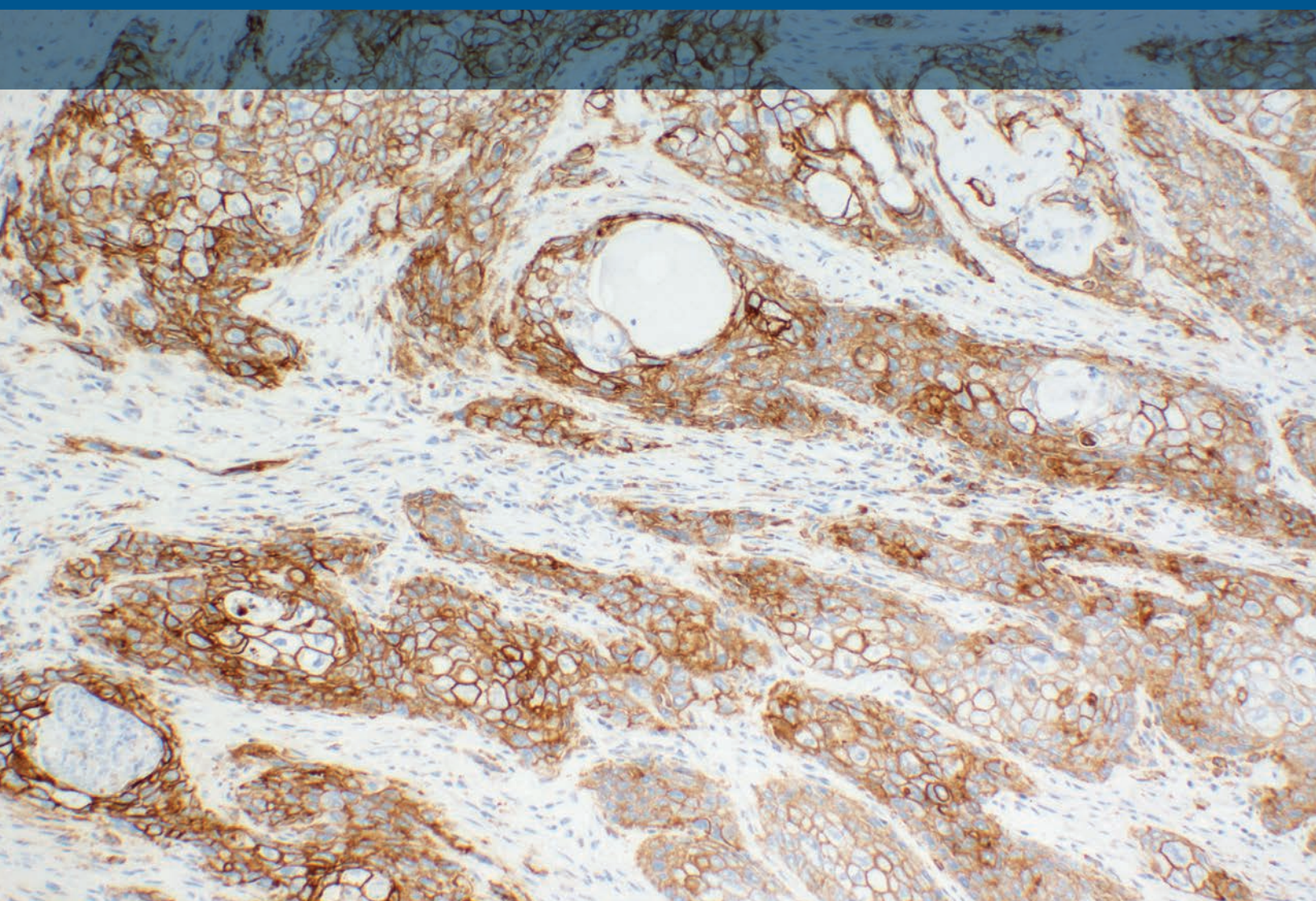


# PD-L1 IHC 28-8 pharmDx Interpretation Manual – Urothelial Carcinoma Including Muscle Invasive Urothelial Carcinoma (% Tumor Cell Expression)

IVD marked for in vitro diagnostic use (Worldwide version)

Rx



**Agilent**

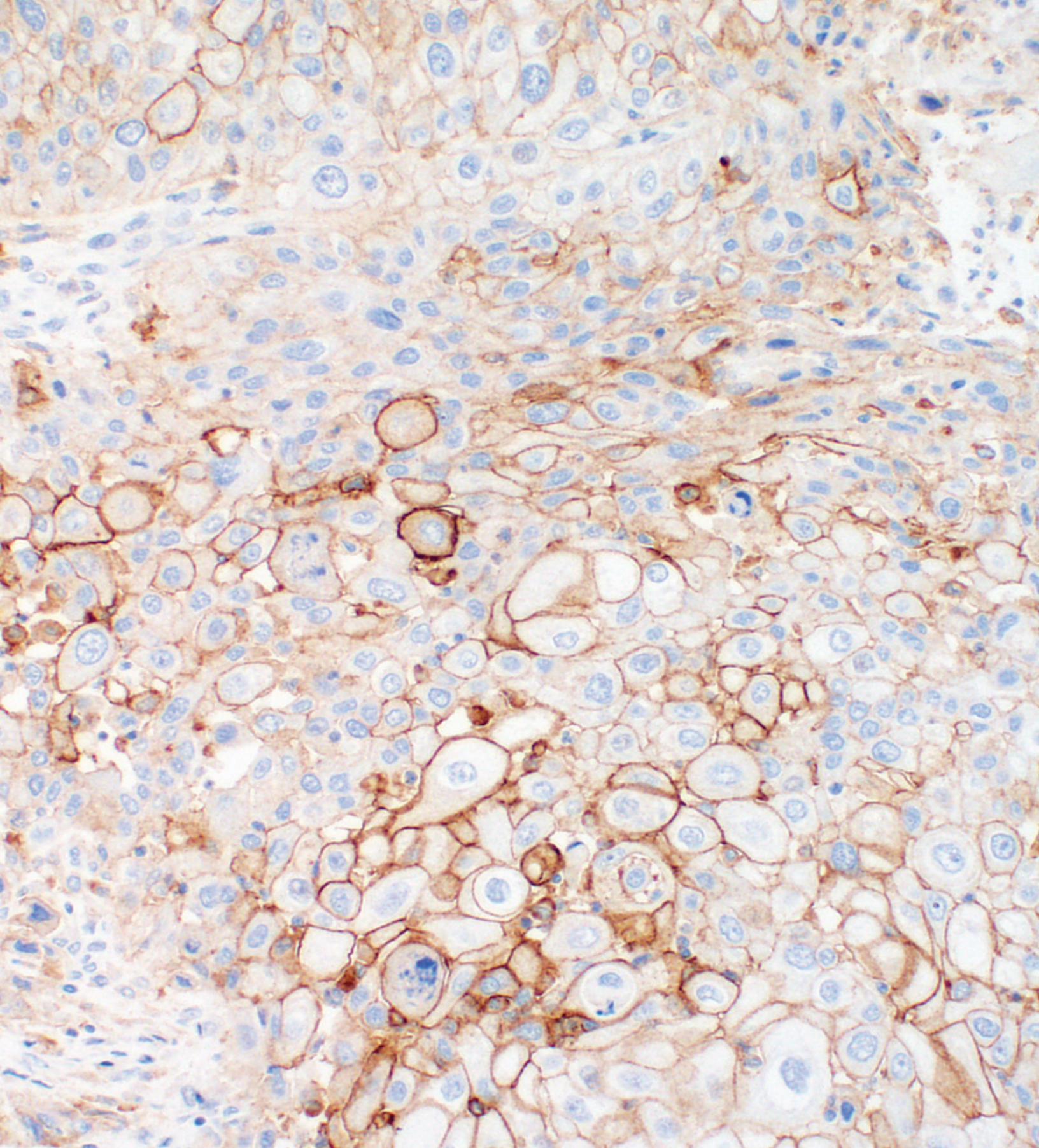
Trusted Answers



# Table of Contents

<b>Introduction</b>	<b>5</b>
Intended Use	5
PD-L1 IHC 28-8 pharmDx Interpretation Manual - Overview	5
<b>Acknowledgment</b>	<b>6</b>
<b>PD-L1 Overview</b>	<b>7</b>
<b>The Role of the PD-1/PD-L1 Pathway in Cancer</b>	<b>8</b>
<b>PD-L1 IHC 28-8 pharmDx Overview</b>	<b>9</b>
<b>Technical Considerations for Optimal Performance of PD-L1 IHC 28-8 pharmDx</b>	<b>11</b>
Specimen Collection and Processing	11
Tissue Processing	11
Control Tissues	12
PD-L1 IHC 28-8 pharmDx Staining Procedure	12
Reagent Storage	12
Reagent Preparation	12
Controls to Assess Staining Quality	13
Staining Protocol	13
Deparaffinization, Rehydration, and Target Retrieval	13
Staining and Counterstaining	14
Mounting	14
<b>PD-L1 IHC 28-8 pharmDx Technical Checklist</b>	<b>15</b>
<b>Guidelines for Scoring PD-L1 IHC 28-8 pharmDx in UC and MIUC</b>	<b>16</b>
<b>Slide Evaluation Flowchart</b>	<b>17</b>
<b>Recommendations for Interpretation of PD-L1 IHC 28-8 pharmDx in UC and MIUC</b>	<b>18</b>
Patient Specimen Stained with H&E	18
PD-L1 IHC 28-8 pharmDx Control Slide	18
Positive Control Tissue Slides	19
Negative Control Tissue Slides	19
Patient Specimen Stained with Negative Control Reagent	19
Patient Specimen Stained with Primary Antibody	20
<b>Tips and Considerations</b>	<b>20</b>
Non-evaluable Specimen	20
Indeterminate Specimen	20
<b>PD-L1 IHC 28-8 pharmDx Suggested Scoring Methods for Calculating PD-L1 Percent Tumor Cell Expression</b>	<b>21</b>
<b>Reporting Results</b>	<b>23</b>
<b>PD-L1 IHC 28-8 pharmDx Immunostaining Examples in UC and MIUC</b>	<b>25</b>
<b>Challenging Cases for UC and MIUC PD-L1 IHC 28-8 pharmDx</b>	<b>30</b>
<b>Troubleshooting Guide for PD-L1 IHC 28-8 pharmDx</b>	<b>35</b>
<b>Bibliography</b>	<b>37</b>





PD-L1 IHC 28-8 pharmDx  
Interpretation Manual - UC and MIUC (% Tumor Cell Expression)



# Introduction

## Intended Use

### For In Vitro Diagnostic Use

PD-L1 IHC 28-8 pharmDx is intended for use in specific types of cancer, which may vary by region. Please refer to the local Instructions for Use (IFU) for a description of the types of cancer in which PD-L1 IHC 28-8 pharmDx is intended for use.

PD-L1 IHC 28-8 pharmDx is a qualitative immunohistochemical assay using Monoclonal Rabbit Anti-PD-L1, Clone 28-8 intended for use in the detection of PD-L1 protein in formalin-fixed, paraffin-embedded (FFPE) urothelial carcinoma (UC) and muscle invasive urothelial carcinoma (MIUC) tissues using EnVision FLEX visualization system on Autostainer Link 48.

PD-L1 protein expression in UC and muscle invasive urothelial carcinoma (MIUC) is determined by using % tumor cell expression, which is the percentage of evaluable tumor cells exhibiting partial or complete membrane staining at any intensity.

**Note:** For more details regarding the region-specific device intended use and clinical data, please refer to the IFU ([Instructions For Use | Agilent](#)).

## PD-L1 IHC 28-8 pharmDx Interpretation Manual - Overview

This PD-L1 IHC 28-8 pharmDx Interpretation Manual is provided as a tool to help guide pathologists and laboratory technicians to achieve correct and reproducible results in assessing PD-L1 expression in FFPE UC and MIUC specimens. The goal of this manual is to familiarize you with the requirements for scoring UC and MIUC specimens stained with PD-L1 IHC 28-8 pharmDx. Photomicrographs of example cases are provided for reference.

PD-L1 IHC 28-8 pharmDx Instructions for Use (IFU) contains guidelines and technical tips for ensuring high quality staining in your laboratory.

Review of this PD-L1 IHC 28-8 pharmDx Interpretation Manual will provide a solid foundation for evaluating UC and MIUC specimens stained with PD-L1 IHC 28-8 pharmDx. For more details, please refer to the current local version of PD-L1 IHC 28-8 pharmDx IFU, or visit [www.agilent.com](http://www.agilent.com).

## Assay Interpretation

The clinical interpretation of any staining, or the absence of staining, must be complemented by the evaluation of proper controls. An evaluation must be made by a qualified pathologist within the context of the patient's clinical history and other diagnostic tests. This product is intended for in vitro diagnostic (IVD) use.

## Reporting Results

To help understand what information should be reported to the treating physician, please refer to the Reporting Results section of this manual on page 23-24.

# Acknowledgment

## Photomicrographs

Photomicrograph magnification levels may appear different than indicated in respective annotations due to adjustment of image size.

**Note:** Photomicrographs included in this interpretation manual include specimens provided by the following suppliers:

- Tissue samples supplied by BioIVT (Hicksville, NY, USA).
- The data and UC and MIUC specimens used in this project were provided by Centre Hospitalier Universitaire (CHU) de Nice (Nice, France) with appropriate ethics approval and through Trans-Hit Biomarkers Inc.
- The data and UC and MIUC specimens used in this project were provided by US Biolab (Gaithersburg, MD, USA) with appropriate ethics approval and through Trans-Hit Biomarkers Inc.

## PD-L1 Overview

### **The PD-1/PD-L1 Pathway Controls the Immune Response in Normal Tissue**

Programmed death-ligand 1 (PD-L1) is a transmembrane protein that binds to the programmed death-1 receptor (PD-1) during immune system modulation. The PD-1 receptor is typically expressed on cytotoxic T-cells and other immune cells, while the PD-L1 ligand is typically expressed on normal cells. Normal cells use the PD-1/PD-L1 interaction as a mechanism of protection against immune recognition by inhibiting the action of T-cells (Figure a). Inactivation of cytotoxic T-cells downregulates the immune response such that the inactive T-cell is exhausted, ceases to divide, and might eventually die by programmed cell death, or apoptosis.

### **The Tumor Escapes Detection by Utilizing the PD-1/PD-L1 Pathway**

Many tumor cells are able to upregulate the expression of PD-L1 as a mechanism to evade the body's natural immune response. Activated T-cells recognize the PD-L1 marker on the tumor cell, similar to that of a normal cell, and PD-L1 signaling renders the T-cell inactive (Figure b). The tumor cell escapes the immune cycle, continues to avoid detection for elimination and is able to proliferate.

### **Anti-PD-1 Therapy Enables the Immune Response Against Tumors**

Anti-PD-1 therapy works by blocking the PD-1/PD-L1 interaction between tumor cells and activated T-cells, helping to prevent immunosuppression, thereby enabling cytotoxic T-cells to actively remove tumor cells (Figure c).

## The Role of the PD-1/PD-L1 Pathway in Cancer

### Limiting damage to healthy tissue

Inactivation of T-cells limits damage to healthy tissue.

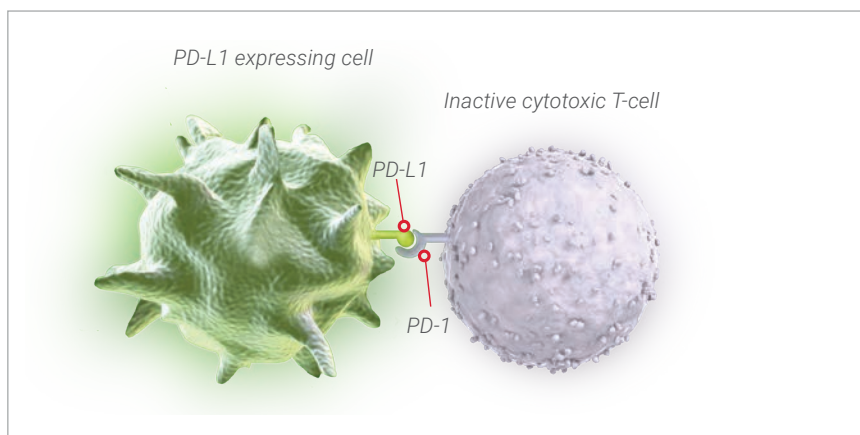


Figure a.

### The tumor escapes detection

Inactivation of T-cells reduces tumor cell killing.

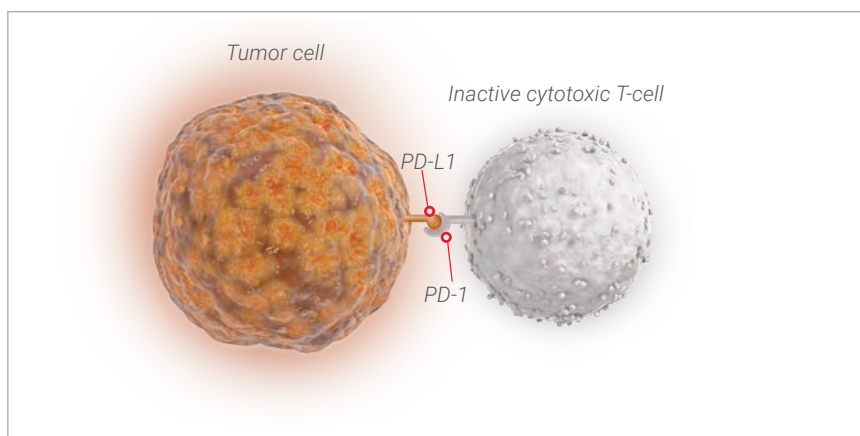


Figure b.

### Immuno-oncology therapies harness the immune response to fight tumors

Blocking PD-1/PD-L1 interaction enables cytotoxic T-cells to actively remove tumor cells.

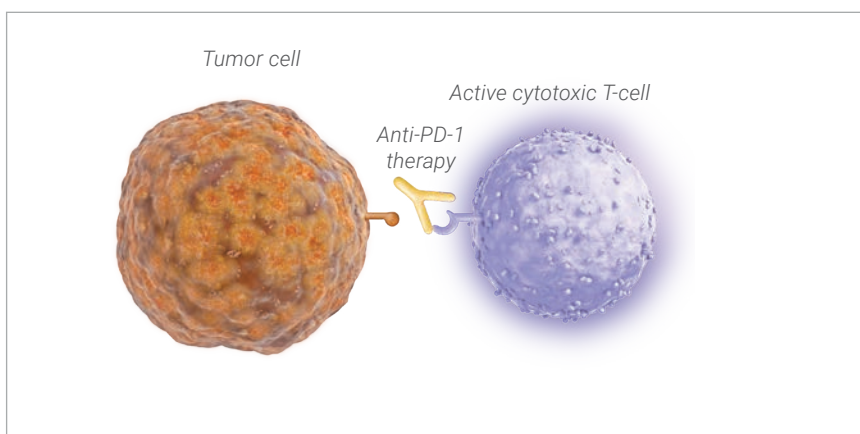


Figure c.



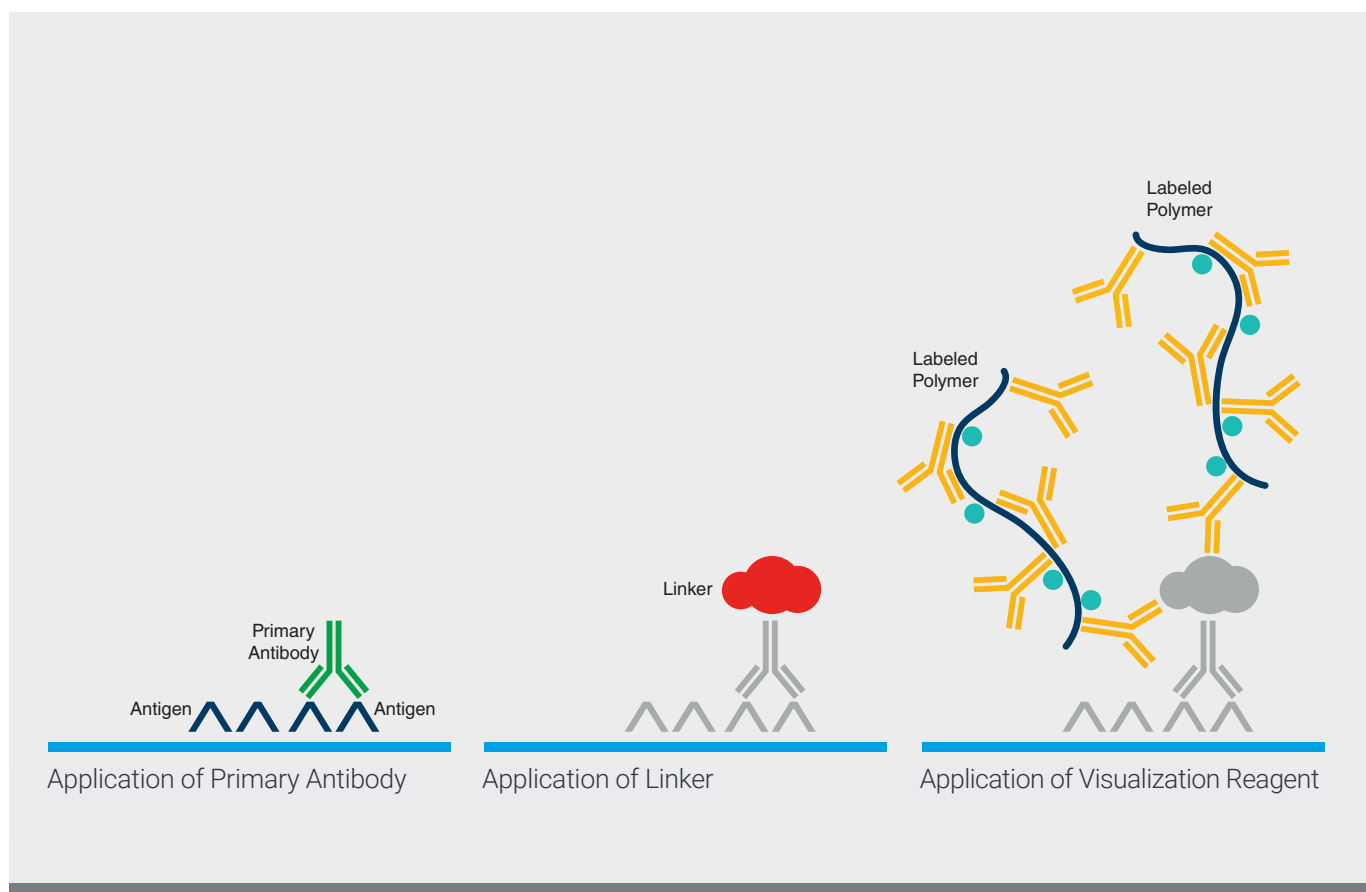
# PD-L1 IHC 28-8 pharmDx Overview

## Code SK005

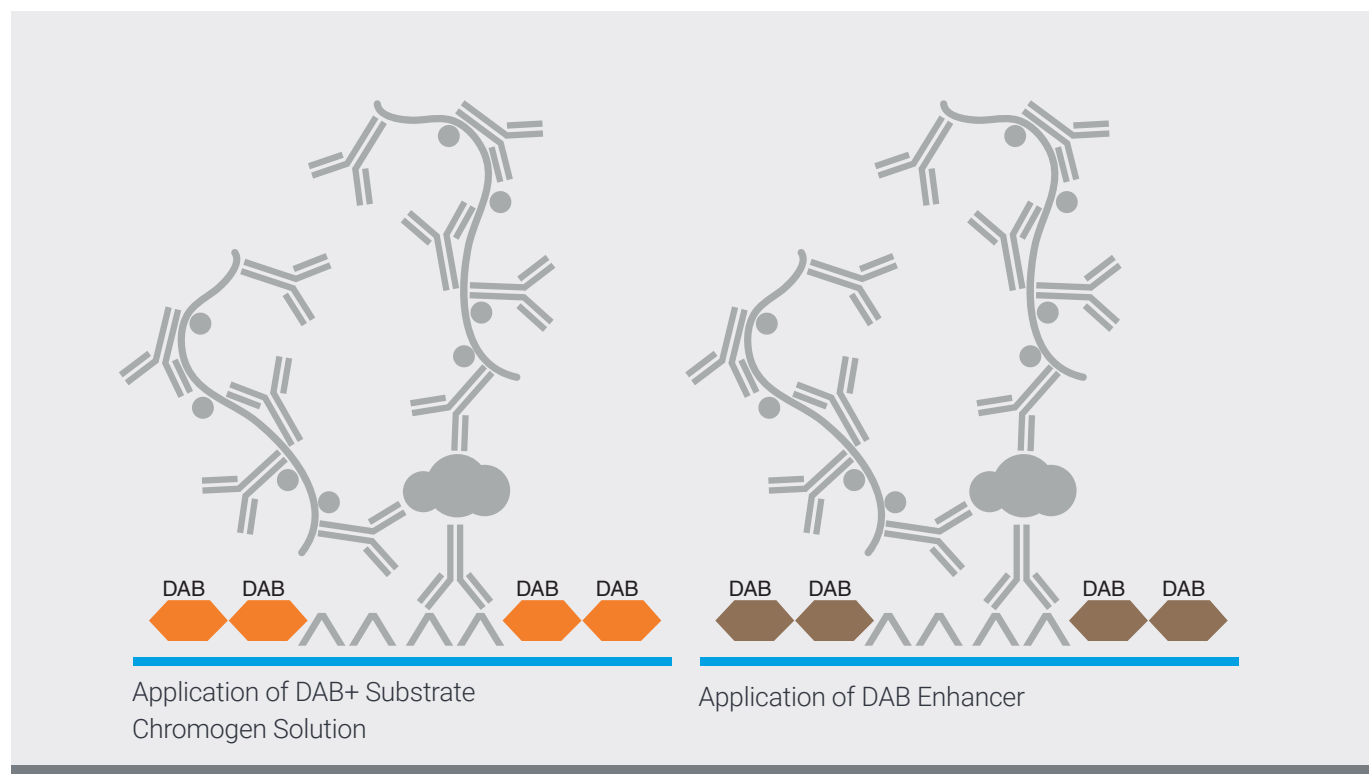
PD-L1 IHC 28-8 pharmDx contains optimized reagents and protocol required to complete an IHC staining procedure of FFPE specimens using PT Link Pre-treatment Module and Autostainer Link 48.

Following incubation with the primary monoclonal antibody to PD-L1 or the Negative Control Reagent (NCR), specimens are incubated with a linker antibody specific to the host species of the primary antibody, and then are incubated with a ready-to-use visualization reagent consisting of secondary antibody molecules and horseradish peroxidase molecules coupled to a dextran polymer backbone (see Figure 1a). The enzymatic conversion of the subsequently added chromogen results in precipitation of a visible reaction product at the site of the antigen. The color of the chromogenic reaction is modified by a chromogen enhancement reagent (see Figure 1b). The specimen may then be counterstained and coverslipped. Results are interpreted using a light microscope. Control Slides containing two FFPE human cell lines are provided to validate staining runs.

## PD-L1 IHC 28-8 pharmDx staining procedure



**Figure 1a.** PD-L1 IHC 28-8 pharmDx staining procedure.



**Figure 1b.** PD-L1 IHC 28-8 pharmDx staining procedure.



**Figure 2.** PD-L1 IHC 28-8 pharmDx staining component.

*All PD-L1 IHC 28-8 pharmDx reagents are to be used on the Autostainer Link 48. All reagents must be used as indicated in the IFU in order for the test to perform as specified.*

PD-L1 IHC 28-8 pharmDx contains reagents to perform 50 tests in up to 15 individual runs (see Figure 2).

- EnVision FLEX Target Retrieval Solution, Low pH, 50x
- Peroxidase-Blocking Reagent
- Primary Antibody: Monoclonal Rabbit Anti-PD-L1, Clone 28-8
- Negative Control Reagent
- PD-L1 IHC 28-8 pharmDx Anti-Rabbit LINKER
- Visualization Reagent-HRP
- DAB+ Substrate Buffer
- DAB+ Chromogen
- DAB Enhancer
- PD-L1 IHC 28-8 pharmDx Control Slides

EnVision FLEX Wash Buffer (20x) (Code K8007) and EnVision FLEX Hematoxylin (Code K8008), are required but not included in the kit. Refer to the IFU for a complete list of required materials and equipment.

# Technical Considerations for Optimal Performance of PD-L1 IHC 28-8 pharmDx

Optimal staining performance is achieved by adhering to the PD-L1 IHC 28-8 pharmDx protocol. Technical problems relating to the performance of PD-L1 IHC 28-8 pharmDx may arise in two areas; those involving specimen collection and specimen preparation prior to performing the test, as well as problems involving the actual performance of the test itself. Technical problems related to the performance of the test are generally related to procedural deviations and can be controlled and minimized through training and a thorough understanding of the product instructions by the user.

## Specimen Collection and Preparation

Specimens must be handled in a way which preserves the tissue for immunohistochemical staining. Tissue should be stained and interpreted as close to the time of biopsy as possible. The stability of PD-L1 immunoreactivity in tissue blocks has not been assessed. Tissue may be susceptible to loss of PD-L1 immunoreactivity with age. Confirm appropriate intact tumor morphology and the presence of sufficient tumor cells for evaluation. Use recommended methods of tissue processing for all specimens.

## Tissue Processing

FFPE tissues are suitable for use. Specimens should be blocked into a thickness of 3 mm or 4 mm, fixed in 10% Neutral Buffered Formalin (NBF), and dehydrated and cleared in a series of alcohols and xylene, followed by infiltration with melted paraffin. The paraffin temperature should not exceed 60 °C. An ischemia time from excision to fixation start time of less than 30 minutes followed by immersion in NBF for 24–48 hours is recommended.

The use of PD-L1 IHC 28-8 pharmDx on decalcified tissue has not been validated and is not recommended.

Cut tissue specimens into sections of 4–5 µm. After sectioning, mount tissues on FLEX IHC microscope slides, Code K8020, or Superfrost Plus charged slides and place in a 58 ± 2 °C oven for 1 hour.

To preserve antigenicity, UC and MIUC tissue sections, once mounted on slides, should be stored in the dark at 2–8 °C, or room temperature up to 25 °C, and stained within 4 months of sectioning. Slide storage and handling conditions should not exceed 25 °C at any point post-mounting to ensure tissue integrity and antigenicity.



## Control Tissues

Differences in processing and embedding in the user's laboratory may produce significant variability in results. Include positive and negative control tissue in each staining run, in addition to the PD-L1 IHC 28-8 pharmDx Control Slide.

Control tissue should be biopsy/surgical specimens, fixed, processed, and embedded as soon as possible in the same manner as the patient sample(s). Control tissue must represent one of the approved tumor indications for PD-L1 IHC 28-8 pharmDx as listed in the Intended Use of the IFU. Control tissues processed differently from the patient specimen validate reagent performance only and do not verify tissue preparation. The ideal positive control tissue provides a complete dynamic representation of weak-to-moderate staining of tumor cells. However, prevalence of PD-L1 expression of immune cells is high, therefore a few staining immune cells is acceptable. Alternatively, negative portions of the positive control tissue may serve as the negative control tissue, but this should be verified by the user.

## PD-L1 IHC 28-8 pharmDx Staining Procedure

The PD-L1 IHC 28-8 pharmDx reagents and instructions have been designed for optimal performance. Further dilution of the reagents, alteration of incubation times, temperatures, or instruments may give erroneous results.

## Reagent Storage

Store all components of PD-L1 IHC 28-8 pharmDx, including Control Slides, in the dark at 2–8 °C when not in use on Autostainer Link 48. Do not use after the expiration date printed on the outside package.

## Reagent Preparation

Equilibrate all components to room temperature (20–25 °C) prior to immunostaining.

### EnVision FLEX Target Retrieval Solution, Low pH

Prepare a sufficient quantity of EnVision FLEX 1x Target Retrieval Solution, Low pH by diluting EnVision FLEX Target Retrieval Solution, Low pH (50x) 1:50 using distilled or de-ionized water; the pH of EnVision FLEX 1x Target Retrieval Solution should be  $6.1 \pm 0.2$ . Do not modify the pH of 1x EnVision FLEX Target Retrieval Solution after preparation under any circumstance. If a problem is suspected with the EnVision FLEX Target Retrieval Solution pH, please refer to the troubleshooting section for more information. One 30 mL bottle of Target Retrieval Solution, Low pH (50x), diluted 1:50 will provide 1.5 L of EnVision FLEX 1x reagent, sufficient to fill one PT Link tank, which will treat up to 24 slides per use. Discard EnVision FLEX 1x Target Retrieval Solution after three uses and no longer than 5 days after dilution. Note the EnVision FLEX Target Retrieval Solution, low pH (50x) is a red-colored solution.

Additional EnVision FLEX Target Retrieval Solution, Low pH (50x), if required, is available as Code K8005.

**EnVision FLEX Wash Buffer (20x)**

Prepare a sufficient quantity of 1x EnVision FLEX Wash Buffer by diluting Wash Buffer (20x) 1:20 using distilled or de-ionized water and mix thoroughly. Store unused 1x solution at 2–8 °C for no more than one month. Discard buffer if cloudy in appearance. Refer to the User Guide for your Autostainer Link 48 for further information. EnVision FLEX Wash Buffer (20x) is available as Code K8007.

**DAB+ Substrate-Chromogen Solution**

Add 9 drops of DAB+ Chromogen per 1 bottle of DAB+ Substrate Buffer and mix. Prepared DAB+ Substrate-Chromogen is stable for 5 days if stored in the dark at 2–8 °C. Mix the DAB+ Substrate-Chromogen Solution thoroughly prior to use. Any precipitate developing in the solution does not affect staining quality.

- When mixing DAB+ Substrate Buffer with DAB+ Chromogen, add 9 drops of DAB+ Chromogen to the bottle of DAB+ Substrate Buffer and mix.
- The color of the DAB+ Chromogen may vary from clear to lavender-brown. This will not affect the performance of the product. Dilute as per the guidelines above. Adding excess DAB+ Chromogen to the DAB+ Substrate Buffer results in deterioration of the positive signal.

**Control to Assess Staining**

Each slide contains sections of two pelleted, FFPE cell lines: NCI-H226\*\* with positive PD-L1 protein expression (originating from human lung squamous cell carcinoma and MCF-7 with negative PD-L1 protein expression (originating from human breast adenocarcinoma).

\*\* Dr. AF Gazdar and Dr. JD Minna at NIH are acknowledged for their contribution in developing NCI-H226 ATCC Number: CRL-5826™.

**Staining Protocol**

Program slides by selecting PD-L1 IHC 28-8 pharmDx staining protocol from the options in the DakoLink drop-down menu. All of the required steps and incubation times for staining are preprogrammed in the DakoLink software. Print and attach slide labels to each slide.

**Deparaffinization, Rehydration, and Target Retrieval**

Use PT Link, Code PT100/PT101/PT200, to perform the Deparaffinization, Rehydration, and Target Retrieval 3-in-1 procedure.

- Set Preheat and Cool to 65 °C, and set Heat to 97 °C for 20 minutes.
- Fill PT Link tanks with 1.5 L per tank of EnVision FLEX Target Retrieval Solution, Low pH, working solution to cover the tissue sections.
- Preheat the Target Retrieval Solution, Low pH to 65 °C.
- Immerse Autostainer racks containing mounted, FFPE tissue sections into the preheated Target Retrieval Solution, Low pH (1x working solution) in PT Link tank. Incubate for 20 minutes at 97 °C.

- As soon as the target retrieval incubation time has been completed, and the temperature has cooled to 65 °C, remove each Autostainer slide rack with slides from the PT Link tank and **immediately** place the slides into a tank (e.g., PT Link Rinse Station, Code PT109) containing room temperature EnVision FLEX Wash Buffer 1x working solution.
- Leave Autostainer rack with slides in room temperature EnVision FLEX Wash Buffer for 5 minutes.

### Staining and Counterstaining

- Place the Autostainer rack with slides on the Autostainer Link 48. Ensure slides remain wet with buffer while loading and prior to initiating the run. Dried tissue sections may display increased nonspecific staining.
- The instrument performs the staining and counterstaining procedures by applying the appropriate reagent, monitoring the incubation time, and rinsing slides between reagents. Counterstaining using EnVision FLEX Hematoxylin, (Code K8008), for 7 minutes, is included in the staining protocol. Do not allow tissue sections to dry after deparaffinization, rehydration, and target retrieval (3-in-1) procedure, or at any time during or after the staining run prior to permanent mounting.

### Mounting

Use nonaqueous permanent mounting media. To minimize fading, store slides in the dark at room temperature (20–25 °C).



# PD-L1 IHC 28-8 pharmDx Technical Checklist

Customer Name/Institution: \_\_\_\_\_

Name and Title: \_\_\_\_\_

Autostainer Link 48 Serial Number: \_\_\_\_\_ Software Version: \_\_\_\_\_

	Yes	No
1. Regular preventive maintenance is performed on the Autostainer Link 48 and PT Link?	<input type="checkbox"/>	<input type="checkbox"/>
2. PD-L1 IHC 28-8 pharmDx is used before the expiration date printed on the outside of the box?	<input type="checkbox"/>	<input type="checkbox"/>
3. All PD-L1 IHC 28-8 pharmDx components, including Control Slides, are stored in the dark at 2–8 °C?	<input type="checkbox"/>	<input type="checkbox"/>
4. All PD-L1 IHC 28-8 pharmDx components, including Control Slides, are equilibrated to room temperature (20–25 °C) prior to immunostaining?	<input type="checkbox"/>	<input type="checkbox"/>
5. Appropriate positive and negative control tissues are identified?	<input type="checkbox"/>	<input type="checkbox"/>
6. Tissues are fixed in 10% neutral buffered formalin?	<input type="checkbox"/>	<input type="checkbox"/>
7. Tissues are infiltrated with melted paraffin, at or below 60 °C?	<input type="checkbox"/>	<input type="checkbox"/>
8. Tissue sections of 4–5 µm are mounted on FLEX IHC Microscope Slides, or Superfrost Plus charged slides?	<input type="checkbox"/>	<input type="checkbox"/>
9. UC and MIUC specimens are stained within 4 months of sectioning when stored in the dark at 2–8 °C or at room temperature up to 25 °C?	<input type="checkbox"/>	<input type="checkbox"/>
10. EnVision FLEX Target Retrieval Solution, Low pH is prepared properly (working solution pH 6.1±0.2)?	<input type="checkbox"/>	<input type="checkbox"/>
11. EnVision FLEX Wash Buffer is prepared properly?	<input type="checkbox"/>	<input type="checkbox"/>
12. DAB+ Substrate-Chromogen Solution is mixed and prepared properly?	<input type="checkbox"/>	<input type="checkbox"/>
13. The Deparaffinization, Rehydration, and Target Retrieval 3-in-1 procedure is followed, using PT Link?	<input type="checkbox"/>	<input type="checkbox"/>
14. Slides remain wet with buffer while loading and prior to initiating run on Autostainer Link 48?	<input type="checkbox"/>	<input type="checkbox"/>
15. The PD-L1 IHC 28-8 pharmDx protocol is selected on Autostainer Link 48?	<input type="checkbox"/>	<input type="checkbox"/>
16. Slides are counterstained with EnVision FLEX Hematoxylin?	<input type="checkbox"/>	<input type="checkbox"/>
17. Do you have all the necessary equipment to perform the PD-L1 IHC 28-8 pharmDx test according to the protocol? If not, specify what is missing in the comments below.	<input type="checkbox"/>	<input type="checkbox"/>

If you answered "No" to any of the above, consult with your local Agilent Technical Support Representative for assistance.

Additional Observations or Comments:

---



---



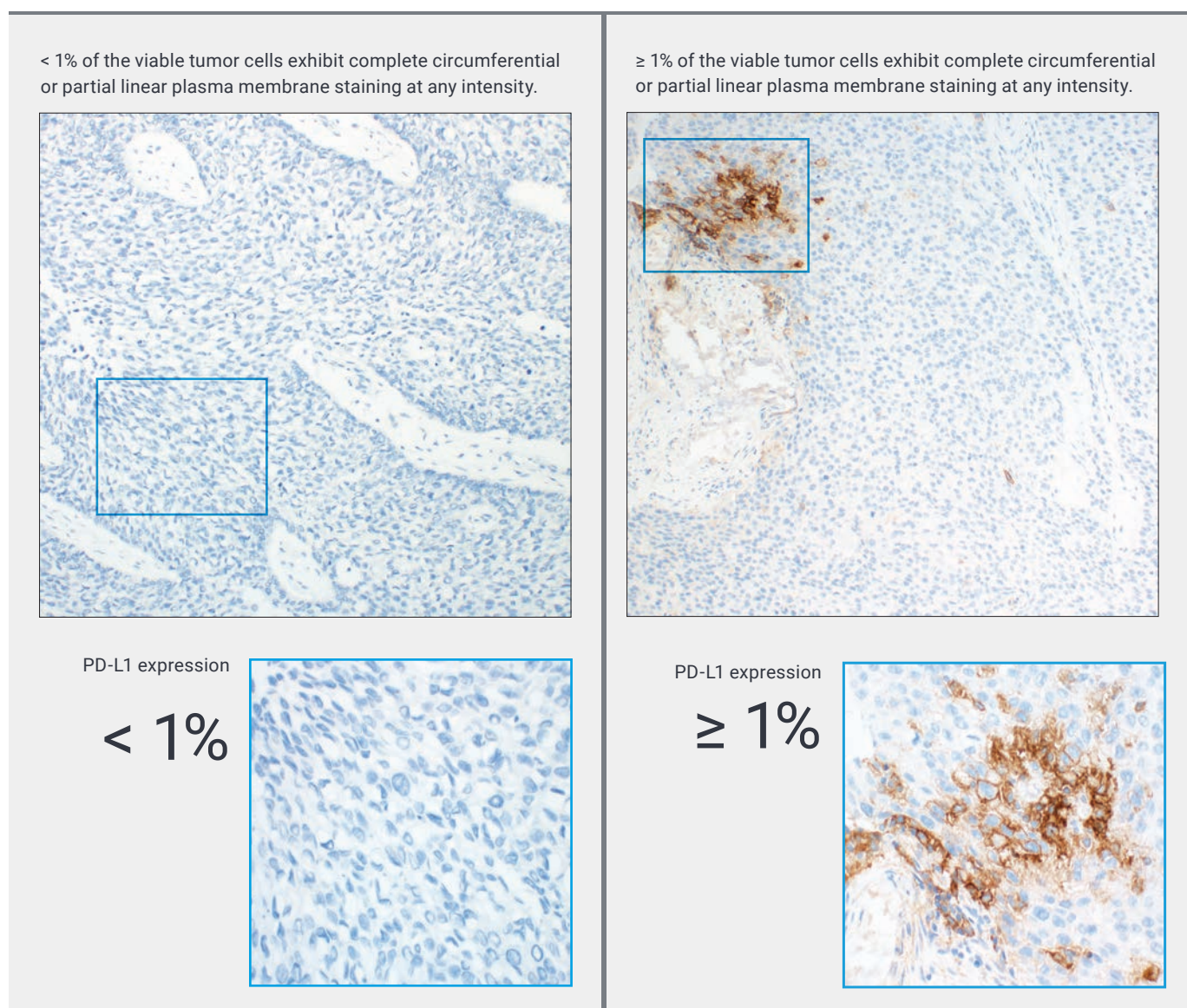
---

# Guidelines for Scoring PD-L1 IHC 28-8 pharmDx in UC and MIUC

Agilent emphasizes that scoring of PD-L1 IHC 28-8 pharmDx must be performed in accordance with the guidelines established in the IFU, within the context of best practices and the pathologist's experience.

This assay was validated for invasive UC including MIUC tissue samples and not for lesions with foci of dysplasia or carcinoma in situ. Hematoxylin and eosin (H&E) stained slides should accompany each PD-L1 stained sample to allow proper assessment of invasive carcinoma, carcinoma in situ, and adjacent normal epithelium.

The percentage of stained viable tumor cells in the specimen determines PD-L1 IHC 28-8 pharmDx result. Scoring guidelines and reporting recommendations are presented in Figure 3. See page 23-24 for an example of a reporting results form for PD-L1 IHC 28-8 pharmDx.



**Figure 3.** Guidelines for scoring and reporting PD-L1 IHC pharmDx results.

# Slide Evaluation Flowchart

The following flow of slide review is recommended when conducting interpretation of PD-L1 IHC 28-8 pharmDx. Refer to the detailed description on pages 18–22.

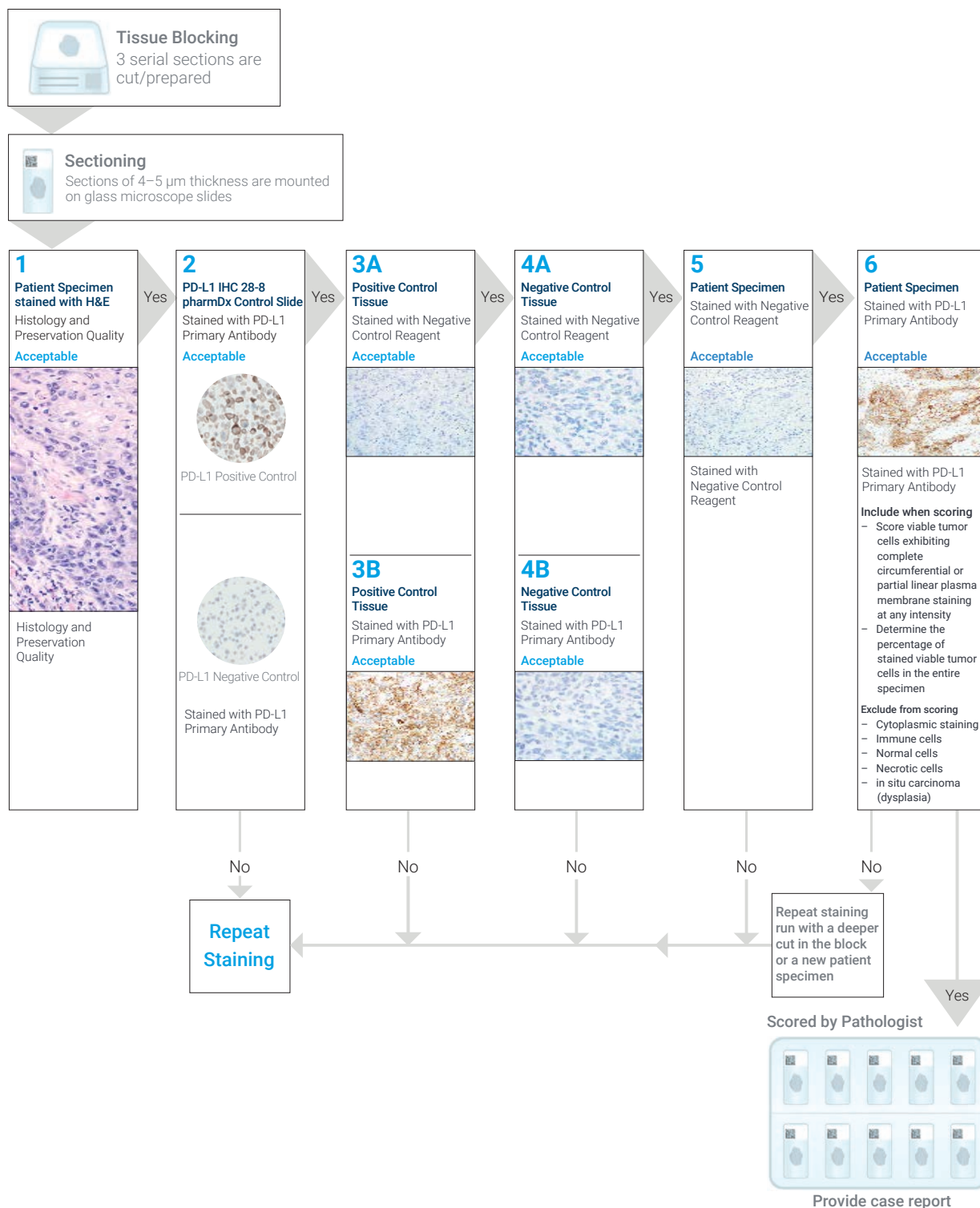


Figure 4. Slide evaluation procedure steps.



# Recommendations for Interpretation of PD-L1 IHC 28-8 pharmDx in UC and MIUC

PD-L1 IHC 28-8 pharmDx evaluation should be performed by a pathologist using a bright field microscope. Before examining the patient specimen for PD-L1 staining, it is important to examine the hematoxylin and eosin (H&E) and controls first, to assess staining quality. Examine a serial section of the patient specimen stained with H&E for histology and preservation quality. Then, examine the PD-L1 IHC 28-8 pharmDx Control Slide, the positive and negative control tissue slides, and the patient specimen slide stained with the Negative Control Reagent (NCR) to assess reagent performance.

Lastly, examine the patient specimen stained with Primary Antibody to assess the staining of viable tumor cells.

PD-L1 staining is defined as complete circumferential or partial linear plasma membrane staining at any intensity. Cytoplasmic staining, if present, is not included in the score. Nonmalignant cells and immune cells (e.g., such as infiltrating lymphocytes or macrophages) may also stain with PD-L1; however, these should not be included in the scoring for the determination of PD-L1 % tumor cell expression.

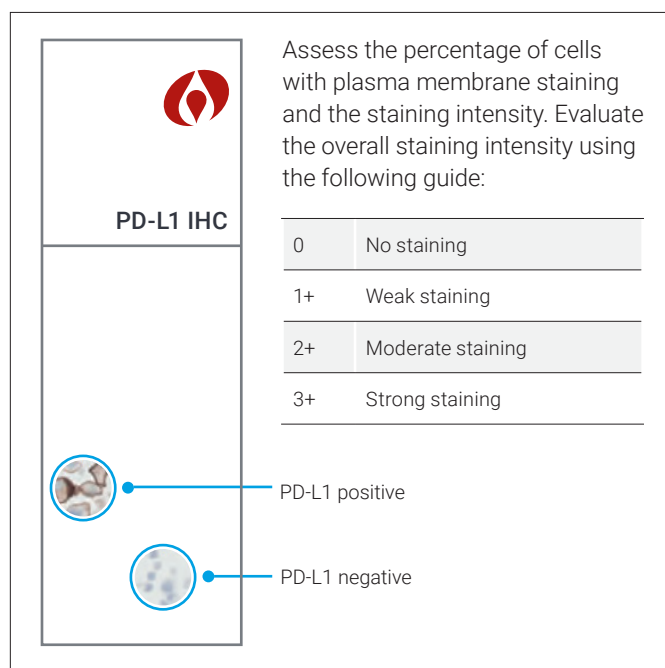
For the determination of PD-L1 % tumor cell expression positive control tissue slides and negative control tissue slides should be supplied by the laboratory. Only the PD-L1 IHC pharmDx Control Cell Line Slide is provided in the kit.

## Patient Specimen Stained with H&E

An hematoxylin & eosin (H&E) stained section is required for the evaluation of histology and preservation quality. PD-L1 IHC 28-8 pharmDx and the H&E staining should be performed on a serial section from the same paraffin block of the specimen.

## PD-L1 IHC 28-8 pharmDx Control Slide

Examine the PD-L1 IHC 28-8 pharmDx Control Slide to ascertain that reagents are functioning properly. Each slide contains sections of cell pellets with positive and negative PD-L1 expression, see Figure 5. The 10x objective is recommended for the evaluation of the cell pellets as a whole. If any staining of the Control Cell Line is not satisfactory, all results with the patient specimens should be considered invalid. For a borderline Control Slide, run outcome should be considered in the context of all run controls, including control tissue. Do not use the Control Slide as an aid in the interpretation of patient results.



**Figure 5.** Each Control Slide contains sections of cell pellets with positive and negative PD-L1 expression.

For the PD-L1 positive cell pellet on the Control Slide, the following staining is acceptable, see Figure 6:

- Plasma membrane staining of  $\geq 80\%$  of cells
- $\geq 2+$  average staining intensity of cells with membrane staining
- Nonspecific staining  $< 1+$  intensity

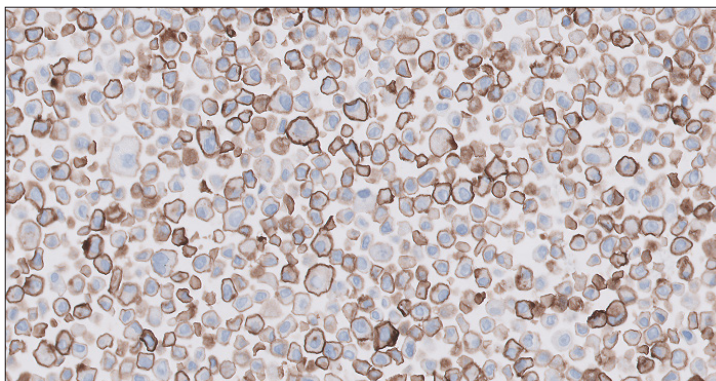


Figure 6. Acceptable Positive PD-L1 Control.

For the PD-L1 negative cell pellet on the Control Slide, the following staining is acceptable, see Figure 7:

- No specific staining
- Nonspecific staining  $< 1+$  intensity

Staining of a few cells in the negative pellet on the Control Slide may occasionally be observed. The presence of  $\leq 10$  cells with distinct plasma membrane staining, and/or cytoplasmic staining with  $\geq 1+$  intensity within the boundaries of the negative cell pellet is acceptable.

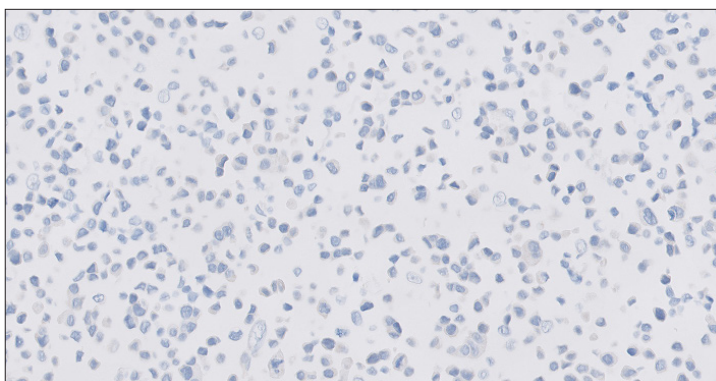


Figure 7. Acceptable Negative PD-L1 Control.

## Positive Control Tissue Slides

Examine the positive control tissue slides

(Primary Antibody, NCR) to ascertain if tissues are correctly prepared, and reagents are functioning properly. Any nonspecific staining should be of  $\leq 1+$  staining intensity. Exclude necrotic or nonviable tumor cells from the evaluation. If the staining of positive control tissues is not satisfactory, all results with the patient specimens should be considered invalid. Do not use control tissue as an aid in the interpretation of patient results.

## Negative Control Tissue Slides

Examine the negative control tissue slides

(Primary Antibody, NCR) to confirm that there is no unintended staining. Any nonspecific staining should be  $\leq 1+$  staining intensity. If plasma membrane staining of malignant cells occurs in the negative control tissue, all results with the patient specimens should be considered invalid. Do not use control tissue as an aid in the interpretation of patient results.

## Patient Specimen Stained with Negative Control Reagent

Examine the patient specimen stained with NCR to ascertain that reagents are functioning properly. The absence of plasma membrane staining of viable tumor cells is satisfactory and nonspecific staining should be  $\leq 1+$  staining intensity. If any staining is not satisfactory, results with the patient specimen should be considered invalid.

The NCR indicates nonspecific staining and allows better interpretation of the patient specimen stained with the Primary Antibody.

## Patient Specimen Stained with Primary Antibody

Staining should be assessed within the context of any nonspecific staining of the patient specimen stained with NCR. A minimum of 100 viable tumor cells must be present in the PD-L1 stained patient slide in order to perform an evaluation.

1	At 4x objective magnification, carefully examine the tumor areas of the entire specimen. All areas with viable tumor cells on the specimen should be evaluated. Exclude nonmalignant cells, necrotic cells, and cellular debris. Cytoplasmic staining, if present, should be disregarded.
2	Use the 10–20x objective magnifications to determine the percentage of viable tumor cells expressing PD-L1 membranous staining. The 40x objective can be used for confirmation if needed. Tumor cells are considered to be PD-L1 positive if they exhibit either partial linear or complete circumferential staining of the plasma membrane at any intensity. Nonmalignant cells and immune cells (e.g., infiltrating lymphocytes or macrophages) may also stain with PD-L1 but must be excluded.
3	Record if the specimen has PD-L1 % tumor cell expression < 1% or ≥ 1%. When determining the percentage of stained tumor cells in the entire specimen, the numerator is the number of stained viable tumor cells and the denominator is the total number of viable tumor cells in the specimen.

$$\% \text{ PD-L1 expression} = \frac{\# \text{ PD-L1 staining tumor cells}}{\text{Total \# viable tumor cells}} \times 100$$

## Tips and Special Considerations

- Include the entire specimen for evaluation of PD-L1 % tumor cell expression
- Use higher magnifications to confirm cell types and areas absent of staining
- Be careful not to overlook weak 1+ staining, which can be missed at 4x and 10x
- Disregard cytoplasmic staining
- Necrotic tissue may stain but should be excluded
- Exclude any nonmalignant cells and immune cells
- Granular staining must demonstrate a perceptible and convincing membrane pattern

## Non-evaluable Specimen

The specimen should be considered non-evaluable if there are fewer than 100 viable tumor cells. A different section from the same block or another block from the same patient may be required to present a sufficient quantity of viable tumor cells for PD-L1 IHC 28-8 pharmDx evaluation.

## Indeterminate Specimen

The tumor cell membrane has been hampered for reasons attributed to the biology of the tumor tissue sample rather than improper sample preparation. For example, high cytoplasmic staining of the tumor cells can hamper the scoring of the membrane staining. An additional cut section or section from another block of the same patient may be required for PD-L1 IHC 28-8 pharmDx evaluation.



# PD-L1 IHC 28-8 pharmDx Suggested Scoring Methods for Calculating PD-L1 Percent Tumor Cell Expression

Agilent offers two different examples of scoring techniques that may be used when assessing stained specimens exhibiting different staining patterns.

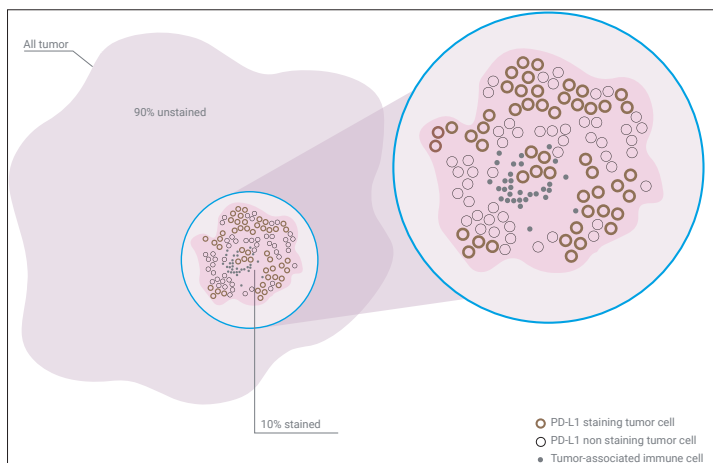
## Example 1: Calculating PD-L1 % tumor cell expression in a specimen with a small PD-L1 staining tumor area

At lower objective magnification, assess the entire specimen for the presence of PD-L1 staining in viable tumor cells at any intensity. Any nonmalignant and immune cells staining PD-L1 positive must be excluded.

- In this example, assume the number of tumor cells is equally distributed in the tumor and that there are a total of 1,000 viable tumor cells in the entire specimen.
- 10% of the tumor area has staining, 90% of the tumor area has no staining.

At a higher objective magnification, carefully examine PD-L1 staining tumor area (blue circle in Figure 8). PD-L1 positive staining is defined as complete circumferential and/or partial linear plasma membrane staining of tumor cells at any intensity.

- 50 out of 100 viable tumor cells are staining PD-L1 positive in the single region of the tumor area (blue circle in Figure 8).
- 50 out of 1000 viable tumor cells are staining PD-L1 positive in the entire tumor specimen (Method 1) which may also be described as 50% PD-L1 positive in a single region representing 10% of the tumor area (Method 2).



**Figure 8.** Example of a tumor with a small PD-L1 staining area.

Determine the overall percentage of PD-L1 staining tumor cells for the entire specimen as shown:

### Method 1

$$\frac{50 \text{ tumor cells staining PD-L1 positive}}{1,000 \text{ viable tumor cells}} \times 100 = 5\% \text{ tumor cell expression}$$

### Method 2

$$\frac{50\% \times 10\%}{100} = 5\% \text{ tumor cell expression}$$

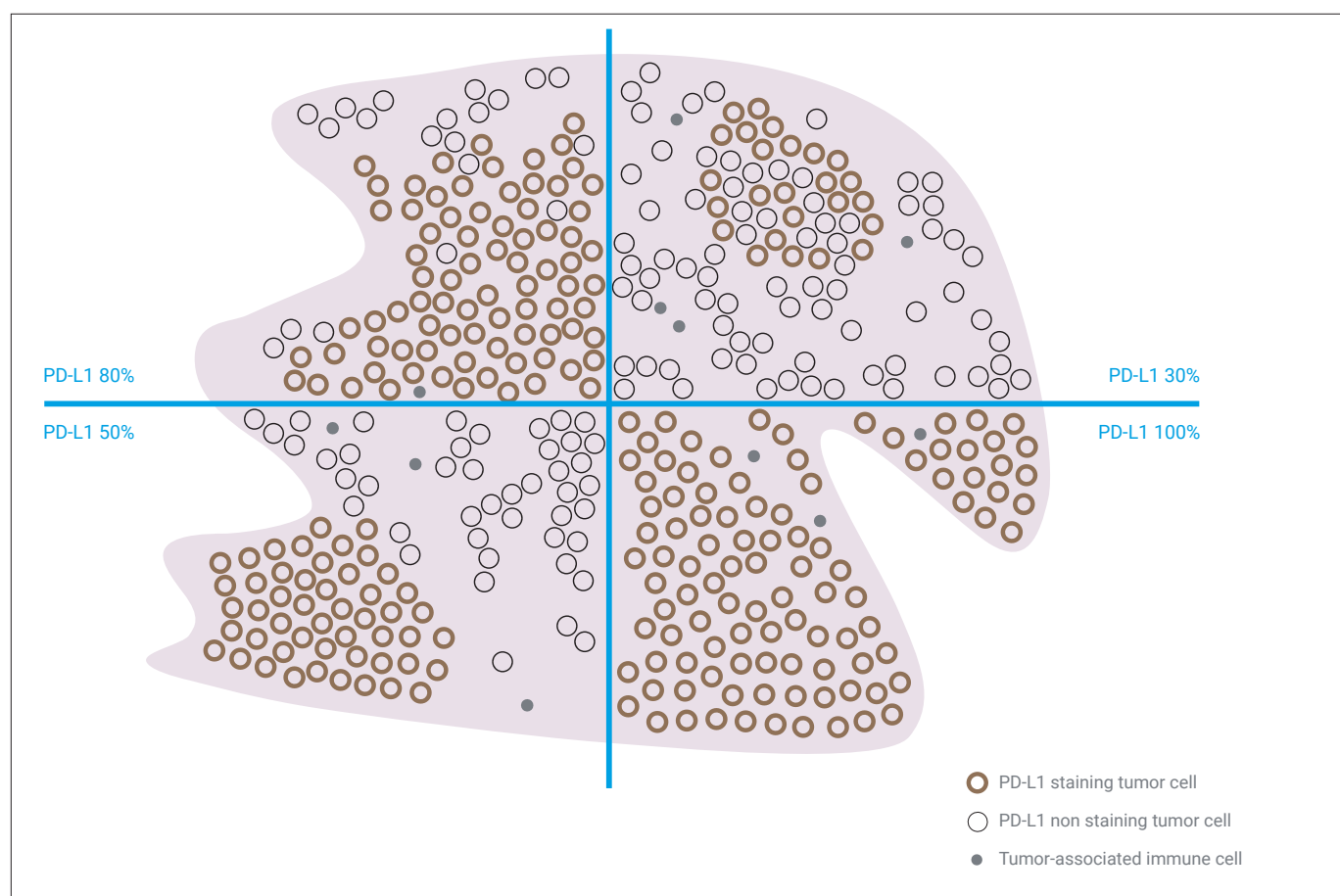
**Example 2: Calculating PD-L1 % tumor cell expression in a specimen with heterogeneous staining**

At lower objective magnification, assess the entire specimen for the presence of PD-L1 staining in viable tumor cells at any intensity. Visually divide the tumor area into regions. Any nonmalignant and immune cells staining PD-L1 positive must be excluded.

- The tumor area is divided into four equivalent quadrants in Figure 9.

At a higher objective magnification, assess and calculate the percentage of PD-L1 staining tumor cells in each quadrant. PD-L1 positive staining is defined as complete circumferential and/or partial linear plasma membrane staining of tumor cells at any intensity.

- The percentage of PD-L1 staining tumor cells for each of the four respective quadrants are 80%, 30%, 50% and 100%.



**Figure 9.** Example with heterogenous PD-L1 staining area.

Determine the overall percentage of PD-L1 staining tumor cells for the entire specimen:

$$\frac{(80\% + 30\% + 50\% + 100\%)}{4 \text{ quadrants}} = 65\% \text{ tumor cell expression}$$

## Reporting Results - UC

### Suggested information to include when reporting results with PD-L1 IHC 28-8 pharmDx in UC

#### PD-L1 IHC 28-8 pharmDx, Code SK005

##### Summary of Sample Tested:

Date of Run: \_\_\_\_\_ PD-L1 IHC 28-8 pharmDx Lot: \_\_\_\_\_

Staining Run Log ID: \_\_\_\_\_ Specimen ID: \_\_\_\_\_

Patient Identifier: \_\_\_\_\_

Type of Service: IHC Stain with Manual Interpretation

Other: \_\_\_\_\_

Type of Tissue: \_\_\_\_\_

Additional Tests Performed with PD-L1 IHC 28-8 pharmDx: \_\_\_\_\_

##### PD-L1 IHC 28-8 pharmDx Controls Results:

PD-L1 IHC 28-8 Control Slide: Pass ☐ Fail ☐

Positive Control Tissue Slides: Pass ☐ Fail ☐

Negative Control Tissue Slides: Pass ☐ Fail ☐

Patient Specimen, Negative Control Reagent: Pass ☐ Fail ☐

**PD-L1 Results:** Viable Tumor Cells Present: ☐  $\geq 100$  cells ☐ Not evaluable

☐ PD-L1 % tumor cell expression is  $< 1\%$

Percent of UC cells with complete circumferential and/or partial linear membrane PD-L1 staining is  $< 1\%$ .

☐ PD-L1 % tumor cell expression is  $\geq 1\%$

Percent of UC cells with complete circumferential and/or partial linear membrane PD-L1 staining is  $\geq 1\%$ .

Other Comments to Treating Physician: \_\_\_\_\_

*Note: PD-L1 IHC 28-8 pharmDx was validated for invasive UC tissue samples and not for lesions with foci of dysplasia or carcinoma in situ. An H&E stained slide should accompany each PD-L1 stained sample to allow a proper assessment of invasive carcinoma, carcinoma in situ, and adjacent normal epithelium.*

## Reporting Results - MIUC

### Suggested information to include when reporting results with PD-L1 IHC 28-8 pharmDx in MIUC

#### PD-L1 IHC 28-8 pharmDx, Code SK005

##### Summary of Sample Tested:

Date of Run: \_\_\_\_\_ PD-L1 IHC 28-8 pharmDx Lot: \_\_\_\_\_

Staining Run Log ID: \_\_\_\_\_ Specimen ID: \_\_\_\_\_

Patient Identifier: \_\_\_\_\_

Type of Service: IHC Stain with Manual Interpretation

Other: \_\_\_\_\_

Type of Tissue: \_\_\_\_\_

Additional Tests Performed with PD-L1 IHC 28-8 pharmDx: \_\_\_\_\_

##### PD-L1 IHC 28-8 pharmDx Controls Results:

PD-L1 IHC 28-8 Control Slide: Pass ☐ Fail ☐

Positive Control Tissue Slides: Pass ☐ Fail ☐

Negative Control Tissue Slides: Pass ☐ Fail ☐

Patient Specimen, Negative Control Reagent: Pass ☐ Fail ☐

**PD-L1 Results:** Viable Tumor Cells Present: ☐  $\geq 100$  cells ☐ Not evaluable

☐ PD-L1 % tumor cell expression is  $< 1\%$

Percent of MIUC cells with complete circumferential and/or partial linear membrane PD-L1 staining is  $< 1\%$ .

☐ PD-L1 % tumor cell expression is  $\geq 1\%$

Percent of MIUC cells with complete circumferential and/or partial linear membrane PD-L1 staining is  $\geq 1\%$ .

Other Comments to Treating Physician: \_\_\_\_\_

*Note: PD-L1 IHC 28-8 pharmDx was validated for invasive MIUC tissue samples and not for lesions with foci of dysplasia or carcinoma in situ. An H&E stained slide should accompany each PD-L1 stained sample to allow a proper assessment of invasive carcinoma, carcinoma in situ, and adjacent normal epithelium.*



# PD-L1 IHC 28-8 pharmDx

## Immunostaining Examples in UC and MIUC

The following images present examples of UC and MIUC tumor samples stained with PD-L1 IHC 28-8 pharmDx.

An example of urothelial carcinoma of the bladder stained with PD-L1 IHC 28-8 pharmDx. The staining shows a range of PD-L1 expression. This specimen would be appropriate to use as a positive control specimen for detection of subtle changes in assay sensitivity. Note the partial linear (red arrow) and complete circumferential (black arrow) plasma membrane staining.

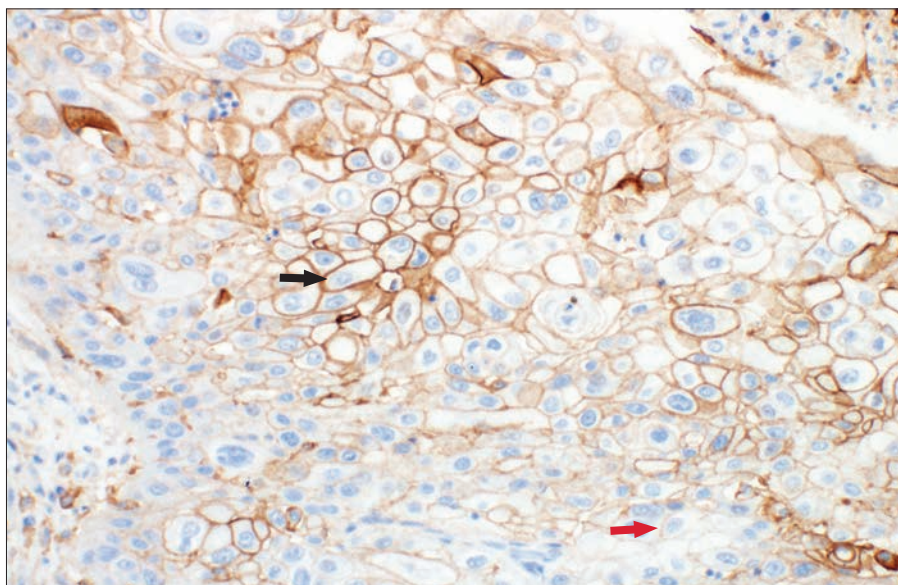


Figure 10. 20x magnification.

Urothelial carcinoma of the bladder  
PD-L1 % tumor cell expression 0%.

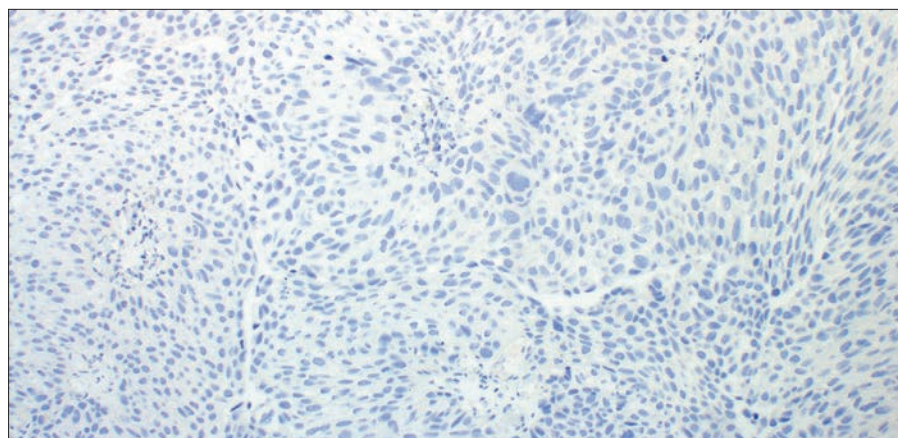
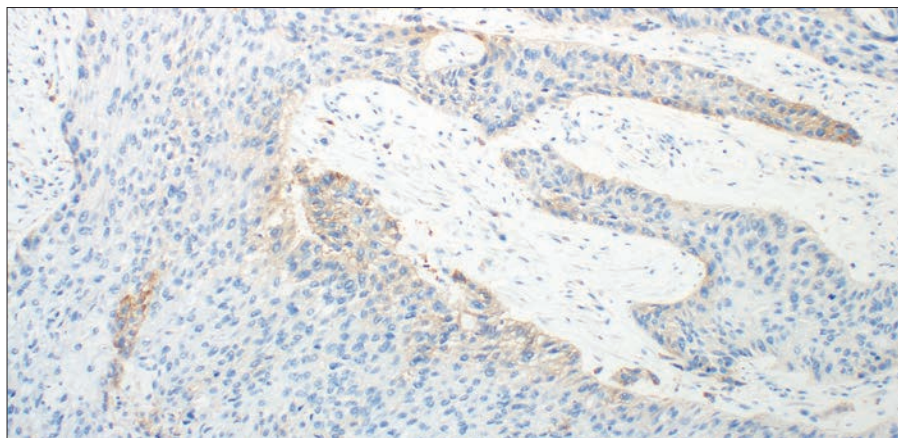


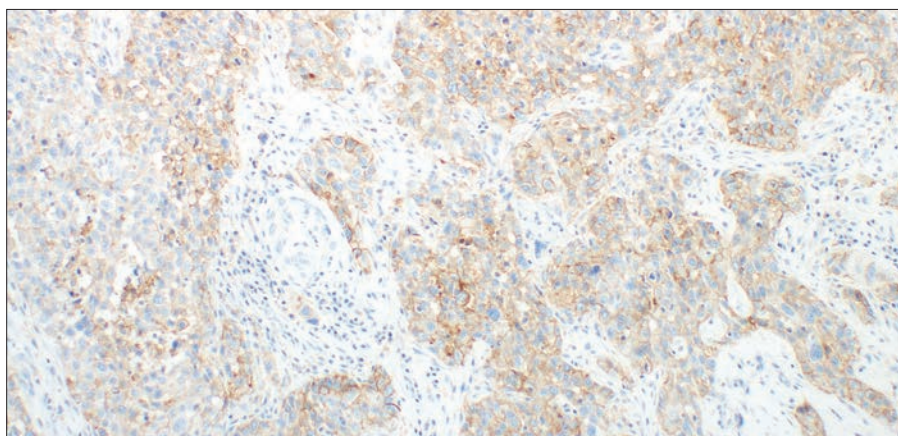
Figure 11. 10x magnification.

Urothelial carcinoma of the bladder  
PD-L1 % tumor cell expression  $\geq 1\%$ .



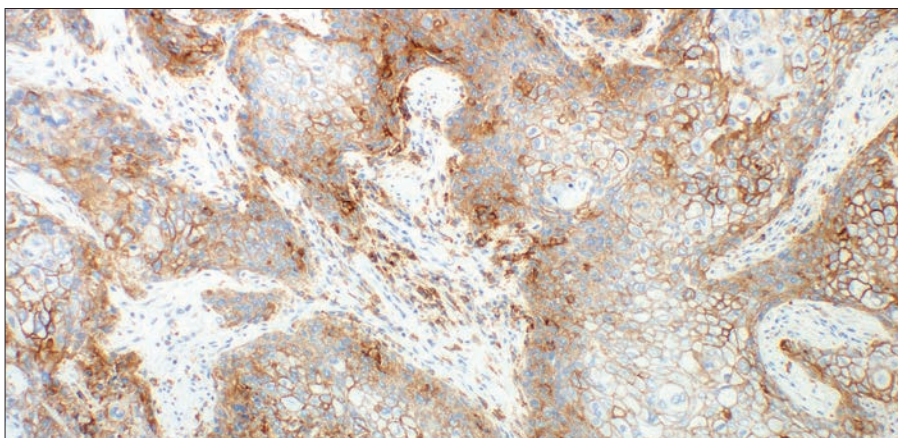
**Figure 12.** 10x magnification.

Urothelial carcinoma of the bladder  
demonstrating  $\geq 1\%$ , moderate  
PD-L1 % tumor cell expression.



**Figure 13.** 10x magnification.

Urothelial carcinoma of the bladder  
demonstrating  $\geq 1\%$ , high PD-L1 % tumor  
cell expression.



**Figure 14.** 10x magnification.



Urothelial carcinoma of the bladder showing strong staining of intra-tumoral immune cells (**red arrows**), while the tumor cells are negative (**black arrows**) for PD-L1. Note the staining of intra-tumoral mononuclear inflammatory cells (histiocytes and lymphocytes) are not included in determining the PD-L1 % tumor cell expression.

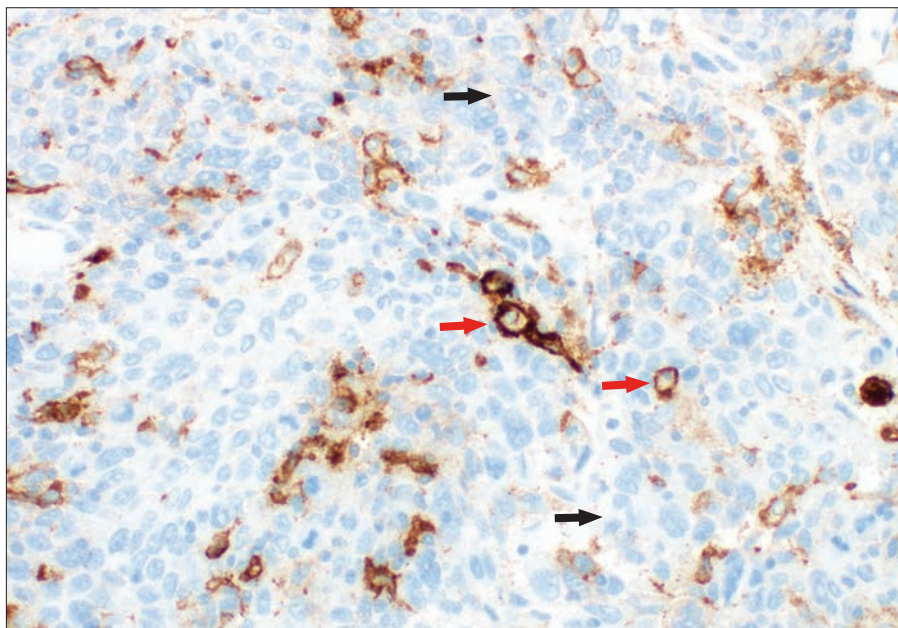


Figure 15. 20x magnification.

Urothelial carcinoma of the bladder showing PD-L1 staining of peri-tumoral immune cells (**red arrows**) and tumor cells (**black arrows**). Note the staining of peri-tumoral mononuclear inflammatory cells (histiocytes and lymphocytes) are not included in determining the PD-L1 % tumor cell expression.

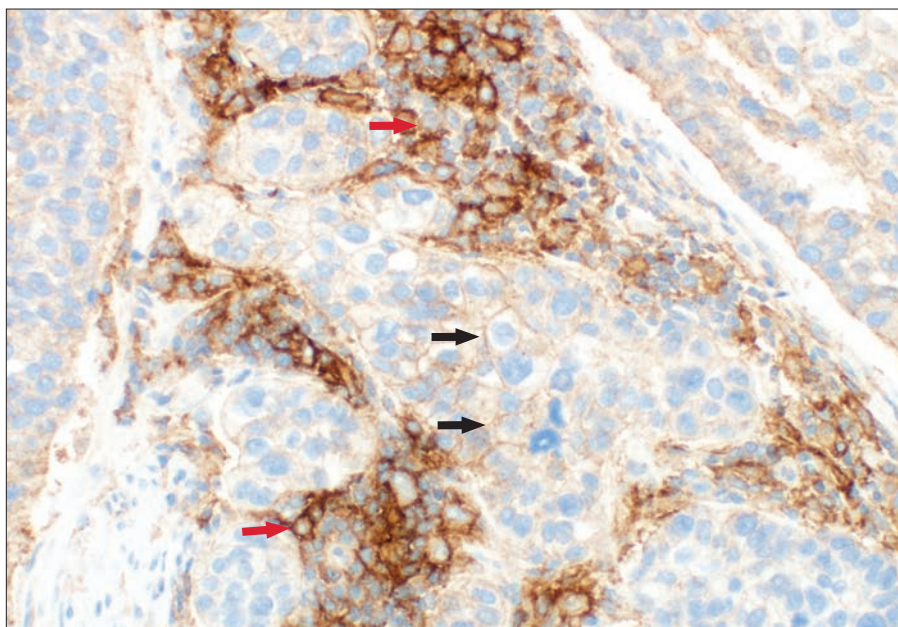


Figure 16. 20x magnification.

H&E stain case of UC of the bladder showing **in situ** component (**red arrow**).

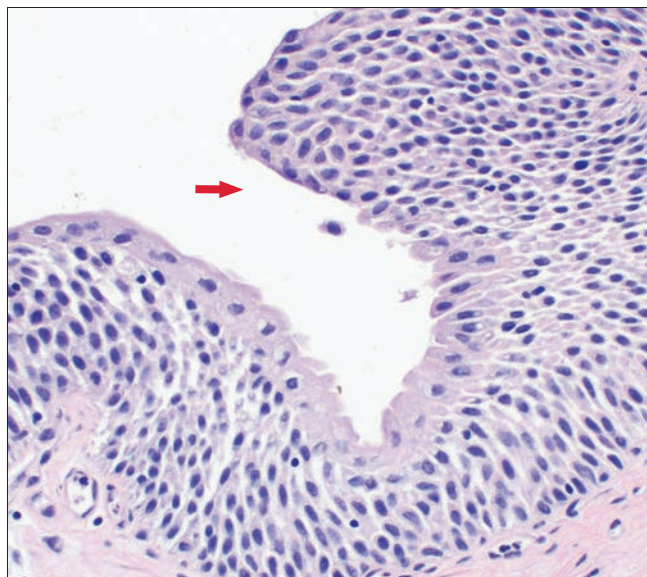


Figure 17a. 20x magnification.

Urothelial carcinoma of the bladder showing negative staining for PD-L1 in the dysplasia/in situ component (**red arrow**). When scoring PD-L1 percent tumor cell expression, the in situ component is not included in the denominator. Only the invasive component is evaluated when determining percent tumor cell expression.

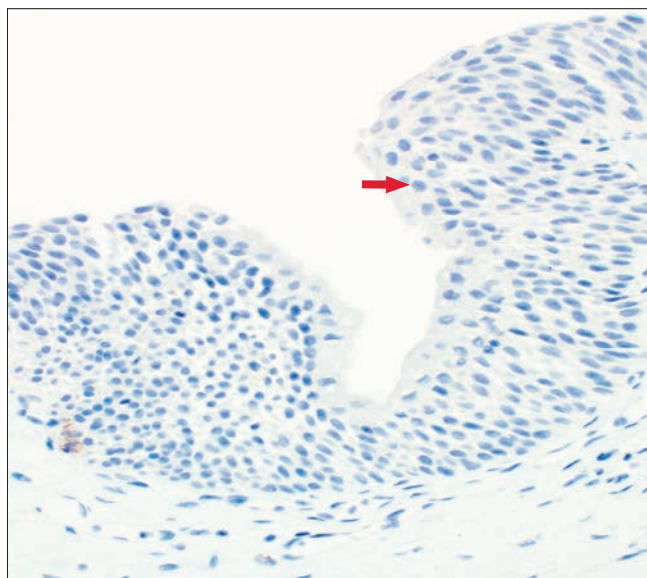


Figure 17b. 20x magnification.

H&E stain of UC of the bladder showing normal urothelium (red arrow).

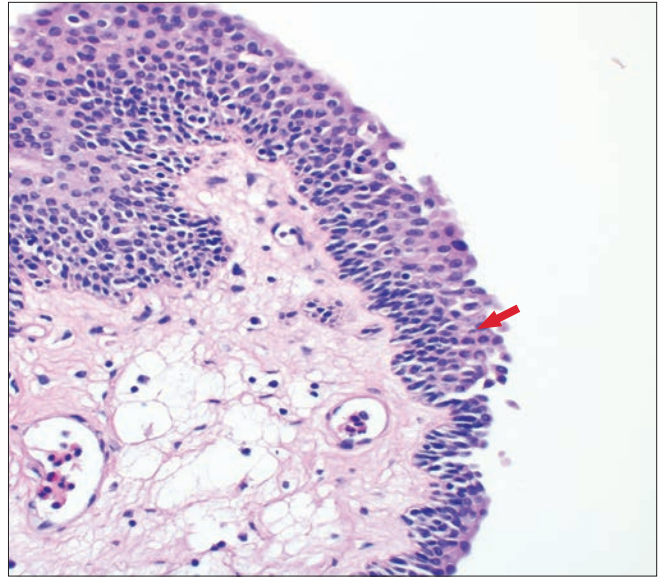


Figure 18a. 20x magnification.

Urothelial carcinoma of the bladder showing normal urothelium (red arrow) staining for PD-L1. When determining the PD-L1 % tumor cell expression for the specimen stained normal urothelium, dysplasia, and in situ carcinoma components are not included in the numerator, and the entire normal urothelium, dysplasia and in situ carcinoma components are not included in the denominator. Only the invasive component is evaluated when determining % tumor cell expression.

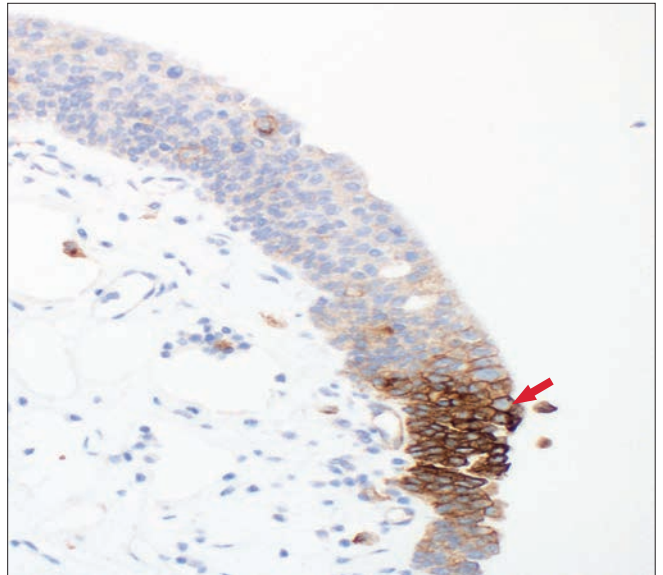


Figure 18b. 20x magnification.



# Challenging Cases for UC and MIUC

## PD-L1 IHC 28-8 pharmDx

### Nonspecific staining

Nonspecific staining is defined as staining that is not related to primary antibody-antigen interaction and represents issues such as reagent trapping, cartilage staining, DAB deposits, tissue folds, and edge drying. Other notable factors include, but are not limited to, pre-analytic fixation and processing of the specimen, incomplete removal of paraffin from sections, and incomplete rinsing of slides.

The use of fixatives other than 10% neutral buffered formalin may be a source of nonspecific staining. If nonspecific staining interferes with interpretation of specific staining, the slide may be considered unacceptable.

### Possible causes of nonspecific staining

- Improper drying of slides; ensure slides remain wet with buffer while loading onto Autostainer Link 48 and prior to initiating run
- Improper deparaffinization procedure
- Incomplete rinsing of reagents from slides
- Improper mixing of wash buffer

The nonspecific staining present on the patient specimen stained with Negative Control Reagent is useful in determining the level of nonspecific staining in the same patient tissue specimen stained with PD-L1. All specimens must have  $\leq 1+$  nonspecific staining.

### Immune cells

Intense staining of inflammatory cell infiltrate in the tumor may occur. Inflammatory cells are not included in determining the % tumor cell expression.

### Necrosis

Necrotic tissue may show nonspecific staining and should not be included in scoring % tumor cell expression.

*Note: If the specimen is excessively necrotic and contains < 100 viable tumor cells in the PD-L1 stained slide, the specimen is considered not evaluable.*

Urothelial carcinoma of the bladder.  
Tumor cells with granular staining  
(**black arrows**) in the cytoplasm and  
no linear membrane staining should not  
be included in scoring.

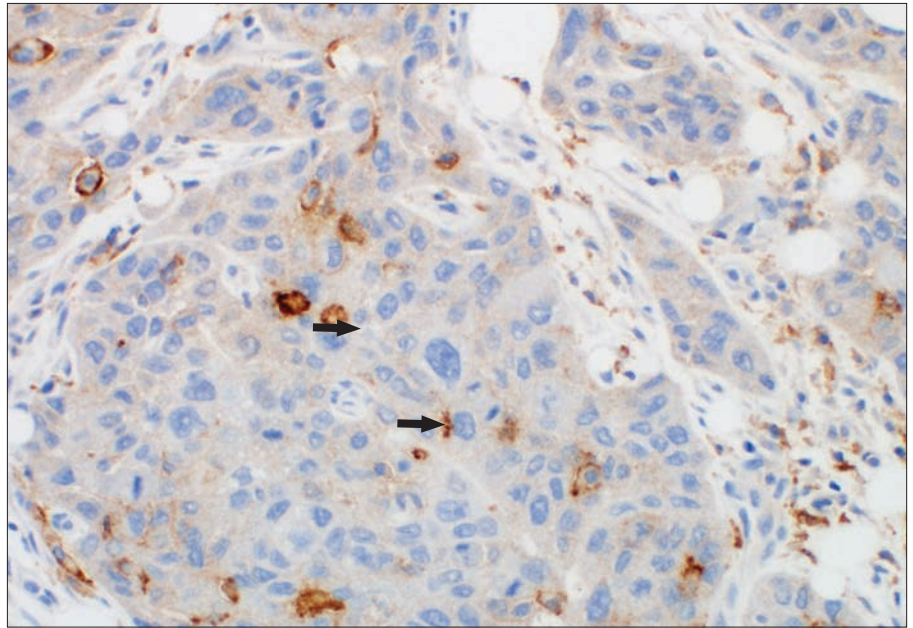


Figure 19. 20x magnification.

Urothelial carcinoma of the bladder.  
Cells with linear membrane staining that is  
distinguishable from cytoplasmic staining  
are included in scoring.

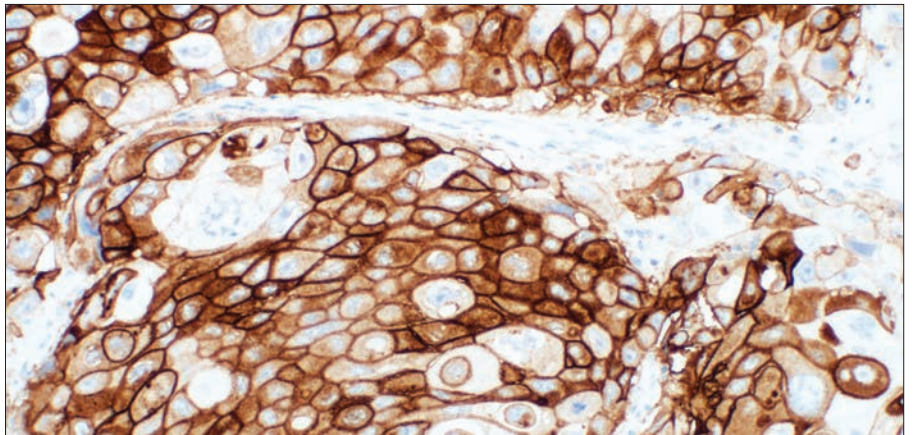


Figure 20. 20x magnification.

Urothelial carcinoma of the bladder.  
This example may be considered an indeterminate case if the excess cytoplasmic staining hampers scoring. Linear membrane staining of the tumor is observed (**black arrow**), however cytoplasmic staining is excessive in much of the specimen (**red arrow**).

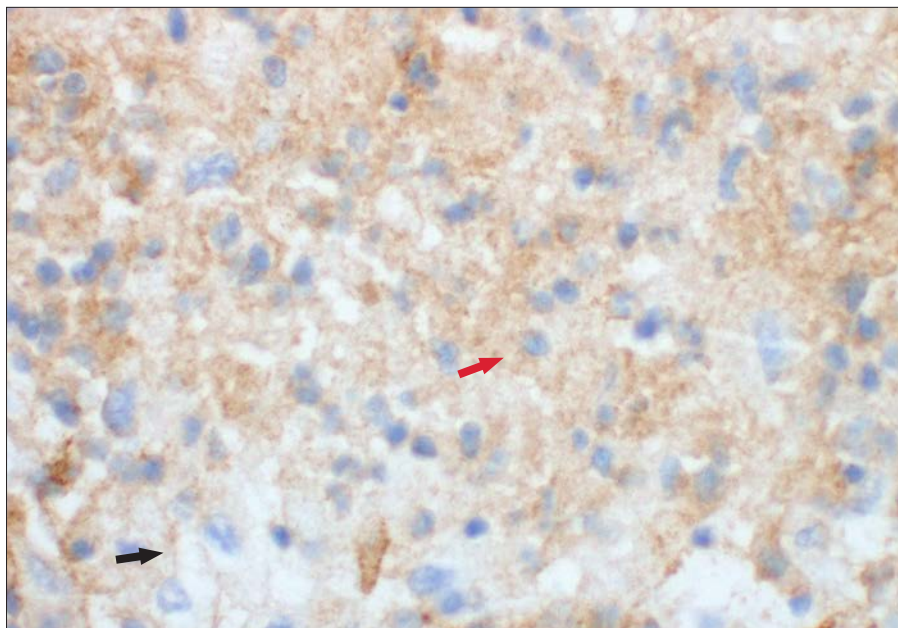


Figure 21. 40x magnification.

Urothelial carcinoma of the bladder.  
Note, plasma membrane staining of foreign body giant cells (FBGC) (**black arrows**) next to tumor cells. FBGCs are created from the fusion of macrophages reacting to foreign material extruded by tumor cells into the stroma. These FBGCs can be misidentified as tumor in the PD-L1 stained slide, but can be recognized as non-tumor in the H&E stained slide. Exclude FBGCs in the scoring of the tumor.

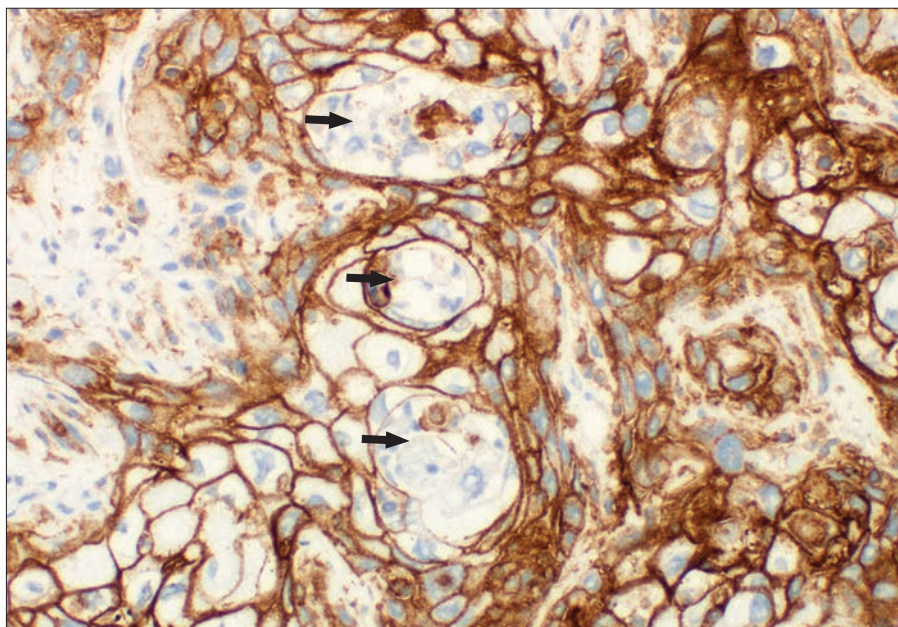


Figure 22. 20x magnification.



Urothelial carcinoma of the bladder. Necrotic tissue may show nonspecific staining and should not be included in scoring % tumor cell expression of the tumor. Care should be taken to only include viable tumor cells for scoring and not necrotic regions. If the specimen is excessively necrotic, the specimen is considered not evaluable. A minimum of 100 viable tumor cells should be present for evaluating the specimen. If the specimen is excessively necrotic and contains < 100 viable tumor cells, the specimen is considered non-evaluable (NE).

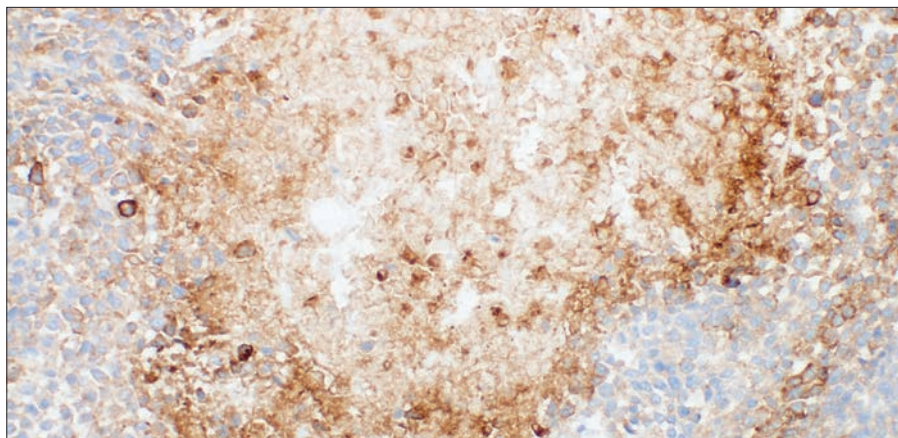


Figure 23. 20x magnification.

Urothelial carcinoma of the bladder with squamous differentiation (**black arrows**). Note, in order to confirm squamous differentiation, intracellular keratin, intercellular bridges, or keratin pearls (**red arrow**) are expected to be present and can be determined by using H&E stain.

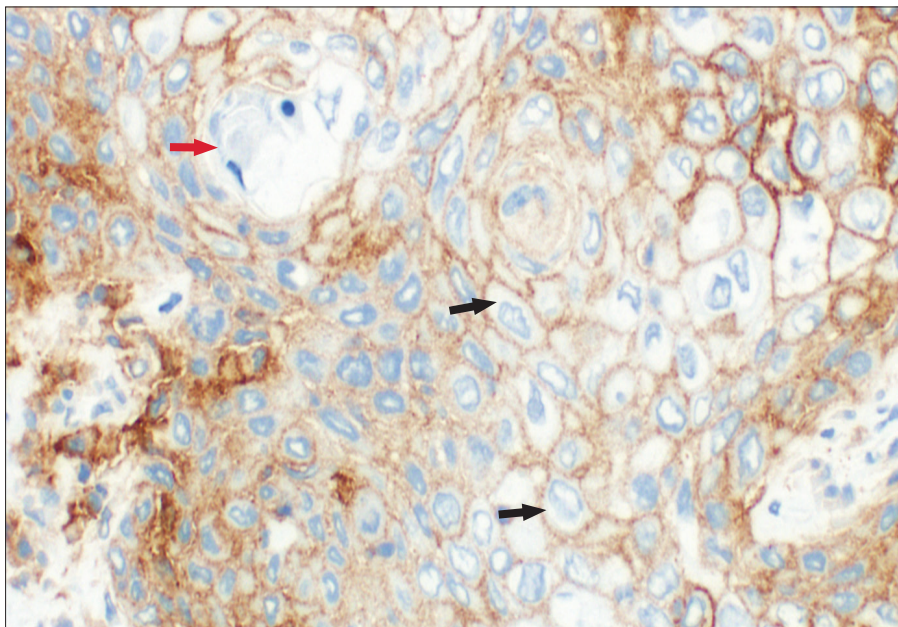
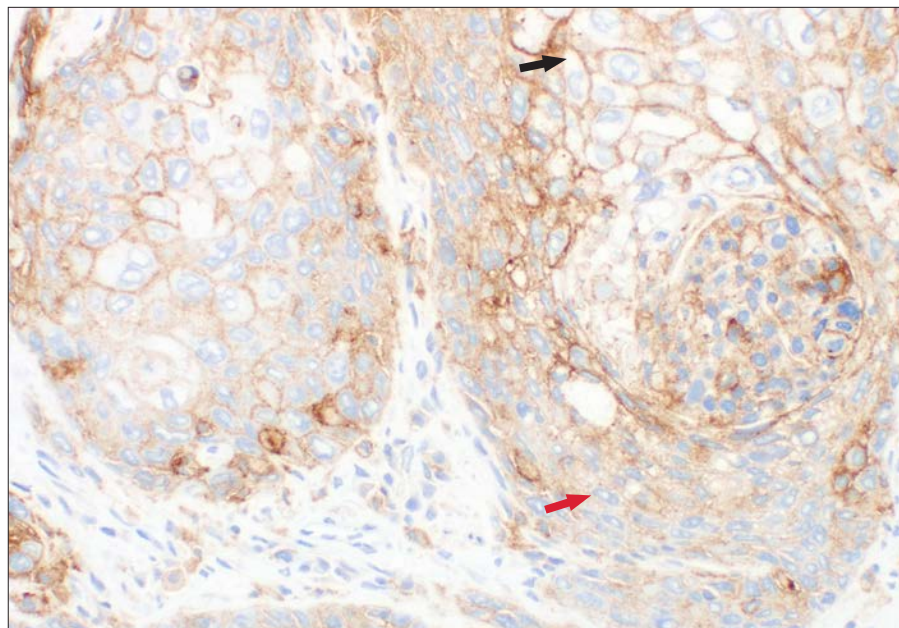


Figure 24. 20x magnification.

PD-L1 staining observed in urothelial carcinoma of the bladder. When scoring for % tumor cell expression, the percentage is determined by the number of stained cells and not area. Note in this example, the presence of staining in smaller tightly packed basaloid cells (**red arrow**) which take up less area than an equal number of well-differentiated staining cells (**black arrow**).



**Figure 25.** 20x magnification.



# Troubleshooting Guide for PD-L1 IHC 28-8 pharmDx

Problem	Probable Cause	Suggested Action
1. No staining of control or specimen slides	1a. Programming error	1a. Verify that the SK005 PD-L1 IHC 28-8 pharmDx program was selected for programming of slides
	1b. Lack of reaction with DAB+ Substrate-Chromogen Solution (DAB)	1b. Verify that DAB+ Substrate-Chromogen Solution was prepared properly
	1c. Sodium azide in wash buffer	1c. Use only EnVision FLEX Wash Buffer, Code K8007
	1d. Degradation of Control Slide	1d. Check kit expiration date and kit storage conditions on outside of package
2. Weak staining of specimen slides	2a. Inappropriate fixation method used	2a. Ensure that only 10% neutral buffered formalin fixative and approved fixation methods are used
	2b. Insufficient reagent volume applied	2b. Check size of tissue section and reagent volume applied
	2c. Inappropriate wash buffer used	2c. Use only EnVision FLEX Wash Buffer, Code K8007
3. Weak staining of specimen slides or the positive cell line on the Agilent-supplied Control Slide	3a. Inadequate target retrieval	3a. Verify that the 3-in-1 pretreatment procedure was correctly performed
	3b. Inappropriate wash buffer used	3b. Use only EnVision FLEX Wash Buffer, Code K8007
4. Excessive nonspecific staining of slides	4a. Paraffin incompletely removed	4a. Verify that the 3-in-1 pre-treatment procedure was correctly performed
	4b. Slides dried while loading onto the Autostainer Link 48	4b. Ensure slides remain wet with buffer while loading and prior to initiating run
	4c. Nonspecific binding of reagents to tissue section	4c. Check for proper fixation of the specimen and/or the presence of necrosis
	4d. Inappropriate fixation method used	4d. Ensure that only neutral buffered formalin fixative and approved fixation methods are used
	4e. Inadequate mixing of wash buffer	4e. Ensure wash buffer is properly mixed
	4f. Increased slide drying	<p>4f. Increase slide quantity per run to extend wash times for slides. Two full racks totaling 24 slides may demonstrate sufficient wash time increases to ameliorate slide drying.</p> <p>If further action is needed, three drop zones may be utilized to increase the amount of reagent on slides which may result in decreased slide drying.</p>
	4g. Improper humidity level maintained during run	<p>4g. Ensuring a complete closure of the instrument lid to create a sealed humidity chamber may decrease slide drying.</p> <p>Loading slide racks from the back to the front of the instrument away from airflow may decrease slide drying.</p> <p>If further action is needed, additional measures to increase humidity can be used.</p>

Problem	Probable Cause	Suggested Action
	4h. Instrument is not properly maintained	4h. Ensure that the instrument is maintained and calibrated per manufacturer's recommendations. Refer to Autostainer Plus Link and Autostainer Link 48 Basic User Guide for proper maintenance to ensure staining quality.
	4i. Autostainer slide racks are not level	4i. Ensure Autostainer slide racks are level prior to initiation of the IHC staining procedure. Please refer to Autostainer Plus Link and Autostainer Link 48 Basic User Guide for additional details.
	4j. Insufficient quality of DI water	4j. Ensure that distilled or de-ionized water (reagent-grade water) is being used and that the water is changed frequently. Refer to 'Materials Required, but Not Supplied' Section 5 for further details.
5. Tissue detached from slides	5a. Use of incorrect microscope slides	5a. Use FLEX IHC Microscope Slides (Code K8020), or Superfrost Plus slides
	5b. Inadequate preparation of specimens	5b. Cut sections should be placed in a $58 \pm 2$ °C oven for 1 hour prior to staining
6. Excessively strong specific staining	6a. Inappropriate fixation method used	6a. Ensure that only approved fixatives and fixation methods are used
	6b. Inappropriate wash buffer used	6b. Use only EnVision FLEX Wash Buffer, Code K8007
7. 1x EnVision FLEX Target Retrieval Solution is cloudy in appearance when heated	7. When heated the 1x EnVision FLEX Target Retrieval Solution turns cloudy in appearance	7. This is normal and does not influence staining
8. 1x EnVision FLEX Target Retrieval Solution does not meet pH specifications	8a. pH meter is not calibrated correctly	8a. Ensure pH meter is calibrated per manufacturer's recommendations. After re-calibration, re-test the pH of 1x EnVision FLEX Target Retrieval Solution. Do not modify the pH of 1x EnVision FLEX Target Retrieval Solution. If the pH is outside the acceptable range ( $6.1 \pm 0.2$ ), discard 1x EnVision FLEX Target Retrieval Solution. Prepare new 1x EnVision FLEX Target Retrieval Solution. Check the pH of the new 1x EnVision FLEX Target Retrieval Solution
	8b. Inferior quality water is used to dilute the EnVision FLEX Target Retrieval Solution concentrate	8b. Ensure that distilled or de-ionized water is used to prepare 1x Target Retrieval Solution
	8c. Incorrect Target Retrieval Solution is used	8c. Ensure that the correct EnVision Flex Target Retrieval Solution specified in "Materials Provided" and "Reagent Preparation" sections of the IFU is used

*Note: If the problem cannot be attributed to any of the above causes, or if the suggested corrective action fails to resolve the problem, please contact Agilent Pathology Support for further assistance. Additional information on staining techniques and specimen preparation can be found in the Education Guide: Immunohistochemical Staining Methods (Taylor C. R. and Rudbeck L. Education Guide: Immunohistochemical Staining Methods – Sixth Edition. Dako, Carpinteria, California. 2013; available from Agilent). Please refer to the latest version of the IFU for the most up to date Troubleshooting Guide: <https://www.agilent.com/en/library/eifu.html?searchTermRedirect=eifu>.*

## Bibliography

- Clinical and Laboratory Standards Institute (CLSI). Protection of Laboratory Workers From Occupationally Acquired Infections; Approved Guideline – Fourth Edition. CLSI document M29-A4 (ISBN 1-56238-962-9). CLSI, 950 West Valley Road, Suite 2500, Wayne, Pennsylvania 19087 USA; Vol.34 No.8; **2014**.
- Clinical and Laboratory Standards Institute (CLSI). Quality assurance for Design Control and Implementation of Immunohistochemistry Assays; Approved guideline – Second Edition. CLSI document I/LA28-A2 (ISBN1-56238-745-6). CLSI, 950 West Valley Road, Suite 2500, Wayne, Pennsylvania 19087 USA; Vol.31 No.4; **2011**.
- Department of Health, Education and Welfare, National Institutes for Occupational Safety and Health, Rockville, MD. Procedures for the decontamination of plumbing systems containing copper and/or leadazides. DHHS (NIOSH) Publ. No. 78-127. **1976**.
- Omata M. D., C-T Liew., et al. Nonimmunologic Binding of Horseradish Peroxidase to Hepatitis B Surface Antigen: A Possible Source of Error in Immunohistochemistry. *Am. J. Clin. Pathol.* **1980**, 73(5), 626-632.
- PD-L1 IHC 28-8 pharmDx Instructions for Use
- Phelps R. M., Johnson B. E., et al. NCI-navy medical oncology branch cell line database. *J. Cell. Biochem.* **1996**, 63(S24), 32-91.
- Therese Phillips M. A., Pauline Simmons B. S., et al. Development of an automated PD-L1 immunohistochemistry (IHC) assay for Non-Small Cell Lung Cancer. *Appl. Immunohistochem. Mol. Morphol.* **2015**, 23(8), 541-549.
- Taylor C. R. and Rudbeck L. Education Guide: Immunohistochemical Staining Methods – Sixth Edition. Dako, Carpinteria, California; **2013**.
- Topalian S. L., Drake C. G., Pardoll D. M. Targeting the PD-1/B7-H1 (PD-L1) pathway to activate anti-tumor immunity. *Curr. Opin. Immunol.* **2012**, 24(2), 207-212.
- Topalian S. L., Hodi F. S., et al. Safety, Activity, and Immune Correlates of Anti-PD-1 Antibody in Cancer. *N. Engl. J. Med.* **2012**, 366, 2455-2465.
- Wang C., Kent B. T., et al. *In Vitro* Characterization of the Anti-PD-1 Antibody Nivolumab, BMS-936558, and *In Vivo* Toxicology in Non-Human Primates. *Cancer Immunol. Res.* **2014**, 2(9), 846-856.

Learn more:

[www.agilent.com/en/product/pharmdx/pd-l1-ihc-28-8-pharmdx](http://www.agilent.com/en/product/pharmdx/pd-l1-ihc-28-8-pharmdx)

Buy online:

[www.agilent.com/chem/store](http://www.agilent.com/chem/store)

Get answers to your technical questions and  
access resources in the Agilent Community:

[community.agilent.com](http://community.agilent.com)

U.S. and Canada

**1-800-227-9770**

**[agilent\\_inquiries@agilent.com](mailto:agilent_inquiries@agilent.com)**

Europe

**[info\\_agilent@agilent.com](mailto:info_agilent@agilent.com)**

Asia Pacific

**[inquiry\\_lsca@agilent.com](mailto:inquiry_lsca@agilent.com)**

D74745\_2.00

This information is subject to change without notice.

© Agilent Technologies, Inc. 2023, 2025  
Published in the USA, September 22, 2025  
29571 2025JUN26

