

Certificate of Analysis

SpectraMax[®] Quant[™] AccuClear[™] Nano dsDNA Assay Bulk Kit

R8357

Lot #	23337366
--------------	----------

	Specification	This Lot
QC Assay 1 AccuClear dsDNA Quantitation Assay	Standard curve $R^2 \geq 0.99$	1.00

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

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

For research use only Store at 2-8 °C

Signed: *Erwin Ramirez*

Date: 04-Jan-2017

Production Chemist