

Measuring the QuantiFluor™ RNA System Using the QuantiFluor™-ST Fluorometer

ABSTRACT

This Application Note describes the protocol for using the QuantiFluor™ RNA System with the QuantiFluor™-ST Fluorometer and the PCR Tube Adapter to detect and quantitate small amounts of RNA.

Publication Date: 11/12.



Download Article

(153 KB)
