

AMPIVIEW® Engrailed 1 (AS) DNP RNA Probes (Xenopus laevis) Set

AMPIVIEW® Engrailed 1 (AS) DNP RNA Probes (*Xenopus laevis*) are DNP-labeled RNA probes targeting the nucleic acid (DNA/RNA or RNA) for Engrailed 1 (EN1) for *in situ* hybridization in tissues and cells. The probes are formulated in a buffered formamide solution with hybridization enhancers.

Ordering Information

[Order Online »](#)

| | |
|-----------------|-----|
| ENZ-GEN346-2000 | 2ml |
|-----------------|-----|

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

| | |
|--------------------|---|
| Use/Stability | Aliquot and store at -20°C. Under these conditions, products are stable until its expiration dates. |
| Handling | Allow contents to warm up to room temperature prior to use. |
| Short Term Storage | -20°C |
| Long Term Storage | -20°C |
| Shipping | Dry Ice |

Regulatory Status

RUO - Research Use Only

Product Details

| | |
|--------------------------------|--|
| Application | ISH (in situ hybridization) |
| Application Notes | AMPIVIEW® Engrailed 1 (AS) DNP RNA probes (<i>Xenopus laevis</i>) have been designed to detect the nucleic acid for Engrailed 1 (EN1) gene in <i>Xenopus laevis</i> or African clawed frogs. AMPIVIEW® Engrailed 1 (AS) DNP RNA probes (<i>Xenopus laevis</i>) are conjugated with dinitrophenol (DNP) labels and have been optimized to produce clear results with Enzo's AMPIVIEW® anti-DNP linker and nanopolymner-based detection systems, POLYVIEW® PLUS, combined with HIGHDEF® chromogens and counterstain to produce clear results that can be visualized with a light microscope. |
| Contents | 2 x 1 mL AMPIVIEW® Engrailed 1 (AS) DNP RNA Probes (<i>Xenopus laevis</i>) (2µg/mL) 1 x 2 mL AMPIVIEW® Hybridization Buffer (1X) |
| Technical Info / Product Notes | AMPIVIEW® Engrailed 1 (AS) DNP RNA Probes (<i>Xenopus laevis</i>) are optimized for detection of the nucleic acid (RNA/DNA or RNA) of African clawed frog's Engrailed 1 in FFPE tissues or cells with AMPIVIEW® anti-DNP linker, POLYVIEW® PLUS detection reagents combined with HIGHDEF® chromogens and counterstains (linker and detection solutions not included). Dilutions and concentration optimization can be done with AMPIVIEW® Hybridization Buffer (1X), included in this set. |



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)