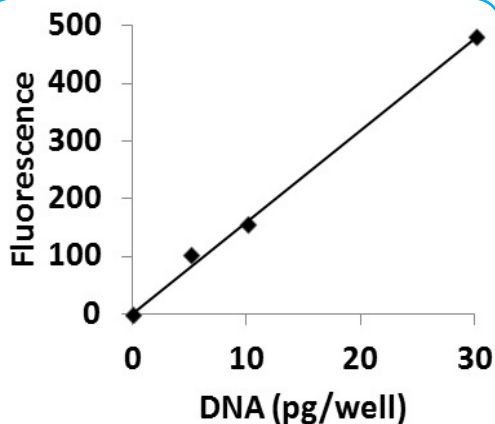


DNA Quantitation

Kits for plate reader or Qubit® reader



Superior DNA quantitation kits for any instrument

AccuBlue®, AccuClear® and AccuGreen™ dsDNA quantitation assays allow precise quantitation of purified dsDNA samples across a wide range of concentrations. Unlike absorbance-based nucleic acid quantitation, fluorescent DNA binding dyes are highly sensitive and selective for double-stranded DNA and provide a more accurate DNA concentration in the presence of contaminating RNA and other common contaminants including free nucleotides, protein, detergents and salts. Biotium offers dsDNA quantitation kits for different sample concentration ranges (see Figure 1).

AccuBlue® and AccuClear® advantages:

- Wide dynamic ranges
- Highly specific for dsDNA
- Compatible with popular plate reader instruments
- Best sensitivity on the market for extremely low-concentration DNA (good for NGS)
- Long working solution stability

Detection limits for microplate reader assays

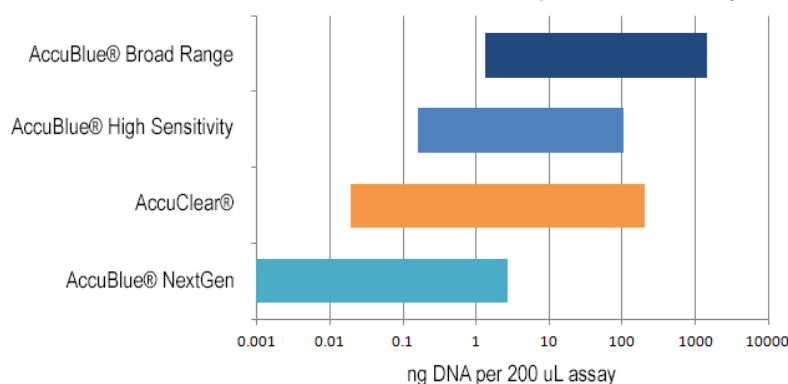


Figure 1. Chart comparing the linear detection ranges for each dsDNA Quantitation Kit for use with fluorescent plate readers. The units are ng of dsDNA per 200 uL reaction. Shown in log scale.

AccuBlue® NextGen

- Linear range: 5 pg-3 ng dsDNA
- Green fluorescence (Ex/Em: 468/507 nm)
- The most sensitive and accurate dsDNA quantitation reagent on the market
- Ideal for quantitation of low-concentration or precious samples

AccuBlue® High Sensitivity

- Linear range: 0.2-100 ng dsDNA
- Green fluorescence (Ex/Em: 500/530 nm)
- Membrane-impermeable dye is non-toxic and non-mutagenic, for safer handling and easy disposal
- Best assay for accurate quantitation of short DNA fragments

AccuClear® Ultra High Sensitivity

- Widest linear range: 0.03-250 ng dsDNA
- Green fluorescence (Ex/Em: 468/507 nm)
- Best choice for most typical DNA samples

AccuBlue® Broad Range

- Linear range: 2-2000 ng dsDNA
- Blue fluorescence (Ex/Em: 350/460 nm)
- Best assay for high-concentration DNA samples, such as minipreps and maxipreps

AccuGreen™ assays for the Qubit® fluorometer

AccuGreen™ for Qubit advantages:

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



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

AccuGreen™ High Sensitivity

- Linear range: 0.1-100 ng dsDNA
- Green fluorescence (Ex/Em: 509/530 nm)
- Direct replacement for the Qubit® dsDNA HS Assay Kit

AccuGreen™ Broad Range

- Linear range: 2-1000 ng dsDNA
- Green fluorescence (Ex/Em: 500/530 nm)
- Direct replacement for the Qubit® dsDNA BR Assay Kit

Ordering Information

Linear Range	Description	Cat. #	Size
Most Sensitive 2.5-3000 pg	AccuBlue® NextGen dsDNA Quantitation Kit with 1 DNA Standard	31060, 31060-T	1000, 200 assays
	AccuBlue® NextGen dsDNA Quantitation Solution	31061, 31061-T	1000, 200 assays
Ultra High Sensitivity 0.03-250 ng	AccuClear® Ultra High Sensitivity dsDNA Quantitation Kit with 7 DNA Standards	31028	1000 Assays
	AccuClear® Ultra High Sensitivity dsDNA Quantitation Kit with 1 DNA Standard	31029	4000 Assays
	AccuClear® Ultra High Sensitivity dsDNA Quantitation Solution	31027, 31027-T	1000, 200 assays
High Sensitivity 0.2-100 ng	AccuBlue® High Sensitivity dsDNA Quantitation Kit with DNA Standard, Trial Size	31006-T	200 Assays
	AccuBlue® High Sensitivity dsDNA Quantitation Kit with 8 DNA Standards	31006	1000 Assays
	AccuBlue® High Sensitivity dsDNA Quantitation Solution	31008, 31008-T	1000, 200 assays
Broad Range 2-2000 ng	AccuBlue® Broad Range dsDNA Quantitation Kit with DNA Standard, Trial Size	31007-T	200 Assays
	AccuBlue® Broad Range dsDNA Quantitation Kit with 9 DNA Standards	31007	1000 Assays
	AccuBlue® Broad Range dsDNA Quantitation Solution	31009, 31009-T	1000, 200 assays
High Sensitivity (for Qubit®) 0.1-100 ng	AccuGreen™ High Sensitivity dsDNA Quantitation Kit (for Qubit®)	31066, 31066-T	500, 100 assays
	AccuGreen™ High Sensitivity dsDNA Quantitation Solution (for Qubit®)	31068, 31068-T	500, 100 assays
Broad Range (for Qubit®) 2-1000 ng	AccuGreen™ Broad Range dsDNA Quantitation Kit (for Qubit®)	31069, 31069-T	500, 100 assays
	AccuGreen™ Broad Range dsDNA Quantitation Solution (for Qubit®)	31070, 31070-T	500, 100 assays
Stand-alone dsDNA standard sets	AccuBlue® Broad Range dsDNA Standards, Set of Nine (0-200 ng/uL calf thymus dsDNA)	31007C	0.5 mL each
	AccuBlue® High Sensitivity dsDNA Standards, Set of Eight (0-10 ng/uL calf thymus dsDNA)	31006C	0.5 mL each