



SpectraMax® Quant™

AccuClear™ Nano 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 AccuClear Nano dye (1:100) in 1X AccuClear Nano buffer and mix.
3. Pipette 200 μL of the working solution into each microplate well.
4. Pipette 10 μL of the SpectraMax Quant AccuClear Nano 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=468 nm, EM=507 nm).

R3650 A
June 2016



Table 1-1: Preparation of dsDNA Standards

Std.	Final Conc.	Std. Volume	1X AccuClear Nano Buffer Volume
A	25 ng/ μ L	100 μ L of 25 ng/ μ L AccuClear Nano Standard	None
B	10 ng/ μ L	40 μ L of 25 ng/ μ L AccuClear Nano Standard	60 μ L
C	3 ng/ μ L	12 μ L of 25 ng/ μ L AccuClear Nano Standard	88 μ L
D	1 ng/ μ L	10 μ L of 10 ng/ μ L of C	90 μ L
E	0.3 ng/ μ L	10 μ L of 3 ng/ μ L of D	90 μ L
F	0.1 ng/ μ L	10 μ L of 1 ng/ μ L of E	90 μ L
G	0.03 ng/ μ L	10 μ L of 0.3 ng/ μ L of F	90 μ L
H	0.01 ng/ μ L	10 μ L of 0.1 ng/ μ L of G	90 μ L
I	0.003 ng/ μ L	10 μ L of 0.03 ng/ μ L of H	90 μ L
J	0 ng/ μ L	None	100 μ L

Complete instructions for using the SpectraMax Quant AccuClear Nano 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.
R3650 A