



Glowing products for science

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

RNA Broad Range Standard, 100 ng/uL

Stand-alone total RNA standard provided with the AccuBlue® Broad Range RNA Quantitation Kit (31073).



Product attributes

Storage Conditions	Store at -10 to -35 °C or below
Shipping condition	Ice pack

Product Description

Stand-alone total mammalian RNA standard provided with the AccuBlue® Broad Range RNA Quantitation Kit (31073).

- Mammalian cell total RNA in 1X Tris-EDTA
- 100 ng/uL
- 300 uL
- Store at -20 °C or below (can be stored at -80 °C for long-term stability)

This datasheet was generated on January 8, 2026 at 12:21:03 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/rna-broad-range-standard-100-ng-ul/>