p16 Monoclonal Antibody	ENZ-ABS377
	p53 Monoclonal Antibody (PAb122)	ADI-KAM-CC002
	Vimentin Monoclonal Antibody	ENZ-ABS694
<b>Detection</b>	MULTIVIEW® PLUS (mouse-HRP/rabbit-AP) IHC Kit	ENZ-KIT181
	MULTIVIEW® (mouse-HRP/mouse-AP) IHC Kit	ADI-950-101
	POLYVIEW® PLUS HRP (Anti-Rabbit) Reagent	ENZ-ACC103
	POLYVIEW® PLUS HRP (Anti-Mouse) Reagent	ENZ-ACC104
	POLYVIEW® PLUS AP (Anti-Rabbit) Reagent	ENZ-ACC110
	POLYVIEW® PLUS HRP-DAB (Anti-Rabbit) Kit	ENZ-KIT159
	POLYVIEW® PLUS HRP-DAB (Anti-Mouse) Kit	ENZ-KIT160
	SAVIEW® PLUS HRP Reagent	ENZ-ACC102
	SAVIEW® PLUS AP Reagent	ENZ-ACC111
<b>HRP Chromogens</b>	HIGHDEF® DAB Chromogen/Substrate Set	ENZ-ACC105
	HIGHDEF® Yellow IHC Chromogen (HRP)	ADI-950-170
	HIGHDEF® AEC (Red) IHC Chromogen (HRP)	ADI-950-210
	HIGHDEF® Blue IHC Chromogen (HRP)	ADI-950-151
	HIGHDEF® Black IHC Chromogen (HRP)	ADI-950-171
	HIGHDEF® Enhancer (DAB)	ADI-950-181
	HIGHDEF® Enhancer (DAB, plus)	ADI-950-182
<b>AP Chromogens</b>	HIGHDEF® Red IHC Chromogen (AP)	ADI-950-140/1
	HIGHDEF® Blue IHC Chromogen (AP)	ADI-950-150
	HIGHDEF® Green AP Chromogen Substitute	ENZ-ACC130
	HIGHDEF® Enhancer (AP)	ADI-950-180
<b>Counterstain</b>	HIGHDEF® Hematoxylin	ENZ-ACC106
<b>Mounting Medium</b>	HIGHDEF® Fluoromount	ADI-950-260
	HIGHDEF® Mount	ADI-950-261

## Accessories

IHC/ISH Peroxidase Block	ENZ-ACC107	IHC Diluent (AP)	ADI-950-241
IHC PAP (Mini) Pen	ADI-950-232	IHC Diluent (Peroxidase I)	ADI-950-242
IHC PAP Pen	ADI-950-233	IHC Diluent (Peroxidase II)	ADI-950-243
IHC Tissue Primer	ADI-950-234	IHC PBS Tablets	ADI-950-240
IHC Universal Negative Control Reagent	ADI-950-231	IHC Fast Red Salt Tablets Kit	ADI-950-191
IHC Wash Buffer	ADI-950-235		

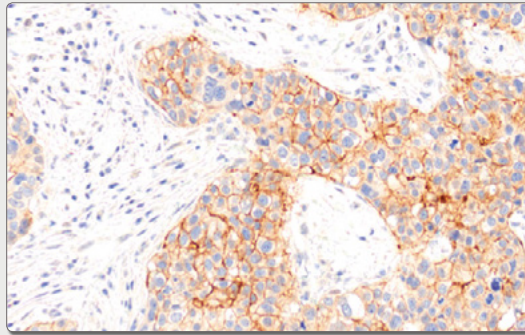
# DETECT WITH CONFIDENCE

## Enzo's POLYVIEW® PLUS Detection System Eliminates False Positive Results Observed with Common Commercially Available Kits

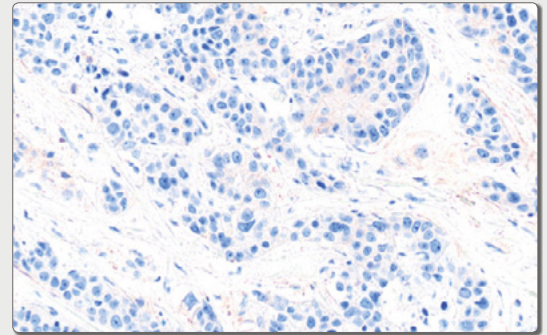
*Comparative analysis of Enzo with a leading market competitor, using samples with clinical relevance, demonstrates elimination of false positive results*

- Breast cancer tissue negative for HER2/Neu
- As expected, tissue stained with POLYVIEW® PLUS HRP Detection System is negative for HER2/Neu; the competitor HRP detection gave a false positive result
- HER2/Neu negative, confirmed by FISH analysis

**Competitor HRP Detection  
(HER2/Neu is false positive)**

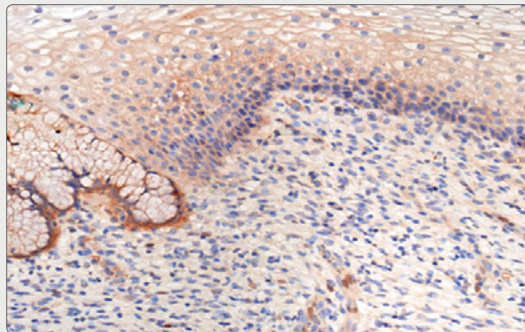


**Enzo's POLYVIEW® PLUS HRP Detection  
(HER2/Neu is negative)**

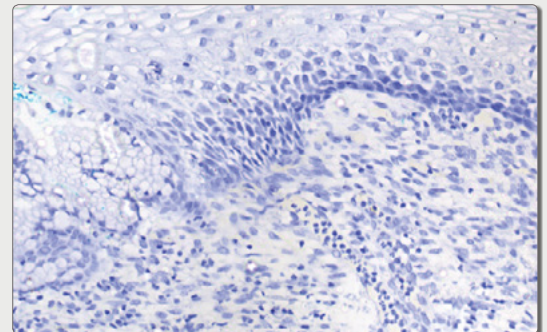


- Cervical cancer tissue negative for p16
- As expected, tissue stained with POLYVIEW® PLUS HRP Detection System is negative for p16; the competitor HRP detection gave a false positive result

**Competitor HRP Detection  
(p16 is false positive)**



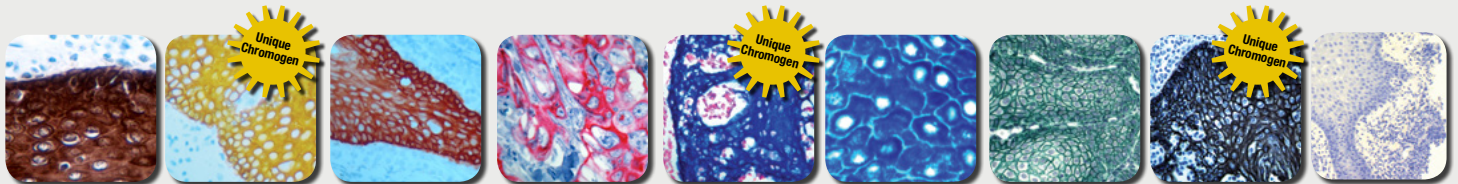
**Enzo's POLYVIEW® PLUS HRP Detection  
(p16 is negative)**



Reference: False-positive results in diagnostic immunohistochemistry are related to horseradish peroxidase conjugates in commercially available assays: G. Nuovo; Ann. Diagn. Pathol. 25, 54 (2016)

## AN UNRIVALED SELECTION OF HIGH DEFINITION CHROMOGENS

Enzo offers the broadest palette of colors to choose from, giving you the ultimate flexibility in planning your multiplex IHC protocol.



**HIGHDEF®  
DAB  
(Brown) HRP**

**HIGHDEF®  
Yellow  
HRP**

**HIGHDEF®  
AEC  
(Red) HRP**

**HIGHDEF®  
Red  
AP**

**HIGHDEF®  
Blue  
HRP**

**HIGHDEF®  
Blue  
AP**

**HIGHDEF®  
Green  
AP**

**HIGHDEF®  
Black  
HRP**

**HIGHDEF®  
Hematoxylin**

ENZ-ACC105

ADI-950-170

ADI-950-210

ADI-950-140  
ADI-950-141

ADI-950-151

ADI-950-150

ENZ-ACC130

ADI-950-171

ENZ-ACC106

[www.enzolifesciences.com/IHC](http://www.enzolifesciences.com/IHC)



# CLEAR AND CONSISTENT RESULTS

## POLYVIEW® PLUS Nanopolymer Peroxidase Detection Solutions

### *Complete Kits with High-Sensitivity and Low Background Deliver Clear Results*

- Complete ready-to-use staining kits and reagents for IHC and ISH
- High sensitivity POLYVIEW® PLUS reagents' biotin-free nanopolymer detection circumvents endogenous background while providing specific antigen localization
- HIGHDEF® DAB (Brown) chromogen provides high intensity color development to deliver sharp, crisp staining
- Compatible with manual and automated platforms

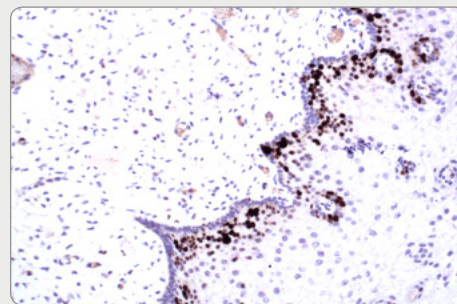
#### **Kits Include:**

Antigen Retrieval Reagent	HIGHDEF® Hematoxylin
IHC Peroxidase Block	Antibody Blocking Buffer
POLYVIEW® PLUS HRP (Anti-Rabbit or Anti-Mouse) Reagent (Ready-to-Use)	HIGHDEF® DAB Substrate/Chromogen

#### **NANOPOLYMER DETECTION KITS**

Product Name	Product #	Size
POLYVIEW® PLUS HRP-DAB (Anti-Rabbit) Kit	ENZ-KIT159	150 Tests
POLYVIEW® PLUS HRP-DAB (Anti-Mouse) Kit	ENZ-KIT160	150 Tests

### **Clear, Consistent Results Eliminate Ambiguity**



Formalin-fixed paraffin-embedded cervical tissue stained using POLYVIEW® PLUS HRP-DAB (Anti-Rabbit) Kit to detect Ki-67 on an autostainer (Leica Bond Rx).

#### **NANOPOLYMER DETECTION REAGENTS**

Product Name	Product #	Size
POLYVIEW® PLUS HRP (Anti-Rabbit) Reagent*	ENZ-ACC103	150 Tests
POLYVIEW® PLUS HRP (Anti-Mouse) Reagent*	ENZ-ACC104	150 Tests
POLYVIEW® PLUS AP (Anti-Rabbit) Reagent*	ENZ-ACC110	150 Tests
POLYVIEW® PLUS AP (Anti-Mouse) Reagent*	ENZ-ACC114	150 Tests

\* Ready-to-Use Reagent

## SAVIEW® PLUS Streptavidin Detection Solutions

### *Streptavidin-Based Detection with Increased Sensitivity and Lower Background*

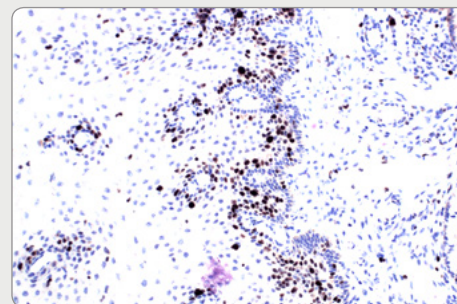
- Reduced carbohydrate and biotin background
- Optimized ready-to-use reagents for IHC and ISH detection
- Compatible with both manual and automated platforms

#### **STREPTAVIDIN DETECTION REAGENTS**

Product Name	Product #	Size
SAVIEW® PLUS HRP Reagent*	ENZ-ACC102	150 Tests
SAVIEW® PLUS AP Reagent*	ENZ-ACC111	150 Tests

\* Ready-to-Use Reagent

### **Strong Signal and Low Background Provides Clear Results**



Formalin-fixed paraffin-embedded cervical tissue stained using SAVIEW® PLUS HRP Reagent to detect Ki-67 on an autostainer (Leica Bond Rx).

## MULTIVIEW® PLUS Multiplex IHC Detection

### *Superior Multiplex Detection*

- Biotin-free nanopolymer-based system for superior multiplex antigen detection
- Complete kit includes HIGHDEF® Chromogens for sensitive, crisp 2-color development
- Preserves precious tissue while reducing reagent and labor costs
- Expandable with additional chromogens (sold separately)

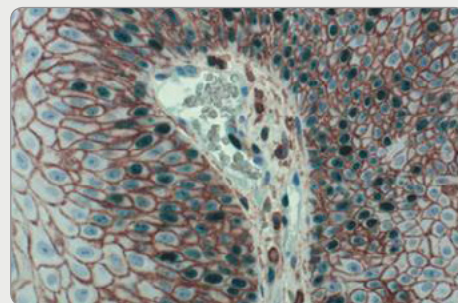
#### **Kit Includes:**

Antigen Retrieval Reagent  
IHC Peroxidase Block  
Antibody Blocker/Diluent  
MULTIVIEW® PLUS HRP & AP Detection Reagents  
HIGHDEF® DAB Chromogen/Substrate  
HIGHDEF® Green AP Chromogen/Substrate  
Hematoxylin  
Empty Mixing Bottle

#### **MULTIPLEX DETECTION KIT**

Product Name	Product #	Size
MULTIVIEW® PLUS (Mouse-HRP/Rabbit-AP) IHC Kit (Brown/Green)	ENZ-KIT181	150 Tests
MULTIVIEW® (Mouse-HRP/Mouse-AP) IHC Kit	ADI-950-101	100 Tests

### **Detection of Multiple Markers on a Single Tissue Section**



Mouse anti-CD44 clone SFF-304 and Rabbit anti-Ki-67 clone MIB-1 was used with MULTIVIEW® PLUS (mouse- HRP/rabbit-AP) IHC Kit (Brown/Green) to perform multiplex IHC.

*You may also be interested in:*



**Tumor Phenotyping and Immune Cell Regulation using Multiplex IHC**



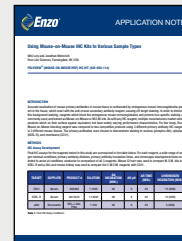
**Three-color IHC Staining on Free- Floating Sections from Human Brain**



**Characterization of Immune Cell Infiltrates in Human Cardiac Allograft Vasculopathy Using MULTIVIEW® IHC Kit**



**Compatibility of POLYVIEW® IHC Detection Reagents and HIGHDEF® IHC Chromogens with Automated IHC Slide Stainers**



**Using Mouse-on-Mouse IHC Kits in Various Sample Types**



**Global Headquarters**  
**ENZO LIFE SCIENCES, INC.**  
10 Executive Blvd.  
Farmingdale, NY 11735  
Ph: 800.942.0430  
Fax: 631.694.7501  
info-usa@enzolifesciences.com

**European Sales Office**  
**ENZO LIFE SCIENCES (ELS) AG**  
Industriestrasse 17  
CH-4415 Lausen, Switzerland  
Ph: +41 61 926 8989  
Fax: +41 61 926 8979  
info-eu@enzolifesciences.com

## LOCAL EUROPEAN OFFICES

**Belgium, The Netherlands & Luxembourg**  
Enzo Life Sciences BVBA  
Avenue Louise 65/Box 11  
1050 Bruxelles  
Belgium  
Ph: +32 3 466 0420  
Fax: +32 3 808 7033  
info-be@enzolifesciences.com

**France**  
Enzo Life Sciences (ELS) AG  
Branch Office Lyon  
13, avenue Albert Einstein,  
F-69100 Villeurbanne, France  
Ph: +33 472 440 655  
Fax: +33 481 680 254  
info-fr@enzolifesciences.com

**Germany**  
Enzo Life Sciences GmbH  
Basler Strasse 57a  
DE-79540 Lörrach  
Germany  
Ph: +49 7621 5500 526  
Fax: +49 7621 5500 527  
info-de@enzolifesciences.com

**UK & Ireland**  
Enzo Life Sciences (UK) Ltd.  
1 Colleton Crescent  
Exeter EX2 4DG  
Ph: 0845 601 1488 (UK customers)  
Ph: +44 1392 825900  
Fax: +44 1392 825910  
info-uk@enzolifesciences.com

For local distributors and detailed product information visit us online:  
**www.enzolifesciences.com**

**Put our experience to work for you!**

Our broad range of scientific expertise and manufacturing capabilities enables us to provide innovative tools for IHC analysis to save you time and money!

MB11-021919