



SpectraMax® Quant™

AccuBlue™ Pico dsDNA Assay Kit

Quick Start Protocol for a 96-Well Microplate

1. Warm assay components to room temperature.
2. Prepare the working solution by diluting the AccuBlue Pico dye (1:400) and AccuBlue Pico Enhancer (1:100) in 1X AccuBlue Pico buffer and mix.
3. Pipette 200 μ L of the working solution into each microplate well.
4. Pipette 10 μ L of the SpectraMax Quant AccuBlue Pico dsDNA standards into the microplate wells and mix.
5. Pipette 10 μ L of each unknown DNA sample into separate microplate wells and mix.
6. Incubate the microplate at room temperature for 5 to 10 minutes in the dark.
7. Measure fluorescence in a fluorescence microplate reader (EX=468 nm, EM=507 nm).

R3648 A
June 2016



Table 1-1: Preparation of dsDNA Standards

Standard	Final Concentration	Standard Volume	1X AccuBlue Pico Buffer Volume
A	300 pg/ μ L	AccuBlue Pico Standard	None
B	50 pg/ μ L	10 μ L of A	50 μ L
C	10 pg/ μ L	12 μ L of B	48 μ L
D	2 pg/ μ L	12 μ L of C	48 μ L
E	0.5 pg/ μ L	15 μ L of D	45 μ L
F	0 pg/ μ L	None	60 μ L

Complete instructions for using the SpectraMax Quant AccuBlue Pico dsDNA Assay Kit can be downloaded from the Knowledge Base on the Molecular Devices support web site: www.moleculardevices.com/support

Click **Knowledge Base**, and then click **Manuals** and search for **SpectraMax Quant AccuBlue Pico dsDNA**.

Contact Us

Phone: +1-800-635-5577

Web: moleculardevices.com

Email: info@moldev.com

Visit our website for a current listing of worldwide distributors.



The trademarks used herein are the property of Molecular Devices, LLC or their respective owners.

Specifications subject to change without notice.

Patents: www.moleculardevices.com/productpatents

FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES

© 2016 Molecular Devices, LLC.

All rights reserved.

R3648 A