

# HIGHDEF® PLUS Red AP Chromogen Kit

Improved, high definition IHC/ISH  
chromogen producing brilliant red color for  
AP-activated development.

HIGHDEF® PLUS Red AP Chromogen Kit is intended to be used in conjunction with alkaline phosphatase (AP)-based immunostaining or in situ hybridization systems.

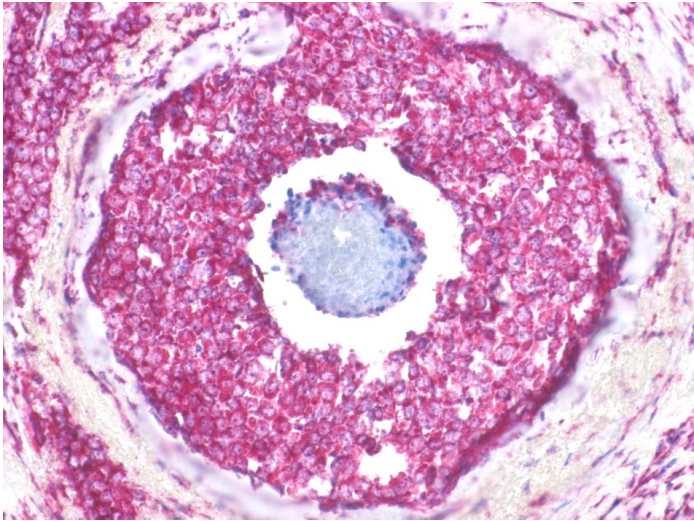
## Ordering Information

[Order Online »](#)

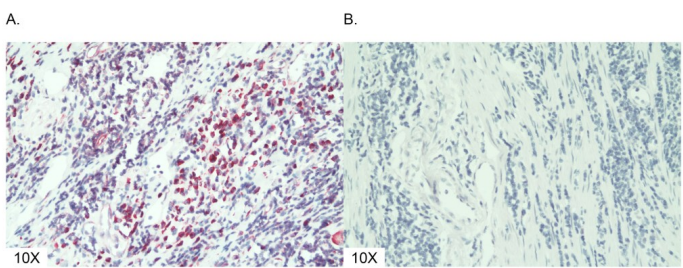
ENZ-ACC158-0030	30ml
-----------------	------

## Manuals, SDS & CofA

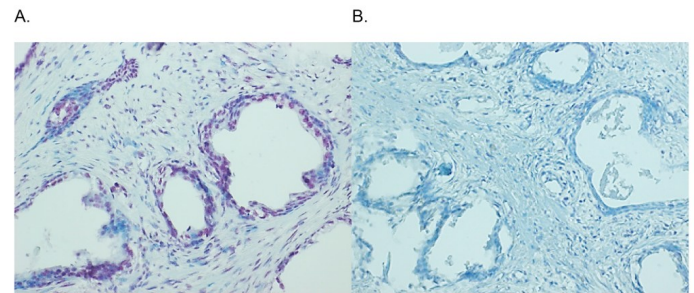
[View Online »](#)



HER-2 protein (red) detected in breast cancer tissue with POLYVIEW® PLUS (anti-rabbit) AP (ENZ-ACC110), combined with HIGHDEF® Red AP chromogen kit (ENZ-ACC158) and hematoxylin (ENZ-ACC106).



In-situ hybridization detection of Ki-67 (red) in ovarian cancer tissue specimens with A. AMPIVIEW® Ki67 (AS) Dig RNA probes (ENZ-GEN271) and B. AMPIVIEW® NSP Dig RNA probes. Tissue samples were pre-treated with antigen retrieval reagent, pH 9 (ENZ-ACC113), Proteinase K (ENZ-33801), and AMPIVIEW® Ki67 (AS) Dig RNA probes (ENZ-GEN271) were amplified with DIGX® mouse anti-digoxigenin linker (RTU) (ENZ-ACC304) and detected with POLYVIEW® PLUS (anti-mouse) AP (ENZ-ACC114), combined with HIGHDEF® Red AP chromogen kit (ENZ-ACC158) and hematoxylin (ENZ-ACC106).



In situ hybridization detection of PTEN (red) in prostate cancer tissue with A. AMPIVIEW® PTEN (AS) DNP RNA Probes (ENZ-GEN131) and B. AMPIVIEW® NSP DNP RNA Probes (no signal). Tissue samples were pre-treated with antigen retrieval reagent, pH 9 (ENZ-ACC113), Proteinase K (ENZ-33801), and AMPIVIEW® PTEN (AS) DNP RNA probes (ENZ-GEN131) were amplified with AMPIVIEW® anti-DNP Linker (RTU) (ENZ-ACC928) and detected with POLYVIEW® PLUS (anti-rabbit) AP (ENZ-ACC110), combined with HIGHDEF® Red AP chromogen kit (ENZ-ACC158) and hematoxylin (ENZ-ACC106).

# Handling & Storage

Long Term Storage            +4°C

Shipping                        Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Application                    IHC, ISH (in situ hybridization)

Contents                        HIGHDEF<sup>®</sup> PLUS Red AP Substrate, HIGHDEF<sup>®</sup> PLUS Signal Enhancer, HIGHDEF<sup>®</sup> PLUS Red AP Chromogen

**Technical Info / Product Notes**      The HIGHDEF<sup>®</sup> PLUS Red AP Chromogen kit is a substrate-enhancer-chromogen system designed to be used for either IHC or ISH when using alkaline phosphatase detection. The system produces a brilliant bright red color and is insoluble in organic solvents; therefore, sections can be dehydrated in alcohol, cleared in xylene (or a xylene-substitute), and permanently mounted. This chromogen substrate system may be used for manual use.



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)