

Product Number:	101351
Lot Number:	501293
Expiration:	17-Apr-2019
Platform(s):	HD-1

Component	Part Number	Lot Number	Concentration
Beads	103206	MF-052418-01	3.23X10 ⁹ beads/mL
Detector/Sample Diluent	101353	MF-052318-11	
Bead Diluent	100458	MF-052318-03	
SBG Dilution Buffer	100376	MF-052318-10	
SBG Concentrate	100439	81072	186nM
Bead Conjugation Buffer	100453	MF-051718-04	
Bead Wash Buffer	101352	MF-051418-09	
Bead Blocking Buffer	100457	MF-051418-07	
Biotinylation Reaction Buffer	100460	MF-051718-08	
EDC, 10mg	100022	TC255240	
EZ-Link NHS-PEG4 Biotin	100348	TC255361	
RGP	103159	N/A ¹	
PSA Capture Antibody	100463	81440	5.1 mg/mL
PSA Detector Antibody	100464	81441	1.07 mg/mL
Calibrator A-H	100758-100765	80230	
Controls 1 and 2	100768 & 100769	80231	

¹ 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	2.55-3.83	
Control 2	57.9-88.1	

Calibrator Levels (pg/mL)	
A	0.000
B	0.100
C	0.327
D	1.14
E	3.44
F	12.1
G	37.1
H	93.4

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

 15 Jun 2018
Signature/Date

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