

AMPIVIEW® PD-L1 (AS) Dig RNA Probes Set

A digoxigenin-labeled RNA probe for the detection of human PD-L1 for *in situ*

hybridization.

AMPIVIEW® RNA probes are uniquely designed with the precision of targeted, sequence-specific RNA probes, powered by Enzo's LoopRNA™ ISH technology to deliver superior sensitivity. AMPIVIEW® PD-L1 (AS) Dig RNA Probes are digoxigenin-labeled probes targeting Human PD-L1 for *in situ* hybridization. The probes are formulated in a buffered formamide solution with hybridization enhancers.

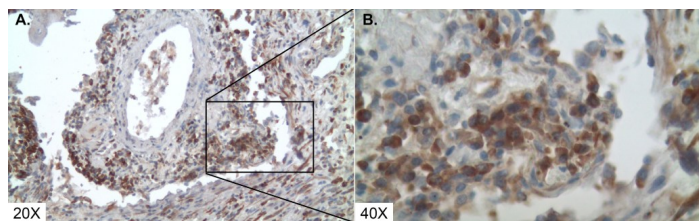
Ordering Information

[Order Online »](#)

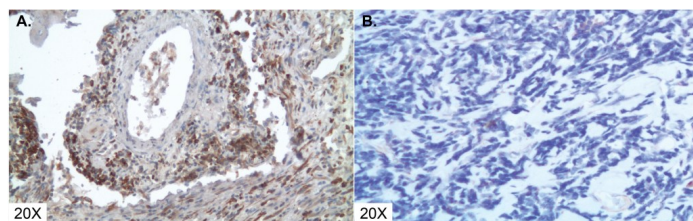
ENZ-GEN283-2000	2ml
-----------------	-----

Manuals, SDS & CofA

[View Online »](#)



PD-L1 (brown) was detected in lung cancer tissue with AMPIVIEW® PD-L1 (AS) 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. A. 20X Image, B. 40X Image.



PD-L1 (brown) was detected in lung cancer tissue with A. AMPIVIEW® PD-L1 (AS) Dig RNA probes and B. 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	Aliquot and store at -20°C. Long term storage recommended at -80°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	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application	ISH (in situ hybridization)
Application Notes	<p>AMPIVIEW® PD-L1 (AS) Dig RNA probes have been designed to target the nucleic acid of CD274 gene which encodes the Programmed Death-Ligand 1 (PD-L1) protein, for its detection in tissues and cells. PD-L1 binds to its receptor Programmed Cell Death 1 (PD-1) on T cells to inhibit T cell activation, suppress immune responses and maintenance of immune tolerance. Many tumors overexpress PD-L1 to evade the immune detection and allow for cancer cells to grow and spread. Immunotherapy drugs block PD-L1, restoring T cell activity and enabling the immune system to attack tumors. AMPIVIEW® PD-L1 (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.</p>
Contents	<p>AMPIVIEW® PD-L1 (AS) Dig RNA Probes (ENZ-GEN282) AMPIVIEW® Hybridization Buffer (ENZ-ACC152)</p>

Technical Info / Product Notes

AMPIVIEW® PD-L1 (AS) Dig RNA Probes are optimized for amplification with DIGX® anti-digoxigenin Linker and detection with POLYVIEW® PLUS detection reagents combined with HIGHDEF® chromogens and counterstains (not included). Dilutions and concentration optimization can be done with AMPIVIEW® hybridization buffer, included in this set.

Last modified: September 30, 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)