

# COMPARING Quantitation Methods

## UV Absorbance w/ Spectrophotometer

Quick, easy and cost-effective



$A_{260}$  of 1 =

5 $\mu$ g/ml dsDNA

40 $\mu$ g/ml RNA

33 $\mu$ g/ml ssDNA

CALCULATE

COST EFFECTIVE:



TIME EFFICIENT:



EASY TO USE:



SENSITIVE:



SPECIFIC:

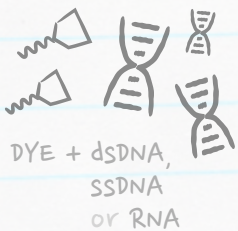


BEST USES:

Cloning

## Fluorescent Dye

Simple, sensitive and nucleic acid-specific



COST EFFECTIVE:



TIME EFFICIENT:



EASY TO USE:



SENSITIVE:



SPECIFIC:

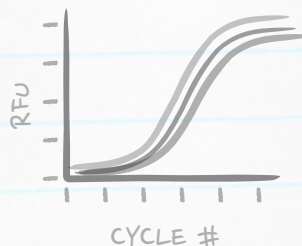
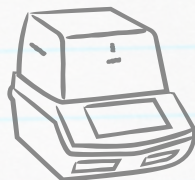
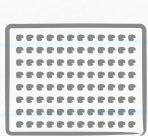


BEST USES:

NGS / STR

## Real-Time Quantitative PCR

Sensitive, specific, great for multiple targets



COST EFFECTIVE:



TIME EFFICIENT:



EASY TO USE:



SENSITIVE:



SPECIFIC:



BEST USES:

NGS / STR