



SpectraMax® Quant™

AccuBlue™ HiRange 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 HiRange dye (1:100) and AccuBlue HiRange Enhancer (1:100) in 1X AccuBlue HiRange buffer and mix.
3. Pipette 200 μL of the working solution into each microplate well.
4. Pipette 10 μL of the SpectraMax Quant AccuBlue HiRange dsDNA standards into the microplate wells and mix.
5. Pipette 10 μL 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=350 nm, EM=460 nm).

R3649 A
June 2016



Table 1-1: Preparation of dsDNA Standards

| Std. | Final Conc. | Std. Volume | AccuBlue HiRange Buffer Volume |
|------|------------------|--|--------------------------------|
| A | 200 ng/ μ L | 100 μ L of 200 ng/ μ L AccuBlue HiRange Standard | None |
| B | 150 ng/ μ L | 75 μ L of 200 ng/ μ L AccuBlue HiRange Standard | 25 μ L |
| C | 100 ng/ μ L | 100 μ L of 200 ng/ μ L AccuBlue HiRange Standard | 100 μ L |
| D | 50 ng/ μ L | 100 μ L of 100 ng/ μ L of C | 100 μ L |
| E | 25 ng/ μ L | 100 μ L of 50 ng/ μ L of D | 100 μ L |
| F | 12.5 ng/ μ L | 100 μ L of 25 ng/ μ L of E | 100 μ L |
| G | 6.25 ng/ μ L | 100 μ L of 12.5 ng/ μ L of F | 100 μ L |
| H | 2 ng/ μ L | 32 μ L of 12.5 ng/ μ L of G | 68 μ L |
| I | 0.2 ng/ μ L | 10 μ L of 2 ng/ μ L of H | 90 μ L |
| J | 0 ng/ μ L | None | 100 μ L |

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

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.
R3649 A