

AMPIVIEW® PD-1 (AS) Dig RNA Probes Set

A digoxigenin-labeled RNA probe for the *in situ* hybridization (ISH) detection of PD-1 RNA or DNA/RNA in cells and tissues

AMPIVIEW® PD-1 (AS) Dig RNA Probes are digoxigenin-labeled RNA probes targeting PD-1 nucleic acid (DNA/RNA or RNA) for *in situ* hybridization in tissues and cells. The probes are formulated in a buffered formamide solution with hybridization enhancers.

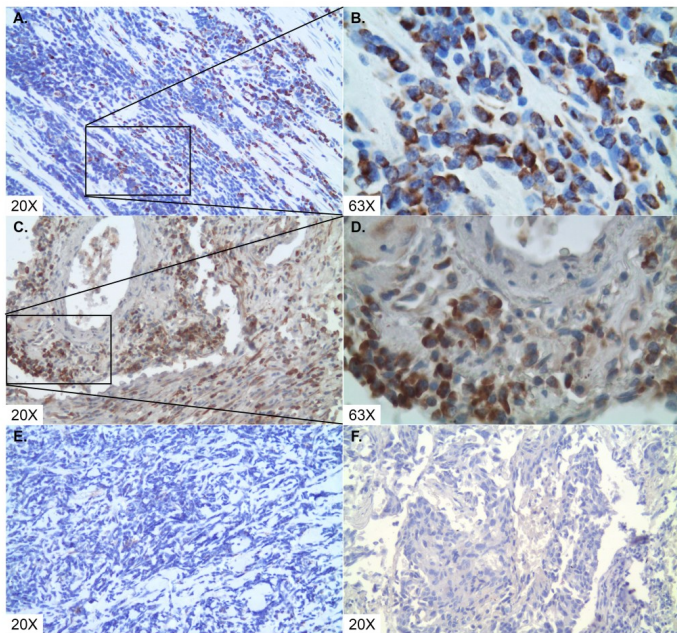
Ordering Information

[Order Online »](#)

| | |
|-----------------|-----|
| ENZ-GEN281-2000 | 2ml |
|-----------------|-----|

Manuals, SDS & CofA

[View Online »](#)



PD-1 (brown) was detected in A-B, E mantic lymphoma and C-D, F lung cancer tissue specimens with A-D AMPIVIEW® PD-1 (AS) Dig RNA probes and E-F AMPIVIEW® NSP Dig RNA Probes, amplified with DIGX® rabbit anti-digoxigenin linker and detected with POLYVIEW® PLUS (anti-r) HRP reagent combined with HIGHDEF® DAB chromogen/substrate and counterstained with HIGHDEF® hematoxylin.

Handling & Storage

| | |
|--------------------|--|
| Use/Stability | Upon receipt, aliquot and store at -20°C. For long-term storage, store at -80°C. These products, under these conditions, are stable until the expiration date. |
| Handling | Allow contents to warm up to room temperature prior to use. |
| Short Term Storage | -20°C |
| Long Term Storage | -80°C |
| Shipping | Dry Ice |

Regulatory Status

RUO - Research Use Only

Product Details

| | |
|--------------------------------|---|
| Application | ISH (in situ hybridization) |
| Application Notes | AMPIVIEW® PD-1 (AS) Dig RNA probes have been designed to detect the nucleic acid that encodes Program Cell Death Protein 1 (PD-1) in tissues and cells. PD-1 is a cell surface receptor on T cells and B cells that plays a role in the regulation of the immune system. AMPIVIEW® PD-1 (AS) Dig RNA probes are conjugated with digoxigenin and have been optimized to be amplified with DIGX® anti-digoxigenin linker and detected with 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® PD-1 (AS) Dig RNA Probes (2µg/mL) 1 x 2 mL AMPIVIEW® Hybridization Buffer (1X) |
| Technical Info / Product Notes | AMPIVIEW® PD-1 (AS) Dig RNA Probes are optimized for detection of PD-1 RNA or RNA/DNA in FFPE tissues or cells with DIGX® anti-digoxigenin 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. |

Last modified: April 16, 2025



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)