

absorbance

