Glowing products for science

AccuBlue® High Sensitivity dsDNA Quantitation Kit with DNA Standards

The state of the s

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

Product Description

The AccuBlue® High Sensitivity dsDNA Quantitation Kit offers sensitive and selective detection of purified dsDNA samples with a linear detection range from 0.2 – 100 ng. The AccuBlue® High Sensitivity assay is also available without any DNA standards (31008), for those that wish to use their own standard.

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 quantitiation 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.

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

Product attributes

Excitation/Emission 500/530 nm (with DNA)

Storage Conditions Store at 2 to 8 °C

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

^{*} Standard assay volume is 200 uL

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 November 8, 2025 at 04:51:56 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/accublue-high-sensitivity-dsdna-quantitation-kit-with-8-dna-standards/

^{**} 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,