



# 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  $\mu$ L of the working solution into each microplate well.
4. Pipette 10  $\mu$ L of the SpectraMax Quant AccuBlue Pico dsDNA standards into the microplate wells and mix.
5. Pipette 10  $\mu$ 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/ $\mu$ L     | AccuBlue Pico Standard | None                           |
| B        | 50 pg/ $\mu$ L      | 10 $\mu$ L of A        | 50 $\mu$ L                     |
| C        | 10 pg/ $\mu$ L      | 12 $\mu$ L of B        | 48 $\mu$ L                     |
| D        | 2 pg/ $\mu$ L       | 12 $\mu$ L of C        | 48 $\mu$ L                     |
| E        | 0.5 pg/ $\mu$ L     | 15 $\mu$ L of D        | 45 $\mu$ L                     |
| F        | 0 pg/ $\mu$ L       | None                   | 60 $\mu$ 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](http://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](http://moleculardevices.com)  
Email: [info@moldev.com](mailto: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](http://www.moleculardevices.com/productpatents)  
FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES

© 2016 Molecular Devices, LLC.  
All rights reserved.  
R3648 A