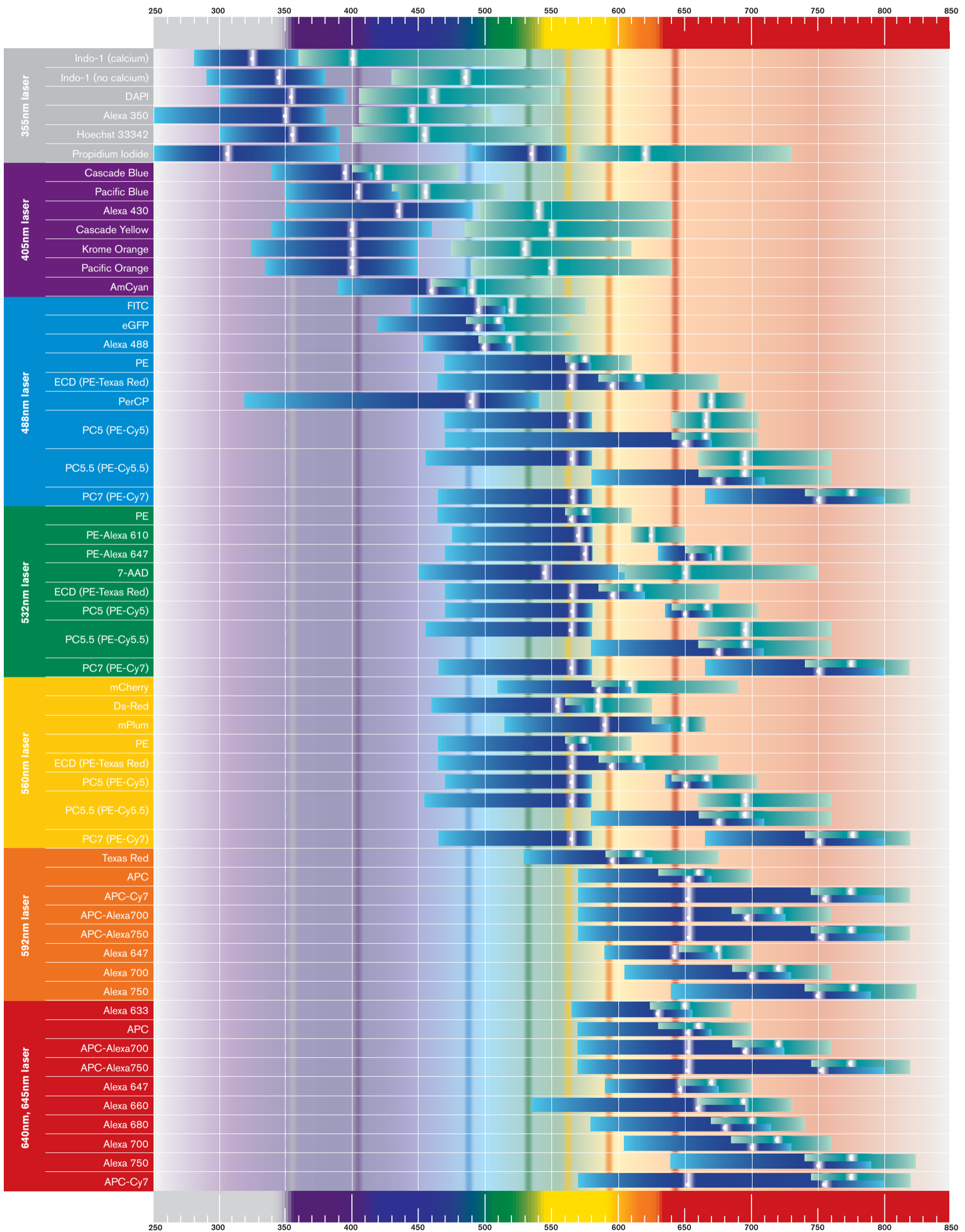
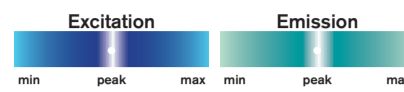


Fluorochrome Guide

MoFlo XDP and MoFlo Astrios



Standard MoFlo XDP Fluorochromes

355nm LASER: Indo-1 (Ca), Indo-1, DAPI, Hoechst, PI **405nm LASER:** Cascade Blue, Pacific Blue, Cascade Yellow **488nm LASER:** FITC, PE, ECD, PE-Cy5, PE-Cy7 **532nm LASER:** PE, PE-Alexa 610, PE-Alexa 647, ECD, PE-Cy5, PE-Cy7 **560nm LASER:** PE, ECD, PE-Cy5, PE-Cy7 **592nm LASER:** APC, APC-Cy7, Alexa 647, Alexa 700, Alexa 750 **640nm LASER:** APC, Alexa 647, Alexa 700, Alexa 750, APC-Cy7

Standard MoFlo Astrios Fluorochromes

355nm LASER: Hoechst (bound), Hoechst (unbound), PI **405nm LASER:** Pacific Blue, Cascade Yellow **488nm LASER:** FITC, PE, ECD, PE-Cy5, PE-Cy5.5, PE-Cy7 **532nm LASER:** PE, PE-Alexa 610, PE-Alexa 647, ECD, PE-Cy5, PE-Cy5.5, PE-Cy7 **560nm LASER:** mCherry, DsRed, mPlum, ECD, PE-Cy5, PE-Cy5.5, PE-Cy7 **592nm LASER:** TR, APC, APC-Cy7, Alexa 700 **645nm LASER:** APC, APC-Cy7, Alexa 700