



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 μ L of the working solution into each microplate well.
4. Pipette 10 μ L of the SpectraMax Quant AccuClear Nano 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).

R3650 B
November 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 B	90 μ L
E	0.3 ng/ μ L	10 μ L of 3 ng/ μ L of C	90 μ L
F	0.1 ng/ μ L	10 μ L of 1 ng/ μ L of D	90 μ L
G	0.03 ng/ μ L	10 μ L of 0.3 ng/ μ L of E	90 μ L
H	0.01 ng/ μ L	10 μ L of 0.1 ng/ μ L of F	90 μ L
I	0.003 ng/ μ L	10 μ L of 0.03 ng/ μ L of G	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 B