

SpectraMax Quant AccuClear Nano dsDNA Assay Kit

The SpectraMax® Quant™ AccuClear™ Nano dsDNA Assay Kit is a part of the SpectraMax Quant family of double-stranded DNA (dsDNA) assay products. Each product is tailored to a specific range of DNA concentrations, and is optimized for use with fluorescence microplate readers such as the SpectraMax® microplate readers. Preconfigured protocols are included in SoftMax® Pro Software to make DNA quantitation analysis with the assay kits part of a simplified workflow.

The SpectraMax Quant AccuClear Nano dsDNA Assay Kit enables quantitation of a wide range of dsDNA concentrations. The assay is linear between 34 pg per well and 250 ng per well in the 96-well microplate format as determined using a fluorescence microplate reader.

Table 1-1: Available Kits

Assay Kit	Explorer Kit	Bulk Kit
SpectraMax® Quant™ AccuClear™ Nano dsDNA Assay Kit	R8356	R8357

Contents

- Chapter 1: About the SpectraMax Quant AccuClear Nano dsDNA Assay Kit** **3**
 - Assay Principles 3
- Chapter 2: Materials and Equipment** **5**
 - Kit Components 5
 - Storage and Handling 5
 - Materials Required but Not Provided 5
 - Compatible Molecular Devices Microplate Readers 5
- Chapter 3: Assay Protocol** **6**
 - Considerations Before you Begin 6
 - Fluorescence Microplate Reader Setup with SoftMax Pro Software 7
 - Example of Assay Protocol 7
 - Assay Setup 7
- Chapter 4: Data Analysis Example** **10**
 - Example Standard Curve 10
- Chapter 5: Obtaining Support** **11**
 - SpectraMax Quant Product Family 11

SpectraMax Quant AccuClear Nano dsDNA Assay Kit

This document is provided to customers who have purchased Molecular Devices equipment, software, reagents, and consumables to use in the operation of such Molecular Devices equipment, software, reagents, and consumables. This document is copyright protected and any reproduction of this document, in whole or any part, is strictly prohibited, except as Molecular Devices may authorize in writing.

Software that may be described in this document is furnished under a non-transferrable license. It is against the law to copy, modify, or distribute the software on any medium, except as specifically allowed in the license agreement. Furthermore, the license agreement may prohibit the software from being disassembled, reverse engineered, or decompiled for any purpose.

Portions of this document may make reference to other manufacturers and/or their products, which may contain parts whose names are registered as trademarks and/or function as trademarks of their respective owners. Any such usage is intended only to designate those manufacturers' products as supplied by Molecular Devices for incorporation into its equipment and does not imply any right and/or license to use or permit others to use such manufacturers' and/or their product names as trademarks.

Each product is shipped with documentation stating specifications and other technical information. Molecular Devices products are warranted to meet the stated specifications. Molecular Devices makes no other warranties or representations express or implied, including but not limited to, the fitness of this product for any particular purpose and assumes no responsibility or contingent liability, including indirect or consequential damages, for any use to which the purchaser may put the equipment described herein, or for any adverse circumstances arising therefrom. The sole obligation of Molecular Devices and the customer's sole remedy are limited to repair or replacement of the product in the event that the product fails to do as warranted.



For research use only. Not for use in diagnostic procedures.

The trademarks mentioned herein are the property of Molecular Devices, LLC or their respective owners. These trademarks may not be used in any type of promotion or advertising without the prior written permission of Molecular Devices, LLC.

AccuBlue and AccuClear are trademarks of Biotium, Inc.

Molecular Devices Patents: <http://www.moleculardevices.com/productpatents>

Product manufactured for Molecular Devices, LLC
1311 Orleans Drive, Sunnyvale, California, United States of America 94089.
Molecular Devices, LLC is ISO 9001 registered.
©2016 Molecular Devices, LLC.
All rights reserved.

Chapter 1: About the SpectraMax Quant AccuClear Nano dsDNA Assay Kit

The SpectraMax Quant AccuClear Nano dsDNA Assay Kit provides a convenient tool for quantitation of a wide range of concentrations of dsDNA samples. The key features of this kit are the following:

- Optimal for a wide range of concentrations of dsDNA samples
- Highly selective for dsDNA over single-stranded DNA or RNA
- Optimized for fluorescence microplate readers, such as SpectraMax® readers
- Simplified data acquisition and analysis with preconfigured protocol in SoftMax® Pro Software

Assay Principles

The SpectraMax Quant AccuClear Nano dsDNA Assay Kit provides an accurate and sensitive method for quantitation of a wide range of DNA concentrations. The kit contains a fluorescent dye that binds to dsDNA. [Figure 1-1](#) shows the fluorescence excitation peak at 468 nm and the emission peak at 507 nm. [Figure 1-2](#) shows that, unlike absorbance-based measurements, AccuClear Nano dye is highly selective for dsDNA over single-stranded DNA (ssDNA) or RNA. As shown in [Example Standard Curve on page 10](#), the assay is linear between 34 pg and 250 ng of dsDNA per well in a 96-well microplate format using microplate readers such as the SpectraMax® i3x Multi-Mode Microplate Reader.

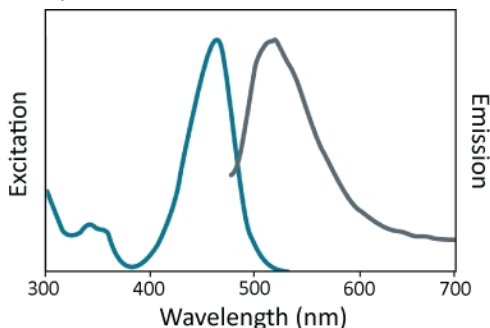


Figure 1-1: SpectraMax Quant AccuClear Nano dsDNA assay dye fluorescent excitation and emission spectra

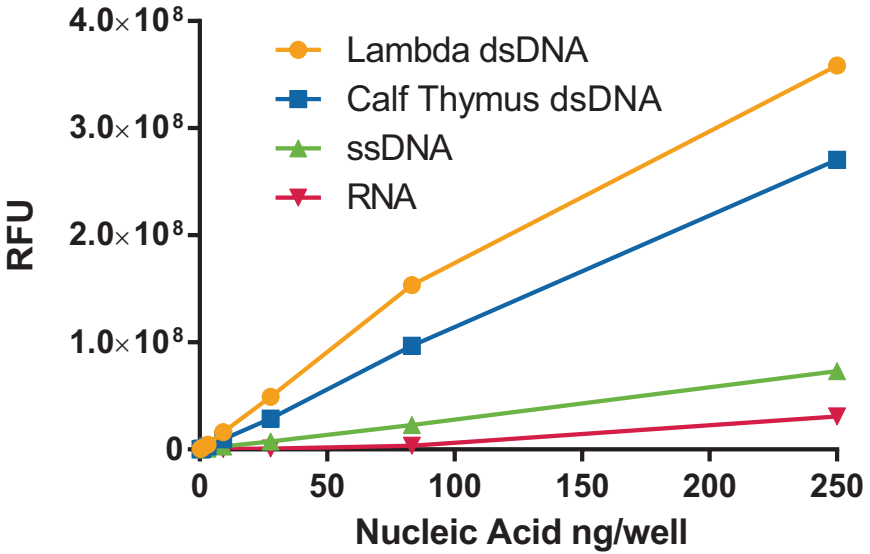


Figure 1-2: SpectraMax Quant AccuClear Nano dsDNA specificity assay

Specificity of SpectraMax Quant AccuClear Nano dsDNA assay is demonstrated by comparing the fluorescent signals collected from the same concentrations of lambda dsDNA, calf thymus dsDNA, ssDNA, and RNA. Plots of relative fluorescence units (RFU) vs. nucleic acid concentration show that the signals are significantly higher for calf thymus dsDNA and lambda dsDNA than for RNA and ssDNA.

Chapter 2: Materials and Equipment

Kit Components

Table 2-1: Components of the SpectraMax Quant AccuClear Nano dsDNA Assay Kit

Item	Explorer Kit (R8356)	Bulk Kit (R8357)
AccuClear Nano Dye, 100X	1 x 0.5 mL	2 x 1 mL
AccuClear Nano Buffer, 20X	2 mL	10 mL
dsDNA standard from calf thymus, 25 ng/μL	1 x 1.0 mL	N/A
dsDNA standards from calf thymus: 0, 0.03, 0.1, 0.3, 1.0, 3.0, 10, and 25 ng/μL (8 vials)	N/A	1 x 0.5 mL each

- The Explorer kit is sufficient for two 96-well microplates.
- The Bulk kit is sufficient for ten 96-well microplates.

The number of microplates is based on the example protocol that is detailed in this document.

Storage and Handling

Store at 4°C. When stored as directed, the kit is stable for at least 6 months from the date it is received.



WARNING! Reagents can contain chemicals that are harmful. Exercise care when handling reagents as described in the related safety data sheet (SDS or MSDS). The safety data sheet is available in the Knowledge Base on the Molecular Devices support web site: www.moleculardevices.com/support

Materials Required but Not Provided

Table 2-2: Reagents and Supplies

Item	Suggested Vendor
Solid black microplates (96-well)	Greiner Bio-One or equivalent
Deionized water for buffer dilution	

Compatible Molecular Devices Microplate Readers

- SpectraMax® i3x Multi-Mode Microplate Reader
- SpectraMax® M2 and M2e Multi-Mode Microplate Readers
- SpectraMax® M3 Multi-Mode Microplate Reader
- SpectraMax® M4 Multi-Mode Microplate Reader
- SpectraMax® M5 and M5e Multi-Mode Microplate Readers
- SpectraMax® Paradigm® Multi-Mode Microplate Reader
- FlexStation® 3 Multi-Mode Microplate Reader
- Gemini™ EM and XPS Microplate Readers
- FilterMax™ F3 and F5 Multi-Mode Microplate Readers

Chapter 3: Assay Protocol

Considerations Before you Begin

Optimal assay conditions for different dsDNA types and microplate formats can vary. Molecular Devices recommends verifying the following to determine the optimal conditions for your assay system.

- Because calf thymus DNA is double-stranded, highly polymerized, and approximately 58% AT (42% GC), it can serve as a reference for most plant and animal DNA. Lambda dsDNA yields similar results. You might want to use a standard similar to your unknown samples in DNA length, structure (for example, linear vs. circular), or GC content. Because GC content varies widely depending on the species, a species-specific standard might be appropriate for bacterial DNA.
- Because of variation among microplate readers, optimization of instrument settings might be needed for best assay performance. Instrument settings and other factors that can affect linearity and the fluorescence signal include the following:
 - Excitation and emission wavelengths and bandwidths
 - Cut-off filters
 - Sensitivity (gain) settings
 - Pipetting accuracy
 - Microplate manufacturer
- Several methods for isolation of DNA from cells or organisms are available. As a result, different types of contaminants can be present in the samples which can cause either an increase or a decrease in the fluorescent signal. Table 3-1 shows the effects of common DNA contaminants on the fluorescent signal in the SpectraMax Quant AccuClear Nano dsDNA assay. Triplicate samples of 250 ng dsDNA per well were assayed in the presence of the contaminants at the indicated final concentrations.

Table 3-1: Effects of common DNA contaminants on the fluorescent signal in the SpectraMax Quant AccuClear Nano dsDNA assay

Compound	Initial concentration in sample	Final concentration in assay (200 μ L)	% Change in signal (+/-)
Sodium chloride	1 M	50 mM	-14%
Magnesium chloride	100 mM	5 mM	-16%
Sodium acetate	600 mM	30 mM	-11%
Ammonium acetate	1 M	50 mM	-14%
Ethanol	20%	1.0%	-21%
Phenol	2%	0.10%	-11%
Chloroform	20%	1%	-34%
SDS	0.20%	0.01%	-31%
SDS	0.02%	0.001%	-9%
Triton X-100	0.20%	0.01%	-36%
Triton X-100	0.02%	0.001%	-20%
Tween-20	0.10%	0.005%	-20%
dNTPs	2 mM	100 μ M	-11%
BSA*	1.0 mg/mL	0.05 mg/mL	-25%

*Not compatible with dsDNA quantitation below 10 ng/well.

Fluorescence Microplate Reader Setup with SoftMax Pro Software

Table 3-2 displays typical fluorescence microplate reader settings. In SoftMax® Pro Software, use the preconfigured SpectraMax Quant AccuClear Nano dsDNA protocol that is available in the protocol library in the software or on the protocol sharing web site (www.softmaxpro.org).

Table 3-2: SpectraMax Quant AccuClear Nano dsDNA assay protocol settings

Parameter	Setting
Read Mode	Fluorescence
Read Type	Endpoint
Wavelengths	Excitation: 468 nm with 9 nm bandwidth Emission: 507 nm with 15 nm bandwidth
PMT and Optics	PMT Gain: Automatic Flashes per read: 10 Read From Top

Example of Assay Protocol

The following method was used for generating a standard curve in a 96-well format.

Assay Setup

Use the following procedure to set up the assay:

1. Allow all components to reach room temperature before use.



Tip: AccuClear Nano dye is provided in DMSO, which can freeze during storage at 4°C. All kit components can be placed in a 37°C water bath for rapid warming. Allow solutions to cool to room temperature before using.

2. To minimize reagent loss in the cap, before removing the required volume, shake or vortex each component well, and then centrifuge vials briefly.
3. AccuClear Nano buffer is supplied at 20X. Dilute the buffer to 1X with deionized water on the day of use.
4. For the Bulk kit, use the dsDNA standards that are provided. To prepare a set of standards for the Explorer kit, dilute the 25 ng/μL dsDNA standard in the 1X AccuClear Nano buffer as shown in [Table 3-3: Preparation of dsDNA Standards on page 8](#).



Note: Prepare the standards fresh the day of the assay. Volumes can be scaled as necessary.

Table 3-3: Preparation of dsDNA Standards

Final Concentration	Standard Volume	1X AccuClear Nano Buffer Volume
25 ng/ μ L	100 μ L of 25 ng/ μ L AccuClear Nano Standard	None
10 ng/ μ L	40 μ L of 25 ng/ μ L AccuClear Nano Standard	60 μ L
3 ng/ μ L	12 μ L of 25 ng/ μ L AccuClear Nano Standard	88 μ L
1 ng/ μ L	10 μ L of 10 ng/ μ L	90 μ L
0.3 ng/ μ L	10 μ L of 3 ng/ μ L	90 μ L
0.1 ng/ μ L	10 μ L of 1 ng/ μ L	90 μ L
0.03 ng/ μ L	10 μ L of 0.3 ng/ μ L	90 μ L
0.01 ng/ μ L	10 μ L of 0.1 ng/ μ L	90 μ L
0.003 ng/ μ L	10 μ L of 0.03 ng/ μ L	90 μ L
0 ng/ μ L	None	100 μ L



Note: When following the assay protocol that is described here, the linear range of the assay is determined to be between 34 pg/well and 250 ng/well. Depending on the microplate reader and assay volume, accuracies of 0.1 pg/well or less might be obtainable.

- Prepare the working solution as follows on the day of the assay: Dilute the AccuClear Nano dye 1:100 in 1X AccuClear Nano buffer in a plastic container (do not use glass), and then vortex or shake to mix well.



Note: Because 200 μ L of working solution is required for each standard and sample that is to be tested, volumes can be scaled as required. Prepare only as much working solution as you plan to use within 24 hours.

- For each dsDNA standard or unknown DNA sample that is to be tested do the following: Add 10 μ L of the standard or sample to a well in a black, 96-well microplate, and then add 200 μ L of the working solution and mix.



Note: To test samples in triplicate, prepare three separate wells for each dsDNA standard and three separate wells for each unknown DNA sample.

- Incubate the microplate at room temperature for 5 minutes in the dark.
- Measure fluorescence in a fluorescence microplate reader with excitation at 468 nm and emission at 507 nm. See [Fluorescence Microplate Reader Setup with SoftMax Pro Software on page 7](#).



Note: In SoftMax[®] Pro Software, use the preconfigured SpectraMax Quant AccuClear Nano dsDNA protocol that is available in the protocol library in the software or on the protocol sharing web site (www.softmaxpro.org).

9. Use the standard curve to calculate the concentrations of the unknown DNA samples. See [Example Standard Curve on page 10](#).
-



Note: Because the fluorescence signal decreases over time after the DNA and dye are combined—an approximately 15% decrease after 3 hours and an approximately 30% decrease after 6 hours—Molecular Devices recommends that you measure fluorescence within 1 hour of setting up the assay.



Note: If the fluorescence of any of the unknown samples exceeds the linear range, then further dilute the sample and use 10 μL of the diluted sample to do the assay. For consistency, use the same volume of sample in all the wells.

Chapter 4: Data Analysis Example

The first step in data analysis is generation of a DNA standard curve that is used to calculate the concentration of unknown DNA samples. A standard curve is generated by plotting the fluorescence values for the DNA standards on the Y-axis and the standard DNA concentrations on the X-axis. A linear curve fit is applied using the drop-down list in the graph section of SoftMax Pro Software. From the standard curve, concentrations of unknown DNA samples are interpolated. A preconfigured assay protocol in the Protocol Library of SoftMax Pro Software enables automatic graphing of standards, and calculation of unknown DNA sample concentrations from the standard curve.

Example Standard Curve

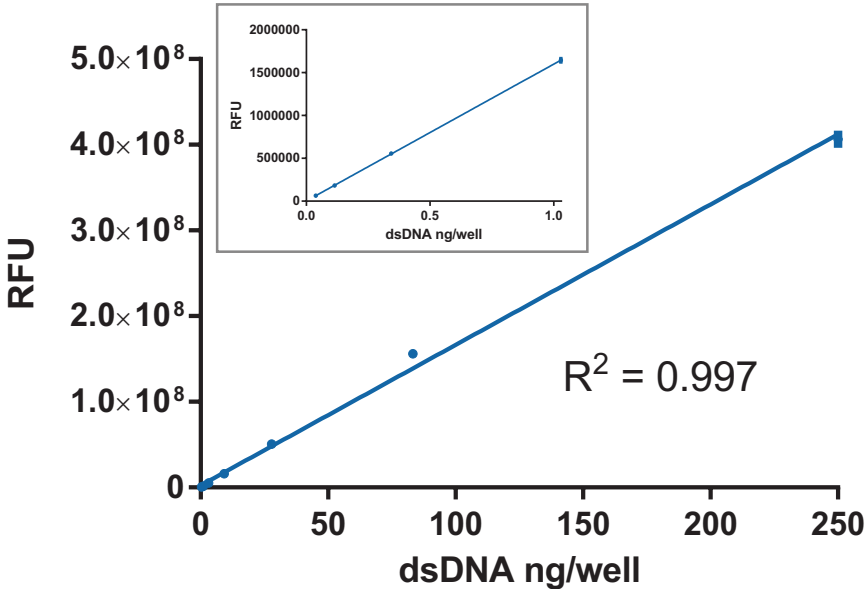


Figure 4-1: SpectraMax Quant AccuClear Nano dsDNA assay standard curve

Linearity of the SpectraMax Quant AccuClear Nano dsDNA assay between 34 ng and 250 ng per well in a 96-well microplate assay with excitation = 468 nm and emission = 507 nm. The inset shows the lower portion of the curve.



Note: The standard curve shown here is for reference only. You must use your own instrument to calculate a standard curve for determining the concentration of DNA in your unknown samples.

Chapter 5: Obtaining Support

Molecular Devices is a leading worldwide manufacturer and distributor of analytical instrumentation, software, and reagents. We are committed to the quality of our products and to fully supporting our customers with the highest possible level of technical service.

Our support web site, www.moleculardevices.com/support, has a link to the Knowledge Base with technical notes, software upgrades, safety data sheets, and other resources. If you do not find the answers you are seeking, follow the links to the Technical Support Service Request Form to send an email message to a pool of technical support representatives.

You can contact your local representative or contact Molecular Devices Technical Support by telephone at 800-635-5577 (U.S. only) or +1 408-747-1700. In Europe call +44 (0) 118 944 8000.

Please have the product name, part number, and lot number available when you call.

SpectraMax Quant Product Family

Table 5-1: SpectraMax Quant Product Family: Available Kits

Assay Kit	Explorer Kit	Bulk Kit
SpectraMax® Quant™ AccuBlue™ Pico dsDNA Assay Kit	R8354	R8355
SpectraMax® Quant™ AccuClear™ Nano dsDNA Assay Kit	R8356	R8357
SpectraMax® Quant™ AccuBlue™ HiRange dsDNA Assay Kit	R8358	R8359

Contact Us

Phone: [800.635.5577](tel:800.635.5577)
Web: www.moleculardevices.com
Email: info@moldev.com

Check our website for a current listing
of worldwide distributors

Regional Offices

USA and Canada
[+1.800.635.5577](tel:+1800.635.5577)
United Kingdom
[+44.118.944.8000](tel:+44.118.944.8000)
Europe*
[00800.665.32860](tel:00800.665.32860)

China (Beijing)
[+86.10.6410.8669](tel:+86.10.6410.8669)
China (Shanghai)
[+86.21.3372.1088](tel:+86.21.3372.1088)
Hong Kong
[+852.2248.6000](tel:+852.2248.6000)

Japan (Osaka)
[+81.6.7174.8831](tel:+81.6.7174.8831)
Japan (Tokyo)
[+81.3.6362.5260](tel:+81.3.6362.5260)
South Korea
[+82.2.3471.9534](tel:+82.2.3471.9534)

Brazil
[+55.11.3616.6607](tel:+55.11.3616.6607)

**Austria, Belgium, Denmark, Finland, France, Germany, Ireland, Netherlands, Spain, Sweden, and Switzerland*