Glowing products for science

AccuBlue® NextGen dsDNA Quantitation Kit

A fluorescent dsDNA quantitation kit with a linear range of 1-3,000 pg DNA, an unprecedented level of sensitivity in dsDNA detection. Designed for use in a fluorescence microplate reader.

Section Sectio

Product attributes

Excitation/Emission 468/507 nm (with DNA)

Storage Conditions Store at 2 to 8 °C

Call us: 800-304-5357 Email: btinfo@biotium.com

Product Description

The AccuBlue® NextGen dsDNA Quantitation assay is the most sensitive and accurate DNA quantitation method for ultra low amounts of DNA, with a linear range from ~1 – 3,000 pg (depending on instrument capability). This assay is designed for use with fluorescence microplate readers. The AccuBlue® NextGen assay is also available without any DNA standards (31061), for those that wish to use their own standard.

With a detection limit as low as 1 pg DNA in a 96-well plate reader assay (see Table 1), AccuBlue® NextGen outperforms rivals such as PicoGreen® in sensitivity and accuracy. Specially designed for scientists who need to quantify extremely dilute or precious DNA, AccuBlue® NextGen is ideal for use in preparing samples for applications such as Next Generation Sequencing (NGS) and digital PCR.

Because the AccuBlue® NextGen assay is so sensitive, results will vary depending on many factors, such as microplate reader (see table below), assay volume, and pipet accuracy.

Plate Reader Detection Limits

Plate reader	Assay format (assay volume)	Lowest DNA detected
BMG CLARIOStar	96-well (200 uL)	1 pg
BMG CLARIOStar	384-well (50 uL)	0.6 pg
Biotek Synergy H1m	96-well (200 uL)	5 pg
SpectraMax Gemini XS	96-well (200 uL)	5 pg
SpectraMax i3	384-well (50 uL)	1.25 pg

Fluorescence-Based dsDNA Quantitation

AccuBlue®, AccuClear® and AccuGreen™ dsDNA quantitation assays allow precise quantitation of purified dsDNA samples across a wide range of concentrations and a variety of fluor 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 concentration in the concentration in the presence of contaminating RNA and other concentration in the concen Biotium offers dsDNA quantitiation kits and solutions for different instruments and sample concentration ranges. See the table below and visit the DNA & RNA Quantitation Technology P

All DNA & RNA Quantitation Kits

Kit	DNA or RNA	Detection range (in assay)*	Dye Ex/Em (nm)	Suggested instrument	Features
AccuGreen™ High Sensitivity DNA	DNA	0.1-100 ng	502/523	Qubit® fluorometer	Compare to the Qubit® dsDNA HS assay from Thermo Fisher
AccuGreen™ Broad Range DNA	DNA	2-1000 ng	500/530	Qubit® fluorometer	Compare to the Qubit® dsDNA BR assay from Thermo Fisher Non-toxic & non-mutagentic DNA quantitation dye
AccuBlue® NextGen DNA	DNA	1-3000 pg**	468/507	Fluorescence microplate reader	Most sensitive assay available for quantitation of precious or dilute samples Optimal for sensitive applications such as NGS or digital PCR
AccuClear® Ultra High Sensitivity DNA	DNA	0.03-250 ng	468/507	Fluorescence microplate reader	Versatile kit with high sensitivity and wide linear range
AccuBlue® High Sensitivity DNA	DNA	0.2-100 ng	485/530	Fluorescence microplate reader	Non-toxic & non-mutagentic DNA quantitation dye
AccuBlue® Broad Range DNA	DNA	2-2000 ng	350/460	Fluorescence microplate reader	Broad linear range with blue fluorescence
AccuBlue® Broad Range RNA	RNA	5-1000 ng	650/670	Fluorescence microplate reader or Qubit® fluorometer	The widest linear range of available RNA quantitation kits Exceptional accuracy, sensitivity, and high RNA selectivity

References

- 1. Analytical Biochemistry (2016) 15 Oct pp74-79 doi:10.1016/j.ab.2016.07.022
- 2. J Assist Reprod Genet (2018) 35, 1497-1501 doi:10.1007/s10815-018-1223-4

This datasheet was generated on November 2, 2025 at 03:43:02 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/accublue-nextgen-dsdna-quantitation-kit/

^{*} Standard assay volume is 200 uL

** AccuBlue® NextGen limit of detection is in the range of 1 pg to 5 pg depending on instrument sensitivityAccuBlue and AccuClear are registered trademarks of Biotium, Inc. AccuBlue,