Glowing products for science

AccuBlue® Broad Range RNA Quantitation Kit

An RNA-selective, fluorescent RNA quantitation kit with the widest linear range (5 ng-1000 ng RNA) of any RNA quantitation assay. Designed for use in a microplate reader or a Qubit® fluorometer.

Product attributes

Call us: 800-304-5357

Excitation/Emission	650/670 nm
Shipping condition	Ice pack
Storage Conditions	See individual components for storage temperature

Email: btinfo@biotium.com

Product Description

The AccuBlue® Broad Range RNA Quantitation Kit boasts the widest linear range (5 ng–1000 ng RNA) of available RNA quantitation kits, while providing exceptional accuracy, sensitivity, and high RNA selectivity. A mammalian RNA standard is included for accurate quantitation of mammalian cell RNA.

Instrument compatibility

The AccuBlue® Broad Range RNA Quantitation assay is designed for use with fluorescence microplate readers equipped with excitation and emission filters for detecting far-red fluorescence. It is also optimized for use in the Qubit® fluorometer, using the pre-programmed RNA Broad Range program. Certain microplate readers are sensitive enough to extend the linear range down to 1 ng, or sample concentrations of 0.1 ng/uL. See product information sheet for more details.

AccuBlue and AccuClear are registered trademarks of Biotium, Inc. AccuBlue, AccuClear and AccuGreen technologies are covered by granted U.S. and international patents. Qubit and Quant-iT are registered trademarks of Thermo Fisher Scientific.

This datasheet was generated on November 17, 2025 at 12:58:10 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/accublue-broad-range-rna-quantitation-kit/