

## Certificate of Analysis

Kit Name:	Simoa™ HIV P24 Kit
Product Number:	102215
Lot Number:	500911
Expiration:	13-Mar-2018

113 Hartwell Ave.  
Lexington MA.  
02421

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

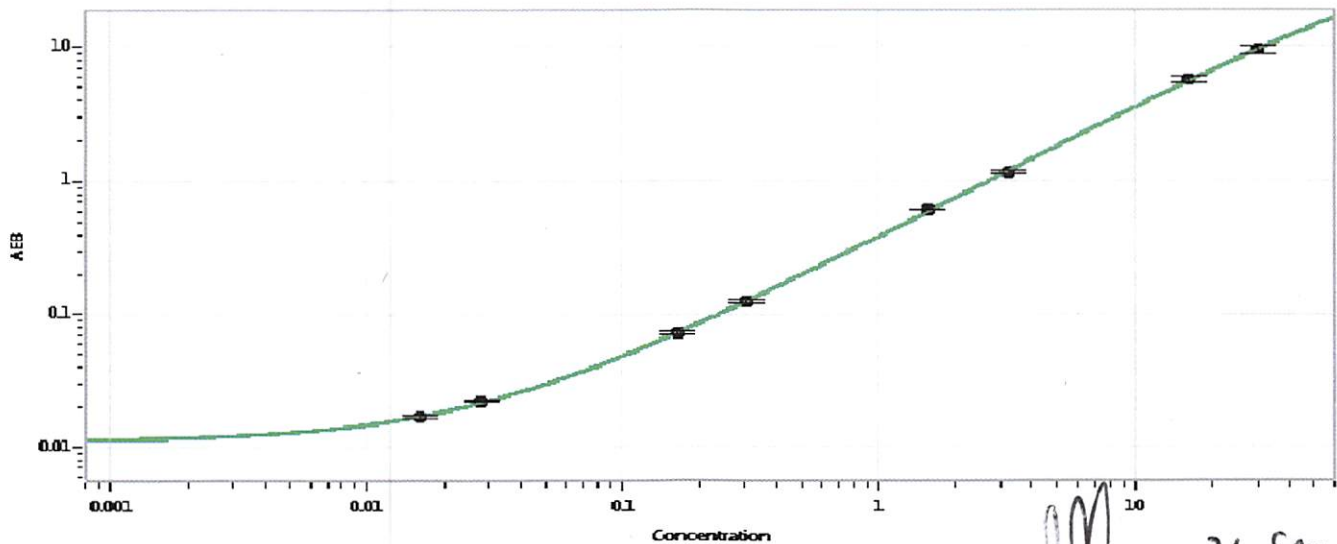
Component	Part Number	Lot Number
HIV P24 Bead Reagent	102211	72613
HIV P24 Detector Reagent	102212	72565
HIV P24 SBG Reagent	102213	72566
HIV P24 Assay Diluent	102214	72564
HIV P24 Calibrators	101310	72150
RGP Reagent	100030	N/A <sup>1</sup>
HIV P24 Controls	100702-100703	62581

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

Calibrators	Assigned Value (pg/mL)	Controls	Mean Range (pg/mL)
A	0.000	Control 1	0.070-0.189
B	0.016	Control 2	7.94-21.1
C	0.028	<b>Controls</b>	<b>Value at Release (pg/mL)</b>
D	0.164	Control 1	0.145
E	0.303	Control 2	15.5
F	1.57		
G	3.24		
H	16.2		
I	30.3		

Patient Panels	Lot Number	Mean Range (pg/mL)	Value at Release (pg/mL)
1	MF-112916-01	0.149-0.261	0.180
2	MF-081616-04	1.77-2.65	2.04
3	MF-081616-05	16.5-26.4	19.5

Calibration Curve:



James Drummey

Senior Director, Manufacturing &  
Operations

*[Handwritten Signature]*  
21-Sep-2017

Review/Approval

Name

Title

Signature/Date

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.