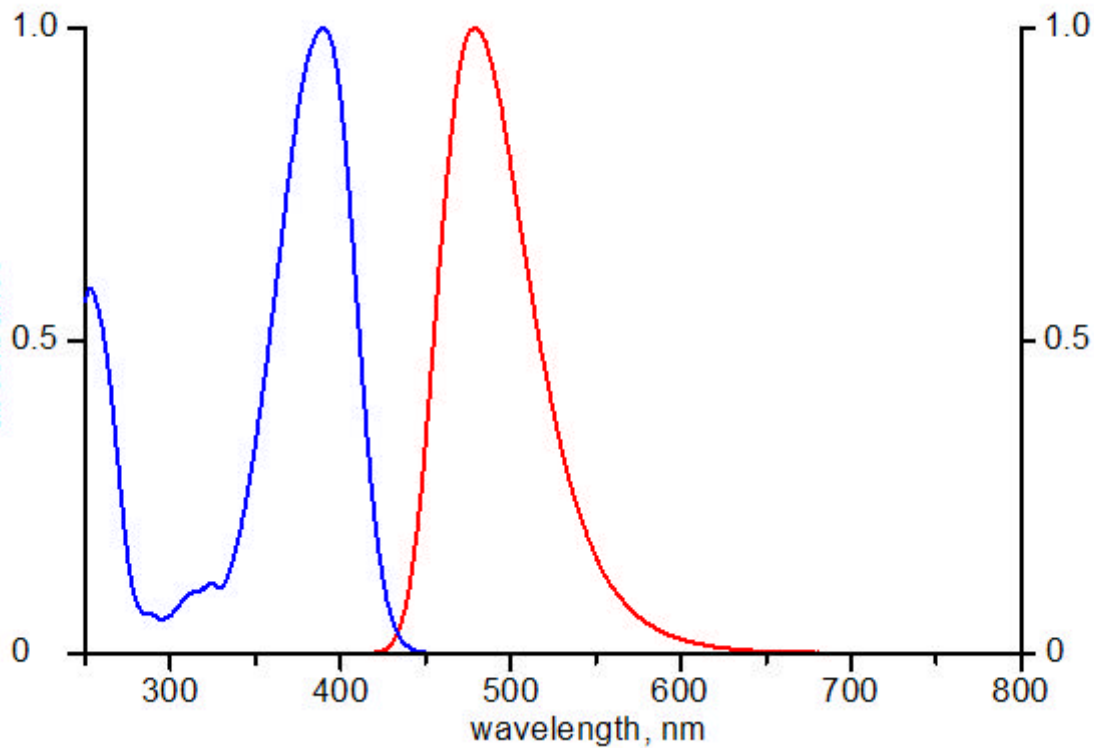


absorbance



fluorescence