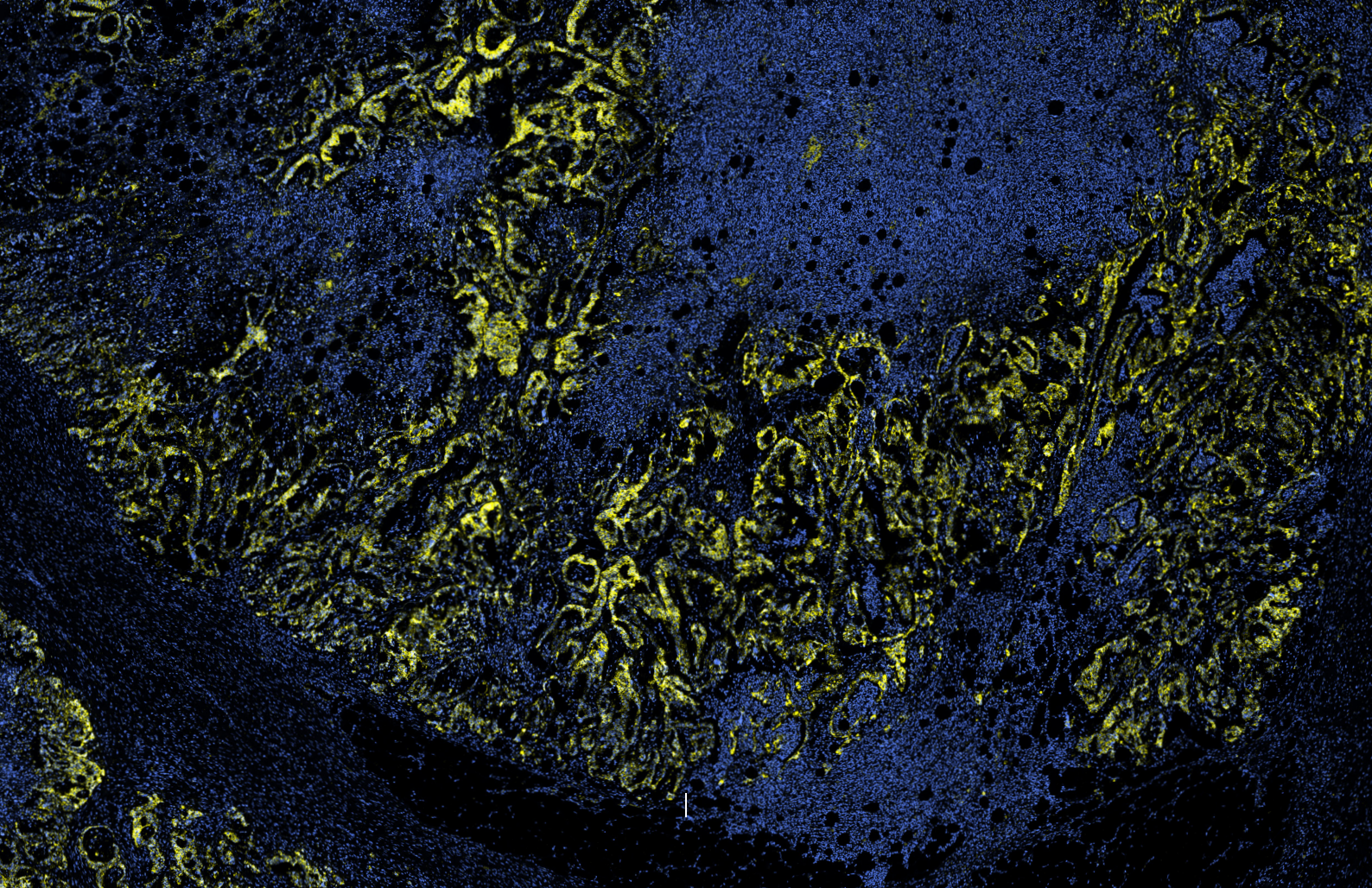




Staining of FFPE Colon Cancer tissue, E-Cadherin/β-Catenin (yellow) and nuclear stain with DAPI (blue).

NaveniFlex™ BOND RX Atto647N/Red

ILLUMINATING FUNCTION IN SPATIAL PROTEOMICS

Item No. 42973

Automated detection of protein-protein interactions and post-translational modifications *in situ*

The NaveniFlex BOND RX Atto647N/Red integrates fluorescence readout automation on the BOND RX/RX™ Fully Automated Research Stainer. These *in situ* kits are meticulously crafted based on our cutting-edge Naveni® *in situ* Proximity Ligation Technology, offering flexibility tailored to your unique primary antibodies and targets.

NaveniFlex BOND RX Atto647N/Red enables you to:

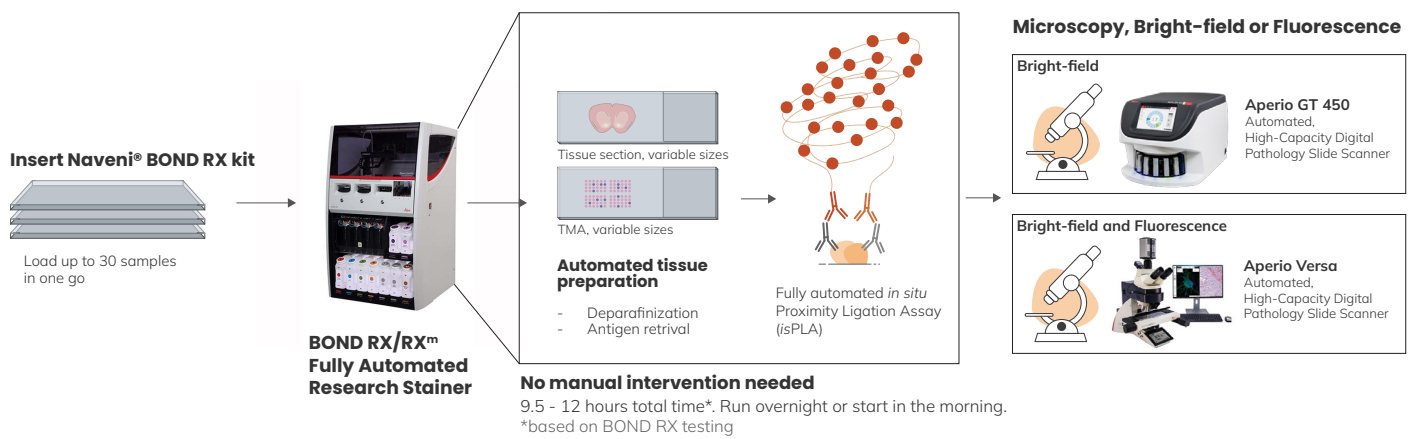
- Detect protein-protein interactions, post-translational modifications, and/or specific protein targets efficiently using the Fully Automated Research Stainer BOND RX/RX™
- Achieve high throughput and enhance your research efficiency and reproducibility through a seamlessly integrated automated workflow on the BOND RX/RX™.
- Save valuable time with the automated workflow, reducing hands-on time, and maximizing overall research productivity.



In partnership with Leica Biosystems

For additional information and images, read more at navinci.se/technology/naveni-bond



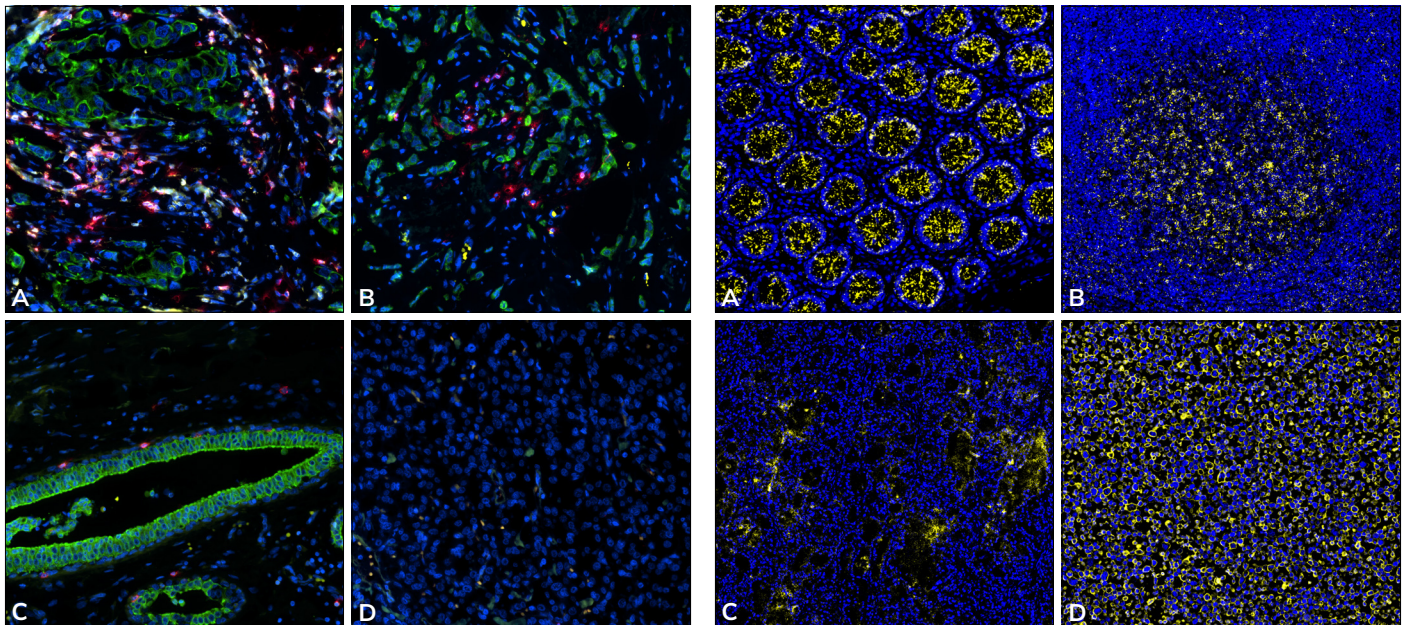


The NaveniFlex BOND RX Atto647N/Red kits include two Navenibodies conjugated to proprietary oligo arms (depicted as orange antibodies in the illustration), directed to the primary antibodies with host origin of mouse and rabbit respectively. Only if the Navenibodies are in close proximity will they generate a rolling circle amplification reaction, leading to a strong and distinct signal.

The kit has undergone thorough validation, including a diverse range of FFPE tissues and five different interaction assays

(PD1/PD-L1, CD8/MHC-I, Mesothelin/Mucin, E-Cadherin/Beta-Catenin and Zap70/Lat) and a dual recognition assay (HER2/HER2). The validation protocol entailed an extensive comparative analysis at three distinct research sites, ensuring the establishment of robust performance characteristics across varied experimental settings.

Please note that the corresponding kit is also available with chromogenic read-out: NaveniBright BOND RX HRP.



Staining of CD8/MHC-I interaction (pink), CD3 (yellow), pan-cytokeratin (green) nuclear stain (DAPI-blue) in a TMA, A) Breast cancer, B) breast cancer, C) healthy breast tissue, D) Adrenal gland (biological negative control) using NaveniFlex BOND RX Atto647N.

A) Staining of B-catenin/E-cadherin interaction in healthy colon tissue, B) Zap 70/LAT interaction in tonsil tissue, C) Mesothelin/Mucin interaction in ovary cancer tissue, D) staining of HER2/HER2 single protein in BT474 cell line.

Available from Navinci

Catalog nr	Kit	Target	Read out	Description
60009/60010	NaveniFlex BOND RX Atto647N/Red	Your choice, use primary antibodies with host origin of mouse and rabbit	Fluorescence	Anti-mouse Navenibody Anti-rabbit Navenibody Buffers for blocking and dilutions and detection reagents Reagents sufficient for 30 FFPE tissue slides, including dead volumes*
60011/60012	Naveni PD1/PD-L1 BOND RX Atto647N/Red	Human PD1/PD-L1 interaction	Fluorescence	Navenibody targeting human PD1 protein based on clone EH33 CST Navenibody targeting human PD-L1 protein based on clone SP142 Abcam RabMAb® Buffers for blocking and dilutions and detection reagents for the PD1/ PD-L1 interaction signal Reagents sufficient for 30 FFPE tissue slides, including dead volumes*

*additional reagents required, read more at navinci.se/technology/naveni-bond
Research use only, not for use in diagnostic procedures