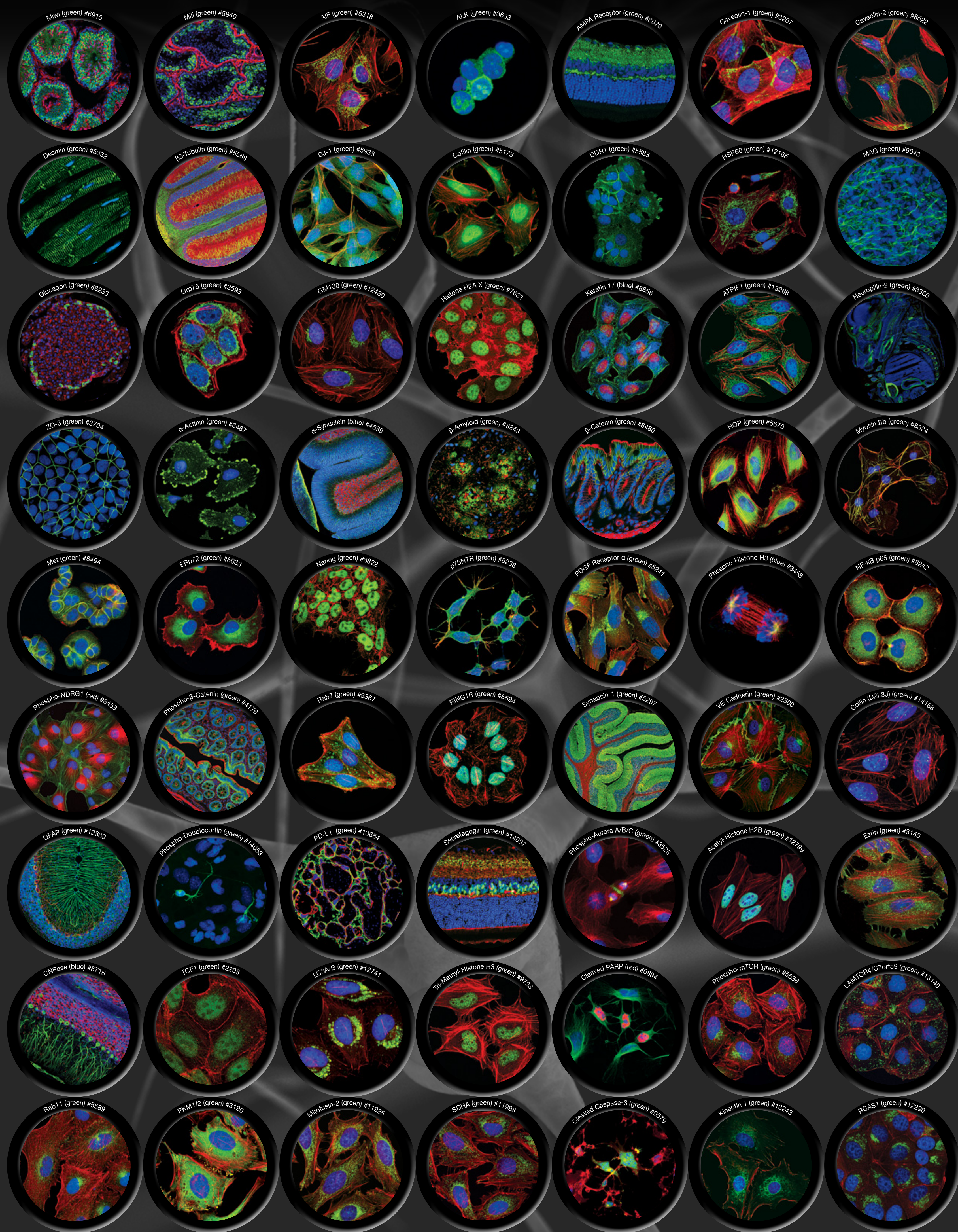


Immunofluorescence



All images were captured at Cell Signaling Technology by our research application scientists.

For more information and comprehensive validation data on our antibodies visit: www.cellsignal.com/ifs

For Research Use Only. Not For Use in Diagnostic Procedures.

© 2014 Cell Signaling Technology, Inc. Cell Signaling Technology® is a registered trademark of Cell Signaling Technology, Inc.



Cell Signaling
TECHNOLOGY®