

DNA & RNA Quantitation

DNA Quantitation



Assays for plate reader

Detection limits for microplate reader assays

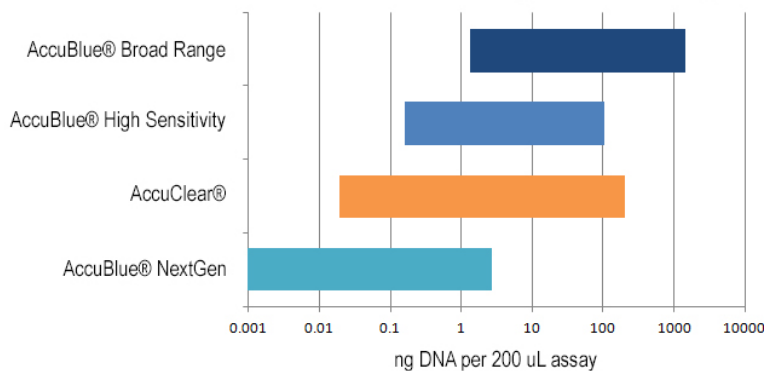


Figure 1. Chart comparing the linear detection ranges for each dsDNA Quantitation Kit for use with fluorescence plate readers. The units are ng of dsDNA per 200 uL assay well. Shown in log scale.

AccuBlue® Broad Range

- Linear range: 2-2000 ng dsDNA
- Best assay for high-concentration DNA samples, such as minipreps and maxipreps

AccuBlue® High Sensitivity

- Linear range: 0.2-100 ng dsDNA
- Membrane-impermeant dye is non-toxic and non-mutagenic, for safer handling and easy disposal

AccuClear® Ultra High Sensitivity

- Widest linear range: 0.03-250 ng dsDNA
- Best choice for most typical DNA samples

AccuBlue® NextGen

- Linear range: 5-3000 pg dsDNA
- The most sensitive and accurate dsDNA quantitation reagent on the market
- Ideal for quantitation of precious samples

Assays for Qubit® fluorometer

Detection limits for Qubit® assays

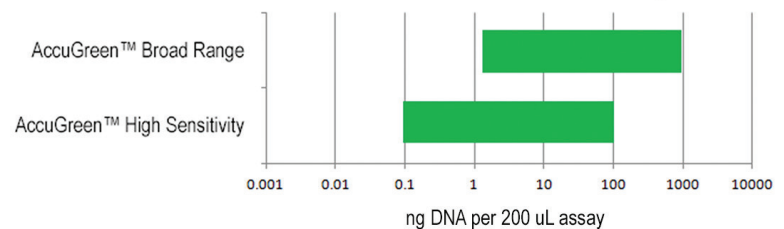


Figure 2. Chart comparing the linear detection ranges for each dsDNA Quantitation Kit for use with the Qubit® fluorometer. The units are ng of dsDNA per 200 uL assay tube. Shown in log scale.

AccuGreen™ (for Qubit®) advantages:

- Precision quantitation of dsDNA
- For use in the pre-programmed assays on the Qubit® fluorometer
- Lower-cost replacements for Qubit® kits

AccuGreen™ Broad Range

- Linear range: 2-1000 ng dsDNA
- Direct replacement for the Qubit® dsDNA BR Assay Kit

AccuGreen™ High Sensitivity

- Linear range: 0.1-100 ng dsDNA
- Direct replacement for the Qubit® dsDNA HS Assay Kit

RNA Quantitation



AccuBlue® RNA Quantitation Kit: for plate reader or Qubit® fluorometer

- Highly specific for RNA over dsDNA
- Superior sensitivity: down to 5 ng on plate reader
- Linear range of 5-1000 ng RNA
- Optimized for use in the Qubit® preprogrammed RNA Broad Range Assay

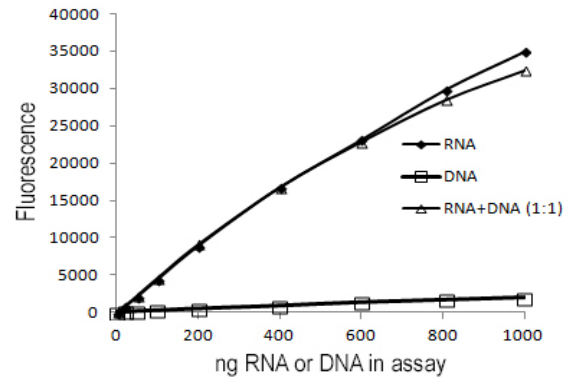


Figure 3. The AccuBlue® Broad Range RNA Quantitation Kit is highly specific for RNA versus dsDNA. The assay can tolerate equimolar contamination of dsDNA in the sample.

RNA Quantitation Kit Ordering Information

Linear Range	Description	Cat. #	Instrument
Broad Range 5-1000 ng	AccuBlue® Broad Range RNA Quantitation Kit	31073, 31073-T	Plate reader & Qubit®

DNA Quantitation Kits Ordering Information

Linear Range	Description	Cat. #	Instrument
Most Sensitive 5-3000 pg	AccuBlue® NextGen dsDNA Quantitation Kit	31060, 31060-T	Plate reader
Ultra High Sensitivity 0.03-250 ng	AccuClear® Ultra High Sensitivity dsDNA Quantitation Kit	31028, 31029	Plate reader
High Sensitivity 0.2-100 ng	AccuBlue® High Sensitivity dsDNA Quantitation Kit	31006, 31006-T	Plate reader
Broad Range 2-2000 ng	AccuBlue® Broad Range dsDNA Quantitation Kit	31007, 31007-T	Plate reader
High Sensitivity 0.1-100 ng	AccuGreen™ High Sensitivity dsDNA Quantitation Kit (for Qubit®)	31066, 31066-T	Qubit®
Broad Range 2-1000 ng	AccuGreen™ Broad Range dsDNA Quantitation Kit (for Qubit®)	31069, 31069-T	Qubit®

DNA Standard Ordering Information

Standard Type	Description	Cat. #	Amount
dsDNA standard sets	AccuBlue® Broad Range dsDNA Standards, Set of Nine (0-200 ng/uL dsDNA)	31007C	0.5 mL each
	AccuBlue® High Sensitivity dsDNA Standards, Set of Eight (0-10 ng/uL dsDNA)	31006C	0.5 mL each
dsDNA single standards	dsDNA Standard, 10 ng/uL	99820-T	1 mL
	dsDNA Standard, 25 ng/uL	31029C	1 mL
	dsDNA Standard, 100 ng/uL	99838-T	1 mL
RNase-free dsDNA	RNase-Free Calf Thymus DNA, 1 mg/mL	31065	1 mL