

Revised: February 19, 2021

Product Information

RNA Broad Range Standard, 100 ng/uL

Catalog Number: 99850

Unit Size: 300 uL

Storage and Handling

The RNA standard should be stored at -20°C or below. Storage at -80°C will extend the shelf life of the RNA standard. The RNA is stable for 6 months from date of receipt when stored as recommended.

Product Description

The RNA Broad Range Standard, 100 ng/uL is purified total mammalian RNA provided in TE buffer. It intended for use in the AccuBlue® Broad Range RNA Quantitation Kit (Cat. # 31073). This RNA standard is identical to the one supplied in the AccuBlue® RNA quantitation kit, but is offered as a stand-alone replacement item in the case that the RNA standard in your kit becomes degraded or otherwise compromised.

The AccuBlue® Broad Range RNA Quantitation Kit has high RNA selectivity and has an extremely wide linear range (5-1000 ng), enabling both high-sensitivity and broad-range quantitation with a single kit. It is optimized for use in fluorescent microplate readers and on Qubit® fluorometers.

Related Products

Catalog number	Product
31073	AccuBlue® Broad Range RNA Quantitation Kit
31065	RNase-Free Calf Thymus DNA, 1 mg/mL
CD201	RNAstorm™ Kit for Isolation of RNA from FFPE Tissue Samples
CD504	RNAstorm™ RNA Isolation Kit
31028	AccuClear® Ultra High Sensitivity dsDNA Quantitation Kit
31066	AccuGreen™ High Sensitivity dsDNA Quantitation Kit (for Qubit®)
41024	Water, Ultrapure Molecular Biology Grade (RNase-, DNase-Free)
41003	GelRed® Nucleic Acid Gel Stain

Please visit our website at www.biotium.com for information on our life science research products including DNA quantitation kits, DNA and RNA extraction kits, environmentally friendly EvaGreen® qPCR master mixes, and other reagents for molecular and cell biology research.

Qubit is a registered trademark of Thermo Fisher.

AccuBlue, AccuClear and AccuGreen are covered by granted and/or pending U.S. and/or international patents.

Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use.