

# SAVIEW<sup>®</sup> PLUS AP

## Reagent

High sensitivity, low background,  
streptavidin-based nanopolymer detection  
reagent for use with HIGHDEF<sup>®</sup> chromogens  
in ISH and IHC applications

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

### Ordering Information

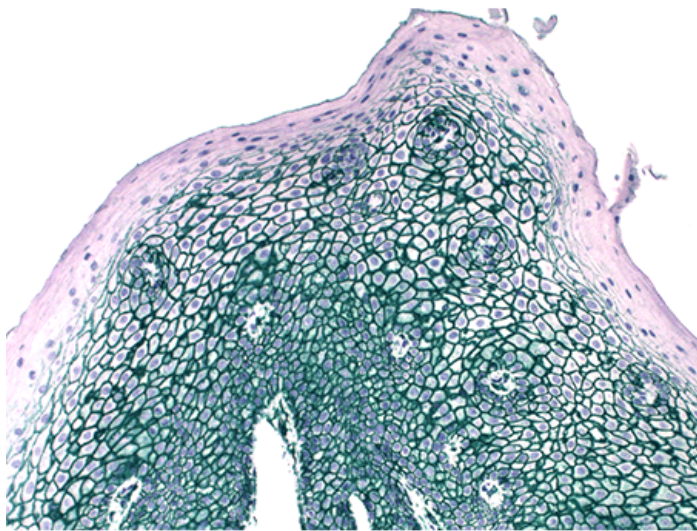
[Order Online »](#)

ENZ-ACC111-0150	150 tests
-----------------	-----------

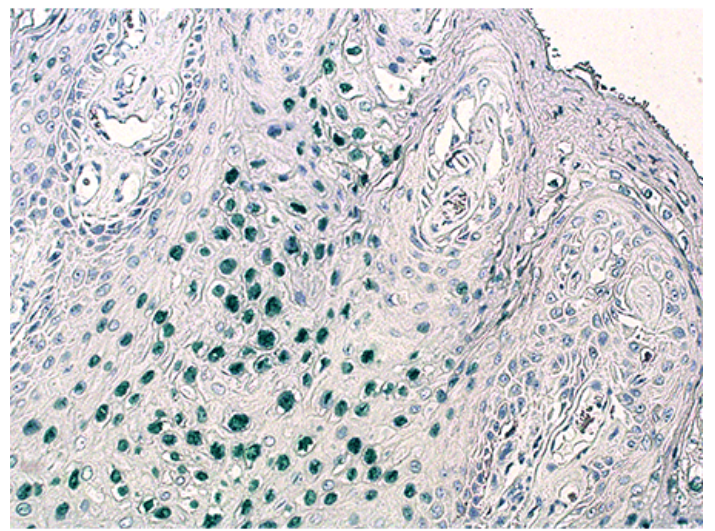
### Manuals, SDS & CofA

[View Online »](#)

- Ready-to-use reagent
- Streptavidin-based detection with reduced carbohydrate and biotin background
- Suitable for both manual and automated platforms
- Streptavidin Blocker/Diluent further reduces background (sold separately)
- 5-10X more sensitive than avidin based detection
- Customize with HIGHDEF<sup>®</sup> chromogens of your choice (sold separately)



Mouse anti-CD44 and SAVIEW® PLUS AP (Enzo Life Sciences) were used to perform IHC: Manual method. Sections were stained with HIGHDEF® Green AP Chromogen/Substrate and counterstained with HIGHDEF® Hematoxylin. Images acquired under 10X objective.



SAVIEW® PLUS AP (Enzo Life Sciences) was used to perform ISH: Manual method. Sections were stained with HIGHDEF® Green AP Chromogen/Substrate and counterstained with HIGHDEF® Hematoxylin. Images acquired under 10X objective.

## 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)
Application Notes	IHC ISH ( <i>in situ</i> hybridization)
Quantity	150 tests × 150 µL (22.5 mL total volume)
Technical Info / Product Notes	<p>The SAVIEW<sup>®</sup> PLUS AP Reagent has been developed by labeling streptavidin molecules with a nanopolymer of alkaline phosphatase (AP) using proprietary technology. This ensures consistent and reproducible detection of biotinylated primary/secondary antibodies or biotinylated nucleic acid probes on tissue sections.</p> <p>The Streptavidin Blocker/Diluent (ENZ-ACC109) further reduces background and is suitable for diluting the SAVIEW<sup>®</sup> PLUS AP Reagent.</p>

Last modified: May 29, 2024



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)