

Streptavidin (AP Conjugate)

As part of the need to increase the sensitivity of your sandwich ELISA, this streptavidin AP conjugate can be used with your sandwich ELISA that utilizes a biotinylated detector antibody and the AMP'D[®] ELISA Signal Amplification kit (Prod. no. ENZ-KIT-100) which requires the presence of the AP as the conjugate.

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

ENZ-PRT119-1000	1ml
-----------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Formulation

Liquid (50 µg/ml). In 0.025M TRIS-HCl, pH 8.0, containing 0.075M sodium chloride, 0.05mM ZnSO₄, 2.45mM MgCl₂, 0.045% sodium azide, 0.05% BSA and 50% Stabilzyme-AP.

Purity Detail

Alkaline phosphatase purified from calf intestine.

Source

Native. Purified from *Streptomyces avidinii* (SA) and calf intestine (AP).

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)