

Certificate of Analysis

Kit Name:	Simoa™ NF-Light® Kit
Product Number:	102258
Lot Number:	501052
Expiration:	15-Sep-2018

113 Hartwell Ave.  
Lexington MA.  
02421

[www.quanterix.com](http://www.quanterix.com)

Component	Part Number	Lot Number
NF-Light® Bead Reagent	102246	73540
NF-Light® Detector Reagent	102248	73497
NF-Light® SBG Reagent	102250	73491
NF-Light® Sample Diluent	102252	73495
NF-Light® Calibrator Diluent	102620	73492
NF-Light® Calibrator Concentrate	102619	73520
RGP Reagent	103159	N/A <sup>1</sup>

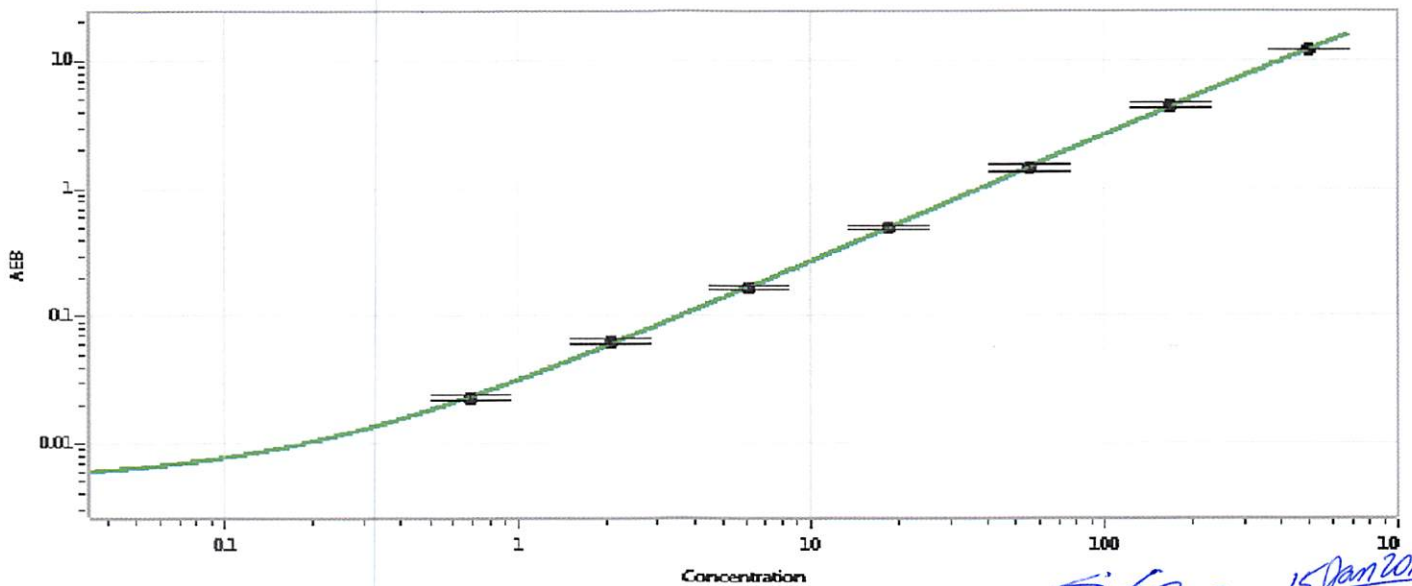
<sup>1</sup> RGP is not Kit Lot Specific

Calibrator Concentrate	Assigned Value (pg/mL)
	7,585

Controls	Lot #s	Range (pg/mL)	Value at Release (pg/mL)
Control 1	72720	3.90-5.86	5.38
Control 2	72720	116-227	197

Patient Panels	Lot Number	Range (pg/mL)	Value at Release (pg/mL)
1	MF-092717-01	4.82-7.28	6.47
2	MF-092717-02	50.6-77.0	70.7
3	MF-092717-03	127-297	242

Calibration Curve:



Review/Approval  
 Name: Sivasankari Murugan Title: QC Supervisor Signature/Date: *[Signature]* 15 Jan 2018

Statement: The above product was tested using Quanterix procedures and conforms to Quanterix's specifications. It is intended for Research Use Only. Not for use in Diagnostic Procedures.