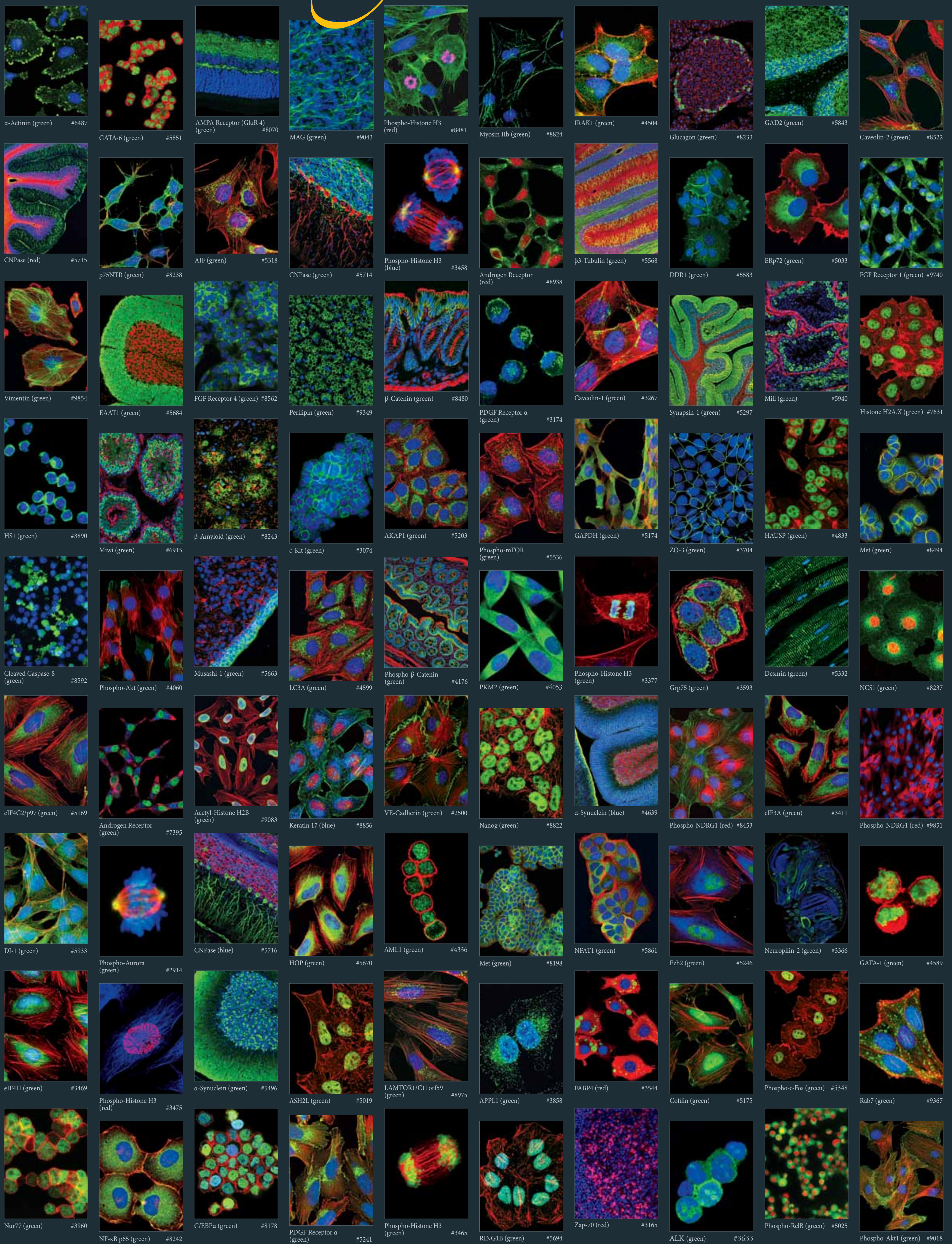


immunofluorescence

featuring **XP[®]** Monoclonal Antibodies



All images were captured at Cell Signaling Technology by our research application scientists. For more information and comprehensive validation data on over 4,000 antibodies visit: www.cellsignal.com

XP[®] Monoclonal Antibodies
one antibody, multiple applications™

XP[®] monoclonal antibodies are a line of high quality rabbit monoclonal antibodies exclusively available from Cell Signaling Technology. Any product labeled with XP has been carefully selected based on superior performance in the most relevant research applications.

exceptional specificity
+ exceptional sensitivity
+ exceptional stability and reproducibility
= exceptional performance™

Our proprietary XMT[™] technology coupled with extensive antibody validation and stringent quality control delivers XP monoclonal antibodies with exceptional performance in the widest range of applications.



Cell Signaling
 TECHNOLOGY[®]