

# Certificate of Analysis

## SpectraMax<sup>®</sup> Quant<sup>™</sup> AccuClear<sup>™</sup> Nano dsDNA Assay Bulk Kit R8357

<b>Lot #</b>	<b>3171190</b>
--------------	----------------

	<b>Specification</b>	<b>This Lot</b>
<b>QC Assay 1 AccuClear dsDNA Quantitation Assay</b>	<b>Standard curve <math>R^2 \geq 0.99</math></b>	<b>1.00</b>

<b>Expiry Date</b>	<b>Up to 6 months from receipt when stored at 2-8 °C</b>
--------------------	--

**This product has been manufactured to ISO9001:2008 Quality Assurance guidelines**

**For research use only Store at 2-8 °C**

**Signed:** *Ervin Ramirez*

**Date:** 19-Jun-2019

**Production Chemist**