

absorbance / intensity

