

Product Number:	101351
Lot Number:	501497
Expiration:	16-Mar-2019

Component	Part Number	Lot Number	Concentration
Beads	103206	MF-052418-01	3.23X10 <sup>9</sup> beads/mL
Detector/Sample Diluent	101353	823608	
Bead Diluent	100458	830608	
SBG Dilution Buffer	100376	829912	
SBG Concentrate	100439	821806	144 nM
Bead Conjugation Buffer	100453	810804	
Bead Wash Buffer	101352	MF-051418-09	
Bead Blocking Buffer	100457	823408	
Biotinylation Reaction Buffer	100460	807503	
EDC, 10mg	100022	TG264930	
EZ-Link NHS-PEG4 Biotin	100348	TC255361	
RGP	103159	N/A <sup>1</sup>	
PSA Capture Antibody	100463	81440	5.1 mg/mL
PSA Detector Antibody	100464	81441	1.07 mg/mL
Calibrator A-H	100758-100765	73180	
Controls 1 and 2	100768 & 100769	73181	

<sup>1</sup> RGP is not Kit Lot specific.


	Mean Range (pg/mL)	* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
Control 1	3.43-5.14	
Control 2	74.2-111	

Calibrator Levels (pg/mL)	
A	0.000
B	0.133
C	0.381
D	1.25
E	3.70
F	12.8
G	39.5
H	98.1

Review/Approval

Sivasankari Murugan  
Name

QC Supervisor  
Title

  
Signature/Date 15 Nov 2018

Statement: The above product was tested using Quanterix procedures and conforms to Quanterix's specifications.