

Naveni® PD1/PD-L1 (HRP) on the BOND RX/RX^m staining instruments

The following pair of antibodies have been optimized on the BOND RX and RX^m on various normal and cancer tissue types.

1. PD-1 antibody based on clone EH33 CST
2. PD-L1 antibody based on clone SP142 Abcam RabMAb®

The kit instruction for the target specific Naveni® PD1/PD-L1 BOND HRP interaction assay will replace **section 2.1 and 2.2** associated with “**Reagent preparation**” in the NaveniBright™ BOND kit instruction (NAB.MR.030.H). Please follow the rest of the NaveniBright Kit instructions to prepare all other reagents and run the assay. Use the default antigen retrieval method ER2 for 40 min at 100°C for optimal detection of PD1/PD-L1 in various tissues.

For research use only.

Kit box 3, Bag:

Storage at -25 to -15°C. FREEZE



Material	Art.no	Amount
Naveni BOND PD-1 antibody 40x	NA.2.030.30	150 µL
Naveni BOND PD-L1 antibody 40x	NA.2.030.31	150 µL

Solution preparation

Primary antibodies

1. Prepare the **Diluent 1** by diluting **Supplement 2, 1:9** in **Diluent 1 (1x)**, see **Table 2** for an example calculation. The calculation is based on 10 slides with a volume of 150 µL per slide. A dead volume of 500 µL is included in the total volume.
2. Use the prepared **Diluent 1** to dilute both the **Naveni PD-1** and **PD-L1** antibodies 1:40 in the same container (**Table 3**). The calculation is based on 10 slides with a volume of 150 µL per slide. A dead volume of 500 µL is included in the total volume.

Table 2. Example calculation: Diluent 1

Kit component	Diluent 1 (µL)
Diluent 1	1778
Supplement 2	222
Total	2000

Table 3. Example calculation: Primary antibody dilution

Kit component	Diluent 1 (µL)
Diluent 1 with Supplement 2	1900
Naveni BOND PD-1 antibody (40x)	50
Naveni BOND PD-L1 antibody (40x)	50
Total	2000