



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

## AccuBlue® High Sensitivity dsDNA Quantitation Solution

A fluorescent dsDNA quantitation kit with a linear range of 0.2-100 ng DNA. Designed for use in a fluorescence microplate reader.



### Product attributes

Excitation/Emission	500/530 nm (with DNA)
Storage Conditions	Store at 2 to 8 °C

### Product Description

The AccuBlue® High Sensitivity dsDNA Quantitation Solution offers sensitive and selective detection of purified dsDNA samples with a linear detection range from 0.2 – 100 ng. This solution does not come with a DNA standard; see the AccuBlue® High Sensitivity Kit ([31006](#)), which contains a set of calf thymus DNA standards.

The AccuBlue® High Sensitivity dsDNA Quantitation Assay uses non-toxic, non-mutagenic EvaGreen® dye (download the [EvaGreen Safety Report](#) for more information), making it safer than PicoGreen® or Quant-iT® kits.

### 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 fluorescence detection instruments. 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 and solutions for different instruments and sample concentration ranges. See the table below and visit the [DNA & RNA Quantitation Technology Page](#) for details on our full line of dsDNA and RNA quantitation kits.

# All DNA & RNA Quantitation Kits

Kit	DNA or RNA	Detection range (in assay)*	Dye Ex/Em (nm)	Suggested instrument	Features
<a href="#">AccuGreen™ High Sensitivity DNA</a>	DNA	0.1-100 ng	502/523	Qubit® fluorometer	Compare to the Qubit® dsDNA HS assay from Thermo Fisher
<a href="#">AccuGreen™ Broad Range DNA</a>	DNA	2-1000 ng	500/530	Qubit® fluorometer	Compare to the Qubit® dsDNA BR assay from Thermo Fisher Non-toxic & non-mutagenic DNA quantitation dye
<a href="#">AccuBlue® NextGen DNA</a>	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
<a href="#">AccuClear® Ultra High Sensitivity DNA</a>	DNA	0.03-250 ng	468/507	Fluorescence microplate reader	Versatile kit with high sensitivity and wide linear range
<a href="#">AccuBlue® High Sensitivity DNA</a>	DNA	0.2-100 ng	485/530	Fluorescence microplate reader	Non-toxic & non-mutagenic DNA quantitation dye
<a href="#">AccuBlue® Broad Range DNA</a>	DNA	2-2000 ng	350/460	Fluorescence microplate reader	Broad linear range with blue fluorescence
<a href="#">AccuBlue® Broad Range RNA</a>	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

\* Standard assay volume is 200 uL  
\*\* AccuBlue® NextGen limit of detection is in the range of 1 pg to 5 pg depending on instrument sensitivity  
AccuBlue and AccuClear are registered trademarks of Biotium, Inc. AccuBlue,

## References

1. Conservation Genet Resour (2011), 3, 753-755 [doi: 10.1007/s12686-011-9450-3](#)
2. Parasites & Vectors (2012), 5:59 [doi:10.1186/1756-3305-5-59](#)
3. PLoS ONE (2013), 8(1) [doi:10.1371/journal.pone.0054743](#)
4. Journal of Chromatography A (2013), [doi:10.1016/j.chroma.2013.11.039](#)
5. Peer J (2013), Nov 19;1:e203 [doi: 10.7717/peerj.203](#)
6. J Biomech (2014), 47(2), 518-525 [doi:10.1016/j.jbiomech.2013.10.036](#)
7. Analytical Biochemistry (2014), [http://dx.doi.org/10.1016/j.ab.2014.03.005](#)
8. J. Bacteriol. (2014), Jul;196(13)2355-66 [doi:10.1128/JB.01493-14](#)
9. Applications in Plant Sciences (2015), [doi: 10.3732/apps.1400077](#)
10. Conserv Genet (2016), 17, 337-353 [doi: 10.1007/s10592-015-0784-3](#)
11. Invertebrate Biology (2015), 134(3) pp 252-259 [doi: 10.1111/ivb.12091](#)
12. Clin Oral Invest (2016) Jan;21(1) 447-452 [doi: 10.1007/s00784-016-1811-6](#)
13. Pharm res (2018), 35, 188 [doi 10.1007/s11095-018-2460-z](#)

This datasheet was generated on January 22, 2026 at 12:37:11 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/accublu-high-sensitivity-dsdna-quantitation-solution/>