

SAVIEW[®] PLUS HRP

reagent

High sensitivity, low background
streptavidin-based nanopolymer detection
reagent for use with HIGHDEF[®] chromogens
in ISH and IHC applications

SAVIEW[®] PLUS HRP reagent is a ready-to-use streptavidin-based nanopolymer detection reagent suitable for immunohistochemical detection of antigens in formalin-fixed paraffin-embedded tissues and frozen sections. The detection reagent may also be used with blood smears, cytosmears, and cell preparations. In addition, the SAVIEW[®] PLUS HRP reagent can be used for *in situ* hybridization detection of biotinylated DNA/RNA probes.

Citations: 1

[View Online »](#)

Ordering Information

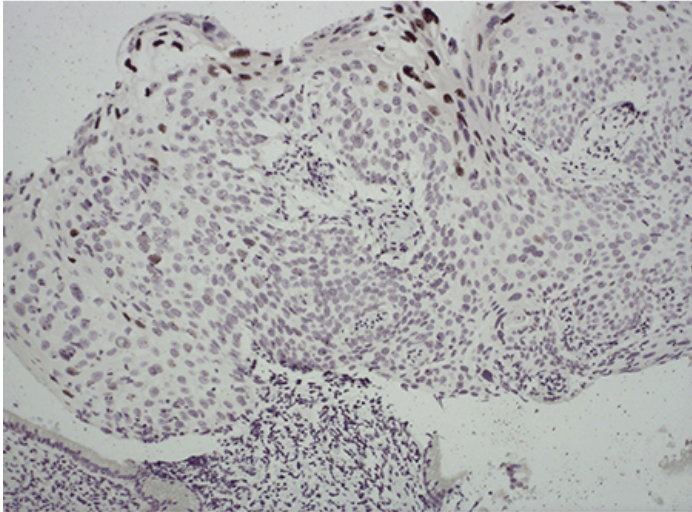
[Order Online »](#)

ENZ-ACC102-0150	150 tests
-----------------	-----------

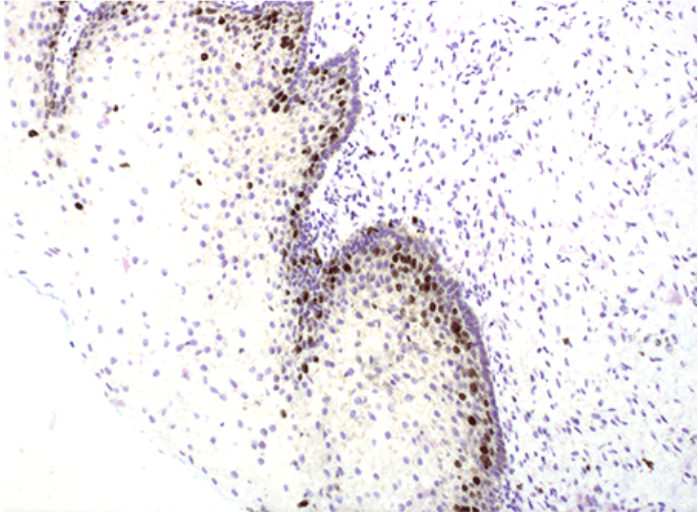
Manuals, SDS & CofA

[View Online »](#)

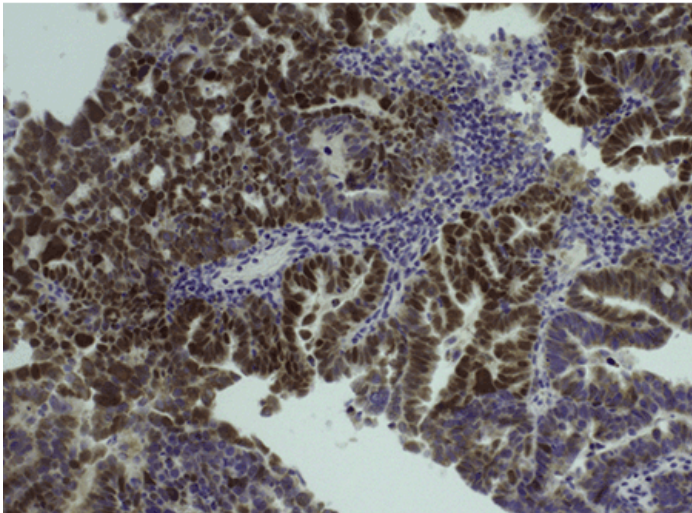
- Ready-to-use reagent
- Streptavidin-based detection with reduced carbohydrate and biotin background
- Suitable for both manual and automated platforms
- Suitable for custom multiplex IHC protocols
- The Streptavidin Blocker/Diluent further reduces background (sold separately)
- Customize with HIGHDEF[®] chromogens of your choice (sold separately)



Cervical tissue screened with PATHO-GENE[®] HPV Type 16/18/31/33/51 (ENZ-32882), followed by SAVIEW[®] PLUS HRP Reagent



Cervical tissue stained with anti-Ki-67 antibody (mouse), then biotinylated anti-mouse secondary antibody, and followed by SAVIEW[®] PLUS HRP reagent



Uterine tissue stained with biotinylated p53 antibody, followed by SAVIEW[®] PLUS HRP reagent

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Do not freeze.
Long Term Storage	+4°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application	IHC, ISH (in situ hybridization)
Quantity	150 tests × 150 µL (22.5 mL total volume)

Technical Info / Product Notes

The SAVIEW[®] PLUS HRP reagent has been developed by labeling streptavidin molecules with a nanopolymer of horse-radish peroxidase (HRP) using proprietary technology. This ensures consistent and reproducible detection of biotinylated primary/secondary antibodies or biotinylated nucleic acid probes on tissue sections.

The Streptavidin Blocker/Diluent (ENZ-ACC109) further reduces background and is suitable for diluting the SAVIEW[®] PLUS HRP reagent.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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