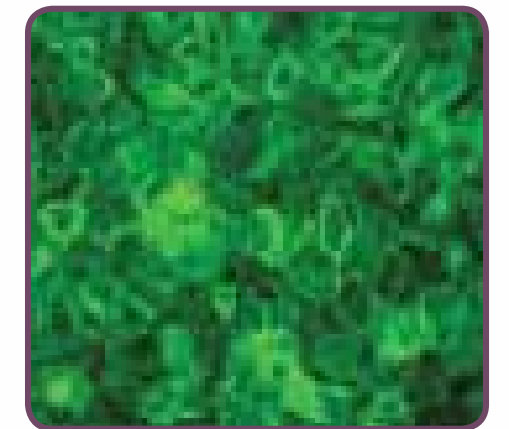
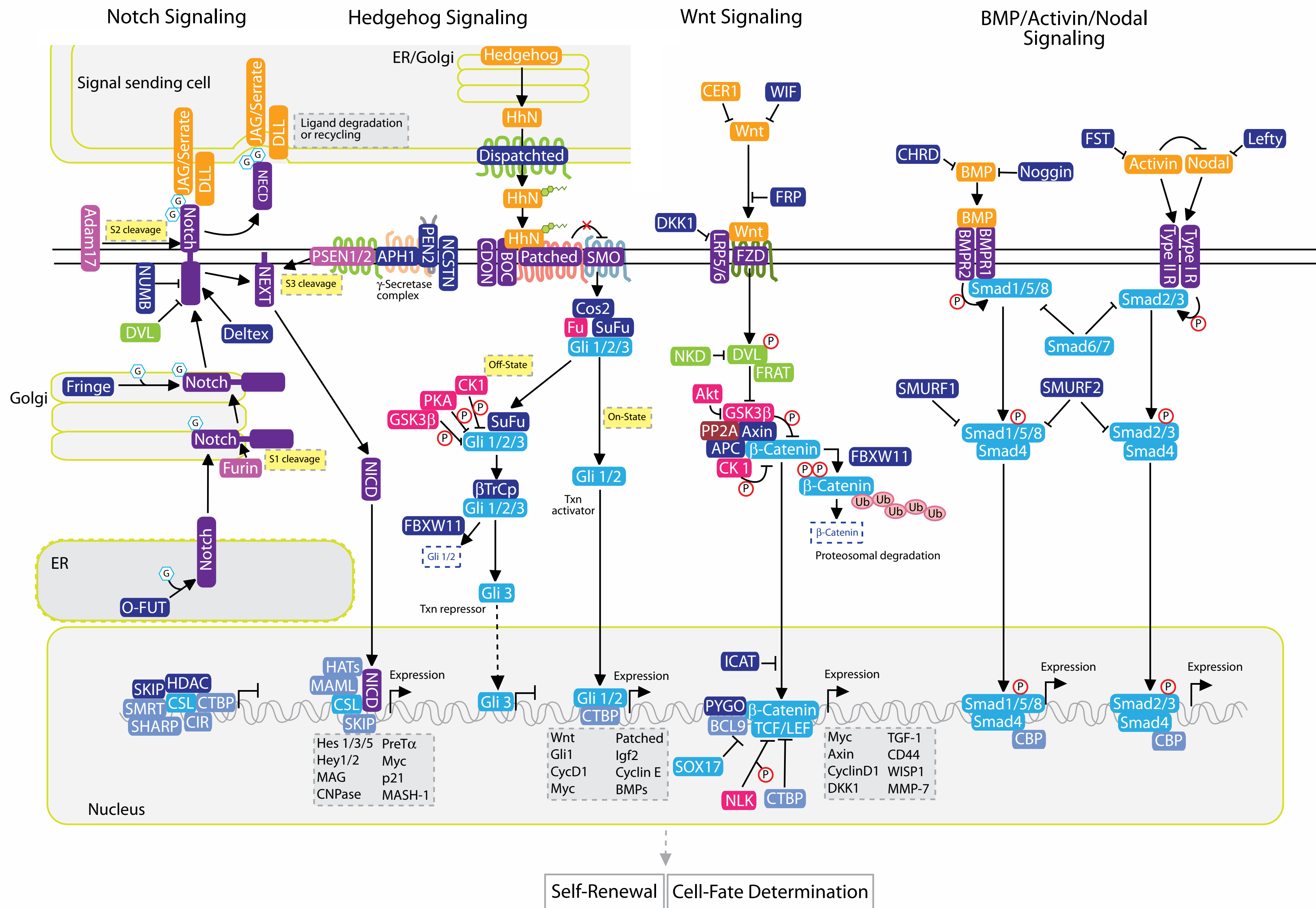
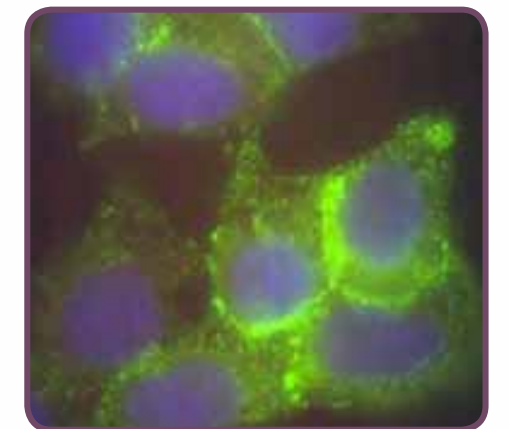


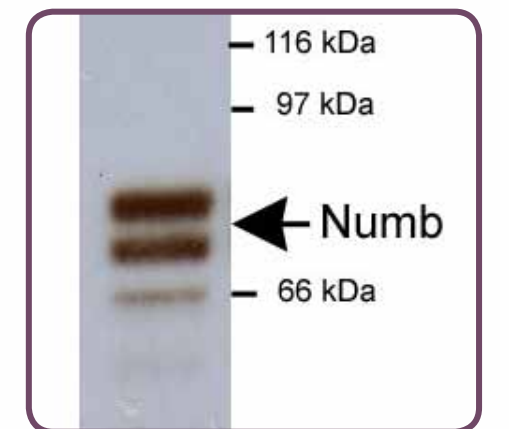
Stem Cell Signaling Pathway



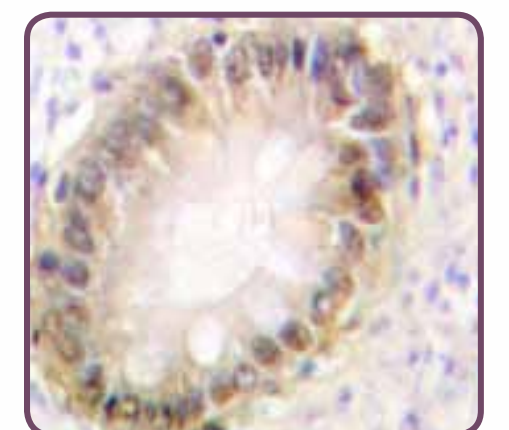
NB110-56924 • Beta Catenin



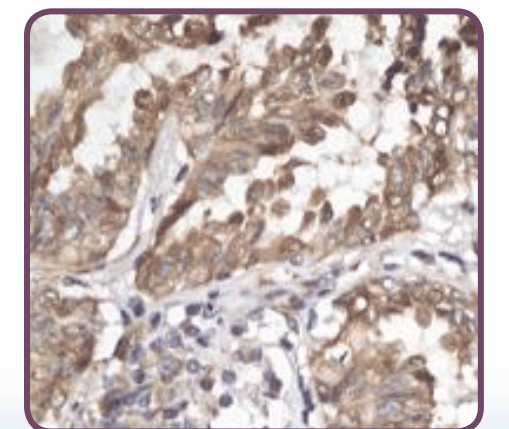
NB600-1161 • Jagged 1



NB500-178 • NUMB



NBP1-19520 • Smad2 / 3



NB110-56923 • Sonic Hedgehog