



Glowing products for science

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

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

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

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 January 17, 2026 at 03:52:32 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/accubroad-range-rna-quantitation-kit/>