Glowing products for science

# AccuGreen™ Broad Range dsDNA Quantitation Kit

The state of the s

A fluorescent dsDNA quantitation kit with a linear range of 2-1000 ng DNA. Designed for use with handheld fluorometers such as the Qubit® fluorometer from Thermo Fisher Scientific.

## **Product Description**

The AccuGreen™ Broad Range dsDNA Quantitation Kit offers sensitive and selective detection of purified dsDNA samples. The kit provides a linear detection range from 2 – 1000 ng and is designed for use with handheld fluorometers such as the Qubit® fluorometers from Thermo Fisher Scientific. The AccuGreen™ Broad Range assay is also available without a DNA standard (31070), for those that wish to use their own standard.

### AccuGreen™ for the Qubit® fluorometer

The AccuGreen™ Broad Range dsDNA Quantitation Kit is designed for use with handheld fluorometers such as the Qubit® fluorometers from Thermo Fisher Scientific.

The linear range of the AccuGreen™ BR assay is 2 to 1000 ng of DNA per tube. DNA samples with concentrations between 0.1 and 1000 ng/uL may be quantified using sample volumes between 1 and 20 uL (for example, 1 uL of 1000 ng/uL is 1000 ng total, and 20 uL of 0.1 ng/uL is 2 ng total, which will both fall within the linear range of the assay). If you use the most common sample volume of 10 uL, the starting sample concentration should be 0.2 to 100 ng/uL.

#### 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 <u>DNA & RNA Quantitation Technology Page</u> for details on our full line of dsDNA and RNA quantitation kits.

#### **Product attributes**

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

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

## **All DNA & RNA Quantitation Kits**

| Kit                                      | DNA<br>or RNA | Detection range (in assay)* | Dye Ex/Em (nm) | Suggested instrument                                 | Features  |
|--|---------------|-----------------------------|----------------|--|---|
| AccuGreen™<br>High Sensitivity DNA       | DNA           | 0.1-100 ng                  | 502/523        | Qubit® fluorometer                                   | Compare to the Qubit® dsDNA HS assay from Thermo Fisher   |
| AccuGreen™ Broad<br>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<br>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<br>Sensitivity DNA | DNA           | 0.03-250 ng                 | 468/507        | Fluorescence microplate reader                       | Versatile kit with high sensitivity and wide linear range   |
| AccuBlue® High<br>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<br>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                      |

This datasheet was generated on November 15, 2025 at 03:38:29 AM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/accugreen-broad-range-dsdna-quantitation-kit/">https://biotium.com/product/accugreen-broad-range-dsdna-quantitation-kit/</a>

<sup>\*</sup> 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,