

DEVELOPMENT OF A GLYCOSYLATION-INDEPENDENT, IHC-VALIDATED RECOMBINANT RABBIT MONOCLONAL ANTIBODY AGAINST THE CANCER STEM CELL MARKER CD133

Antony W. Wood, Richard Cotta, Matthew Prokop, and Jessica Simendinger
Cell Signaling Technology, Inc., Danvers MA 01923

T-2147

ABSTRACT

The cancer stem cell hypothesis postulates that a population of self-renewing tumor-initiating cells, termed Cancer Stem Cells (CSCs), may be responsible for driving tumor heterogeneity, metastasis, therapeutic resistance and/or tumor relapse. Tools to identify and characterize putative CSCs are therefore of significant value for the cancer research community. CD133 is a 5-transmembrane (5-TM) cell surface glycoprotein that shows elevated expression in putative CSCs from multiple tumor types. Numerous studies have used antibodies directed against CD133 to isolate putative CSCs for characterization, *in vitro* culture, transplantation and drug discovery studies. However, the most commonly used antibodies used to study CD133⁺ CSCs are raised against glycosylated CD133 epitopes; this is problematic because the glycosylation status of CD133 varies in response to environmental conditions (e.g., hypoxia) or cell differentiation status. Furthermore, available anti-CD133 antibodies have not been rigorously validated for immunohistochemistry, which is critical for understanding CSC biology *in situ*. To address this problem, we have developed a recombinant rabbit monoclonal antibody that targets an extracellular, glycosylation-independent epitope of CD133. This reagent [CD133 (D4W4N) XP[®] Rabbit mAb] has been rigorously validated in Western immunoblot and immunohistochemistry, where it demonstrates robust and specific staining of CD133 protein across diverse cell and tissue types. It is hoped that this reagent will enable accelerated progress in elucidating the role of putative cancer stem cells in the development, metastasis, therapeutic resistance and relapse of tumors.

METHODS

Monoclonal Antibody Generation

CD133 (D4W4N) XP[®] Rabbit mAb #86781 is a recombinant rabbit monoclonal antibody, generated at Cell Signaling Technology, Inc. using patented XMT[®] Technology.

Western Blot

Western blot analyses were performed using extracts from cell or tissue extracts, as described in the figure legends. Western blot protocol details can be found at <http://www.cellsignal.com/wbprotocol>

Immunohistochemistry (IHC)

Paraffin-embedded tissue sections were deparaffinized and rehydrated, then subjected to antigen retrieval in sodium citrate, pH 6.0. Primary antibodies were incubated overnight at 4°C. Detection was performed using SignalStain[®] Boost IHC Detection Reagent (HRP, Rabbit) #8114 and SignalStain[®] DAB Substrate Kit #8059.

SUMMARY

CD133 (D4W4N) XP[®] Rabbit mAb:

A recombinant rabbit monoclonal antibody that detects human CD133 protein in western blot and IHC with exceptional specificity and sensitivity.

Detection of CD133 is independent of protein glycosylation status (determined by Western blot).

May be used to detect or visualize CD133 in human cell lysates or formalin-fixed paraffin-embedded (FFPE) human normal and tumor tissues.

Exhibits greater sensitivity in IHC when compared to available anti-CD133 antibodies.

Antibody Validation – Cell Model Selection

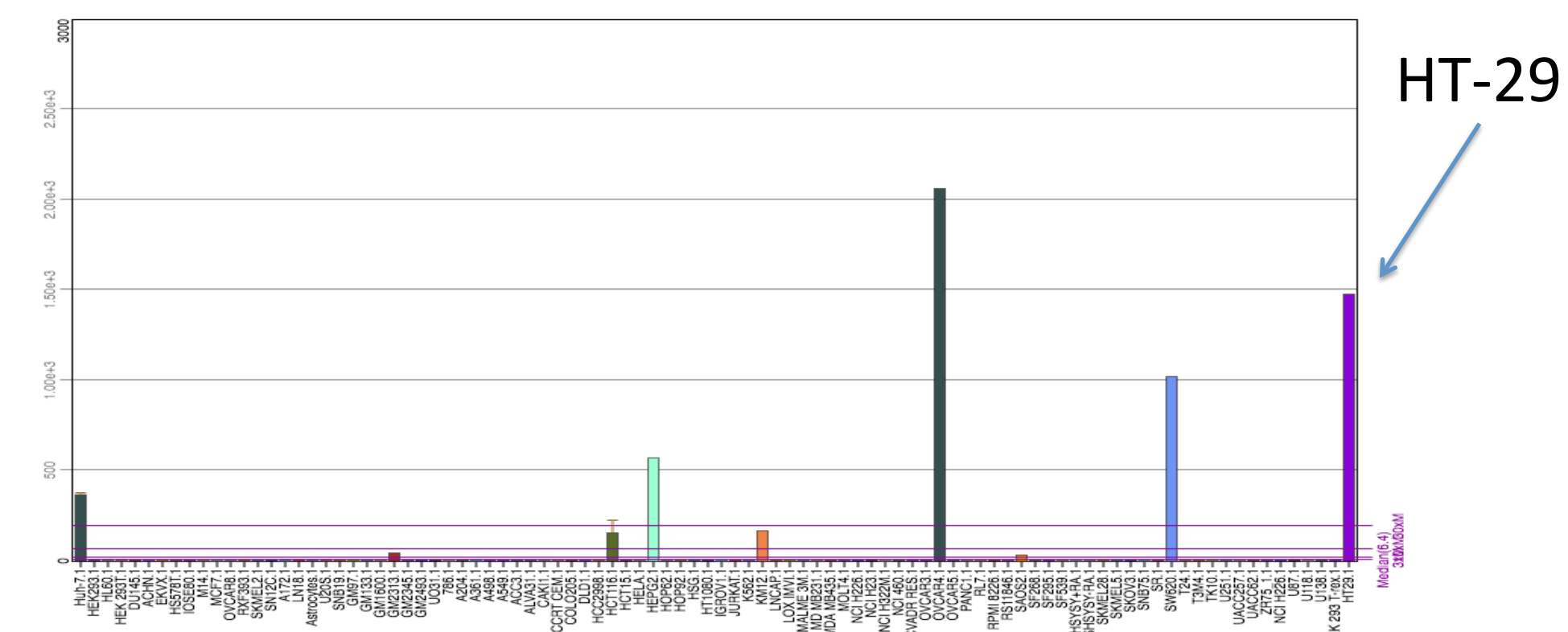


Fig. 1. Messenger RNA expression pattern of human CD133, as reported in BioGPS (www.biogps.org).

Antibody Validation – Western Blotting

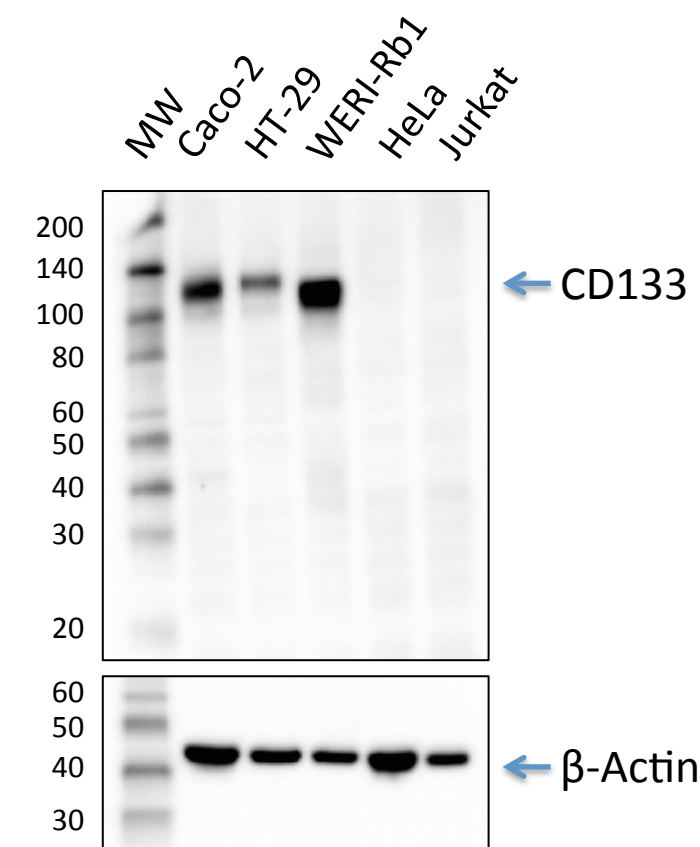


Fig. 2. Western blot analysis of selected human cancer cell lines using CD133 (D4W4N) XP[®] Rabbit mAb #86781 (upper) and β -actin (D6A8) Rabbit mAb #8457 (lower). CD133 protein is robustly detected in confirmed CD133⁺ cancer cell lines, whereas no CD133 protein is detected in CD133⁻ cell lines.

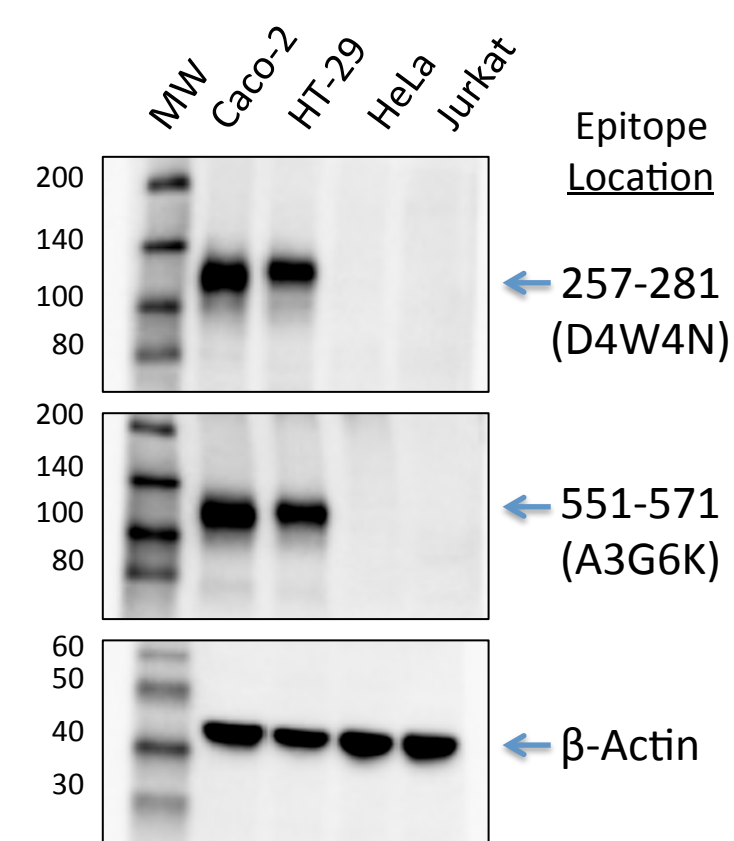


Fig. 3. Western blot analysis of CD133⁺ and CD133⁻ human cancer cell lines, using CD133 (D4W4N) XP[®] Rabbit mAb #86781 (upper), CD133 (A3G6K) Rabbit mAb #5860 (middle) and β -actin (D6A8) Rabbit mAb #8457 (lower). Antibody epitope locations are indicated to the right of respective Western blot images. Detection of CD133 is identical between clones, despite targeting distinct epitopes.

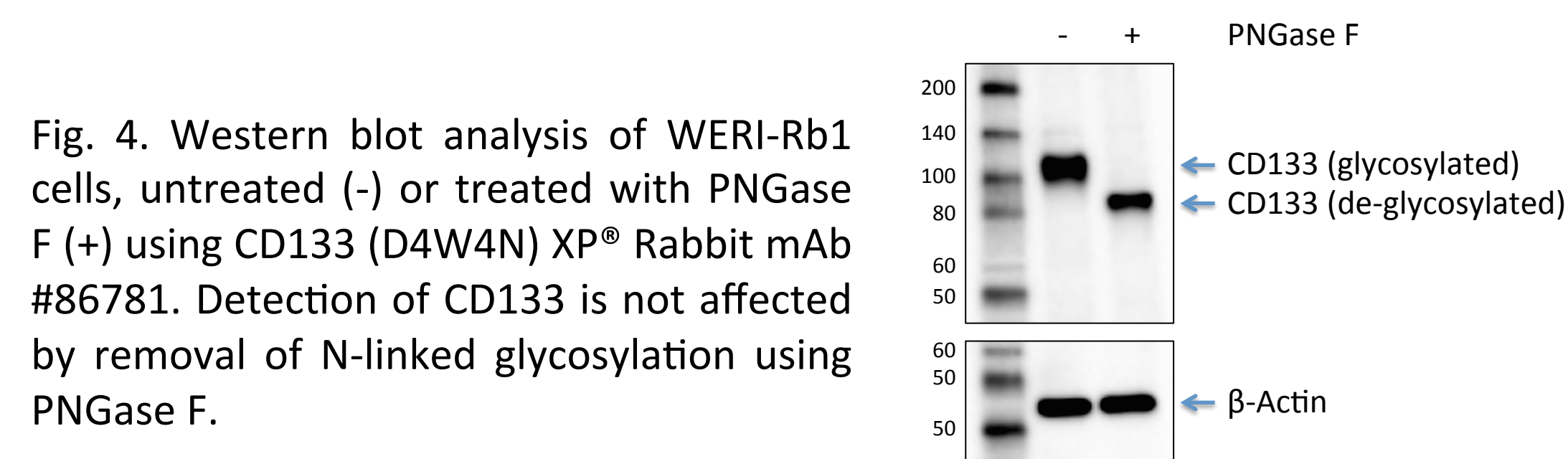


Fig. 4. Western blot analysis of WERI-Rb1 cells, untreated (-) or treated with PNGase F (+) using CD133 (D4W4N) XP[®] Rabbit mAb #86781. Detection of CD133 is not affected by removal of N-linked glycosylation using PNGase F.

Immunohistochemistry – Cell Pellet Validation

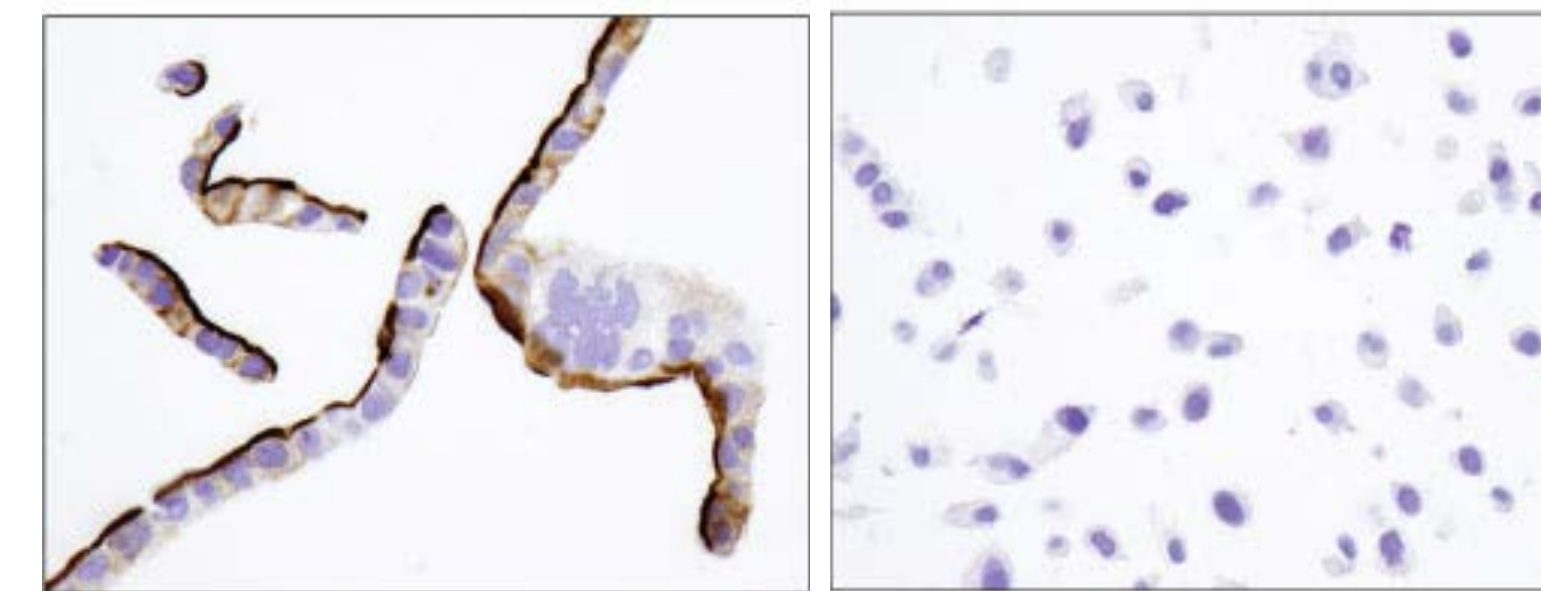


Fig. 5. Immunohistochemical analysis of paraffin-embedded Caco-2 (left) and HeLa (right) cell pellets, using CD133 (D4W4N) XP[®] Rabbit mAb #86781. A strong signal is detected in the CD133⁺ cancer cell line (Caco-2), but signal is completely absent in the CD133⁻ cell line (HeLa).

Immunohistochemistry – FFPE human normal tissues

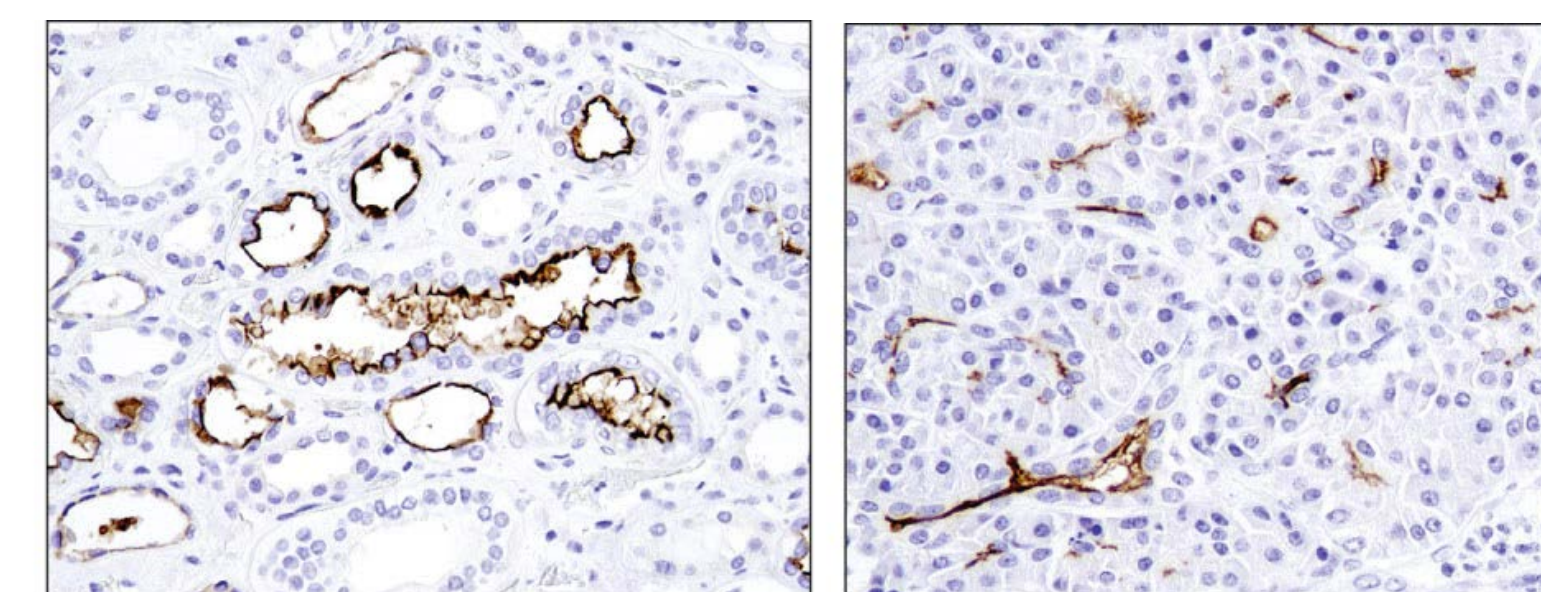


Fig. 6. Immunohistochemical analysis of paraffin-embedded human normal kidney (left) and human normal pancreas (right), using CD133 (D4W4N) XP[®] Rabbit mAb #86781. CD133 protein is detected in kidney tubules and pancreatic acinar cells.

Immunohistochemistry – FFPE human tumor tissues

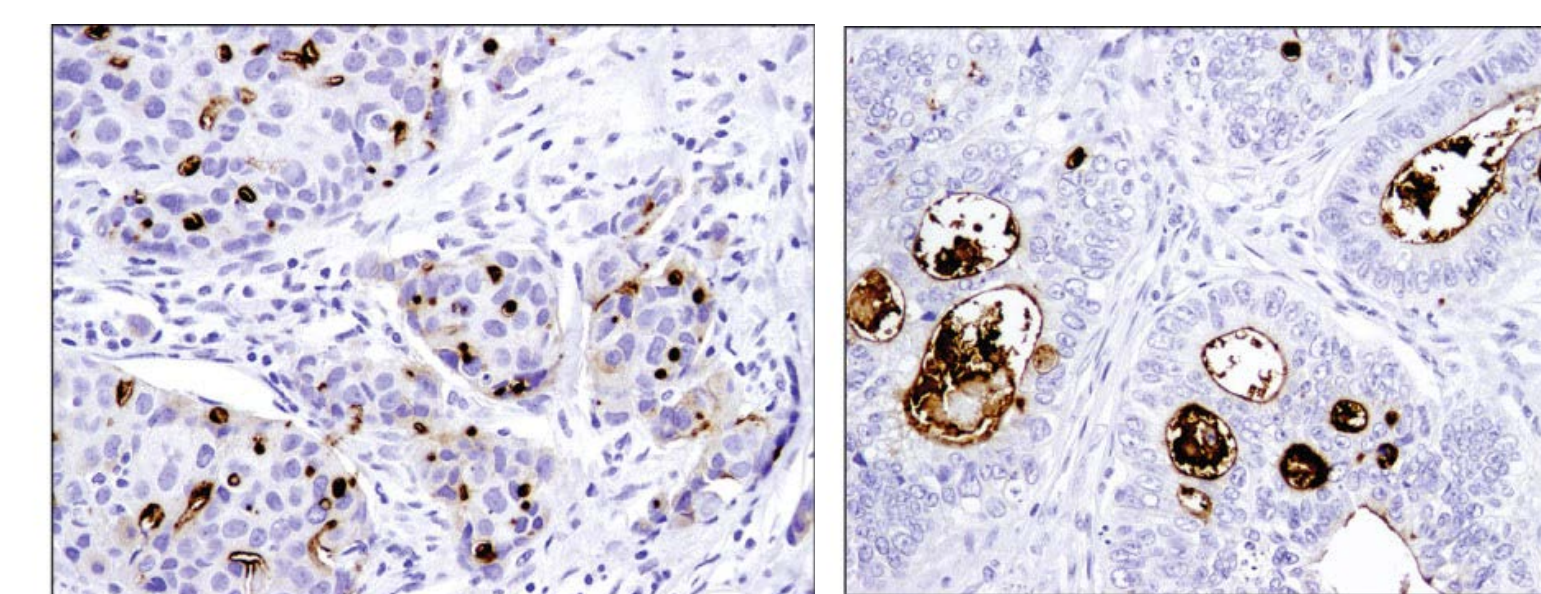


Fig. 7. Immunohistochemical analysis of paraffin-embedded human ductal carcinoma of the breast (left) and human colon carcinoma (right), using CD133 (D4W4N) XP[®] Rabbit mAb #86781.

Immunohistochemistry – Performance Comparison

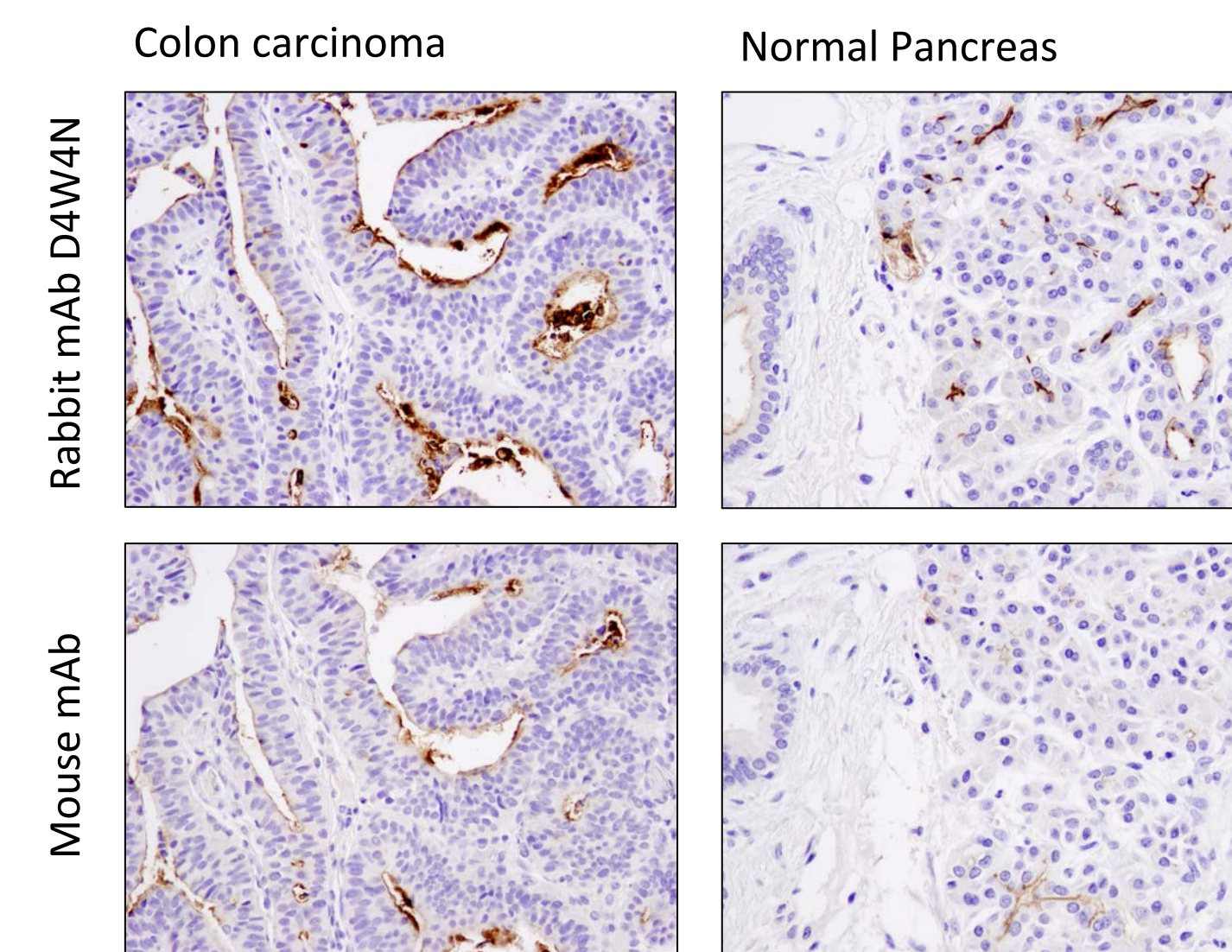


Fig. 8. Immunohistochemical analysis of paraffin-embedded human colon carcinoma (left panels) and paraffin-embedded normal human pancreas (right panels), using CD133 (D4W4N) XP[®] Rabbit mAb (upper panels), and an anti-CD133 mouse monoclonal antibody (lower panels). CD133 (D4W4N) XP[®] Rabbit mAb exhibits more robust detection of CD133 protein in both normal and tumor tissues when compared to the mouse monoclonal antibody.