

AMP'D[®] ELISA Signal Amplification Kit

AMP'D[®] ELISA Signal Amplification Kit provides up to 50-fold increase in sensitivity over traditional ELISAs while detecting lower concentrations of target in samples.

The AMP'D[®] ELISA Signal Amplification kit is designed to replace traditional alkaline phosphatase (AP) substrates, such as pNPP (p-Nitrophenyl phosphate), with a combination substrate and amplifier system that results in greater sensitivity when compared to a classic substrate enzyme linked-immunosorbent assay (ELISA).

In the AMP'D[®] ELISA Signal Amplification kit, bound AP converts a substrate that is utilized in a second enzyme reaction system which is initiated by addition of the amplifier reagent. It is this amplification step that allows for greater (amplified) color production at lower analyte concentrations resulting in an increase in assay sensitivity.

Ordering Information

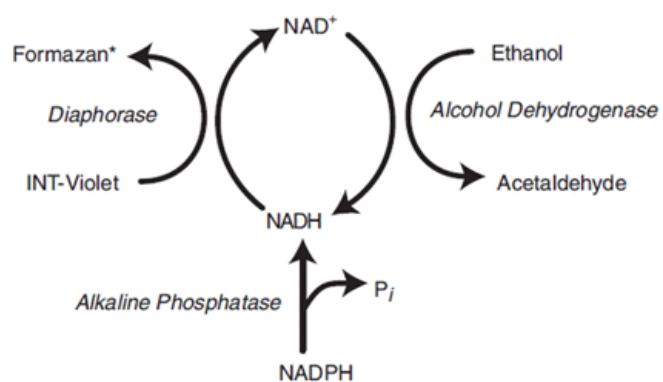
[Order Online »](#)

| | |
|------------------|-----------------------|
| ENZ-KIT-100-0005 | 5 assays - 5x96 wells |
|------------------|-----------------------|

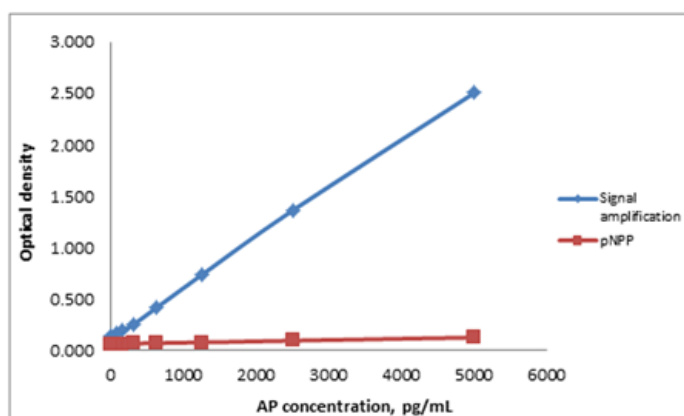
Manuals, SDS & CofA

[View Online »](#)

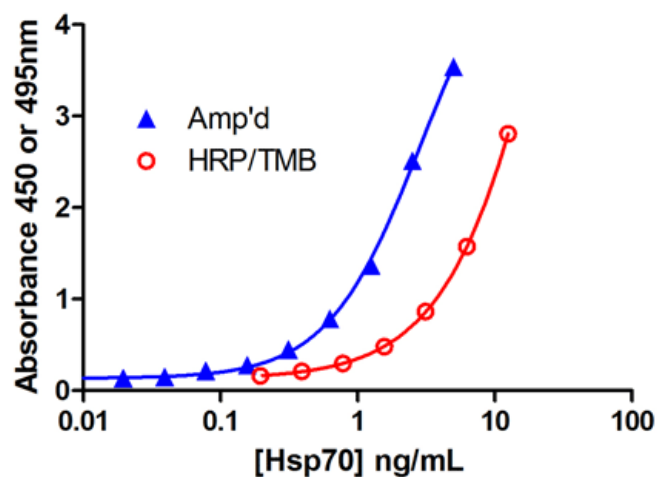
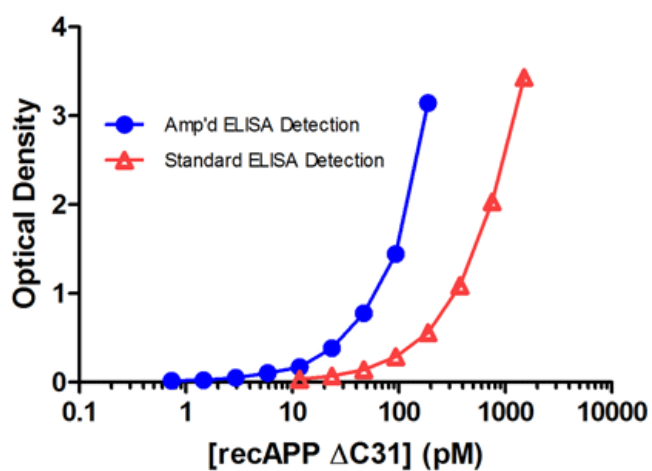
- Designed to amplify sandwich/immunometric ELISA signal
- Increase ELISA sensitivity greater than 10-fold
- Flexibility to work with own targets
- Convenient one or five 96-well plate formats for high throughput analysis
- Easy-to-use, simple procedure with results in just 30 minutes

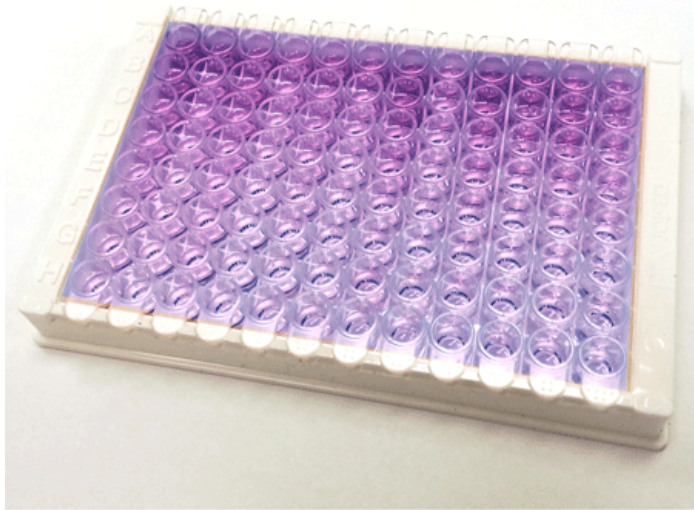


*Colored end product



recAPP Δ C31 Standard Curve





Handling & Storage

| | |
|-------------------|--|
| Handling | Upon receipt, the kit should be stored at 4°C. |
| Long Term Storage | +4°C |
| Shipping | Blue Ice |

Regulatory Status

RUO - Research Use Only

Product Details

| | |
|-------------------|---|
| Application | Colorimetric detection, ELISA |
| Application Notes | <p>For the amplification of ELISA sensitivity designed to replace traditional alkaline phosphatase (AP) substrates.</p> <p>Assay should be read at a wavelength of 495nm.</p> |
| Assay Time | 30 minutes |
| Compatibility | This product is compatible with the Absorbance 96 Plate Reader . |
| Contents | Signal Amplification Substrate, Amplifier, Substrate Diluent, and Amplifier Diluent |
| Sensitivity | Improves ELISA sensitivity ~10 to 50-fold. |
| Wavelength | 495 nm |

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-
fr@enzolifesciences.com](mailto:info-fr@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)