

# AMPIVIEW® PD-L1 (AS) DNP RNA Probes Set

**A DNP-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) DNP RNA Probes are DNP-labeled probes targeting Human PD-L1 for *in situ* hybridization. The probes are formulated in a buffered formamide solution with hybridization enhancers.

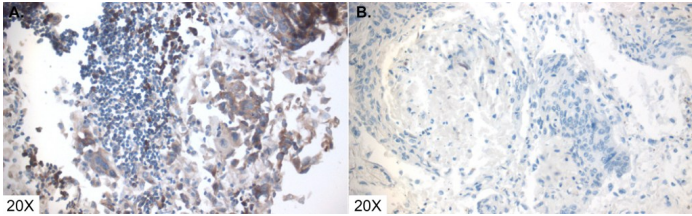
## Ordering Information

[Order Online »](#)

ENZ-GEN291-2000	2ml
-----------------	-----

## Manuals, SDS & CofA

[View Online »](#)



PD-L1 (brown) was detected in lung cancer tissue with A. AMPIVIEW® PD-L1 (AS) DNP RNA probes and B. AMPIVIEW® NSP DNP RNA Probes, amplified with AMPIVIEW® rabbit anti-DNP 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. 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	<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) DNP RNA probes are conjugated with dinitrophenol (DNP) and have been optimized to be amplified with AMPIVIEW® Rabbit anti-DNP 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) DNP RNA Probes (ENZ-GEN290)</p> <p>AMPIVIEW® Hybridization Buffer (ENZ-ACC152)</p>

## Technical Info / Product Notes

AMPIVIEW<sup>®</sup> PD-L1 (AS) DNP RNA Probes are optimized for amplification with AMPIVIEW<sup>®</sup> Rabbit anti-DNP Linker and detection with POLYVIEW<sup>></sup><sup>®</sup> PLUS detection reagents combined with HIGHDEF<sup>®</sup> chromogens and counterstains (not included). Dilutions and concentration optimization can be done with AMPIVIEW<sup>®</sup> 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)