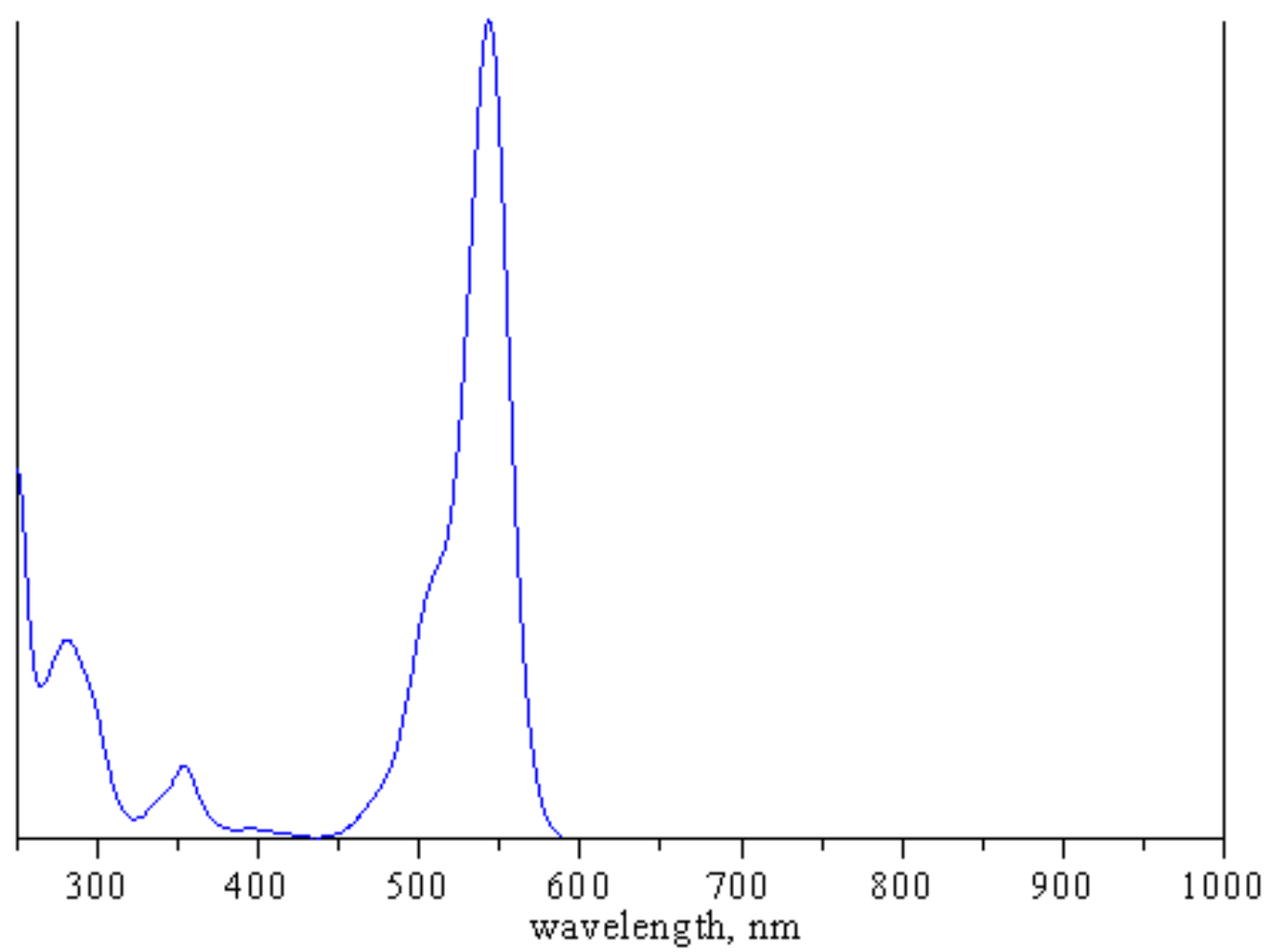


absorbance



wavelength, nm