

Product Number:	101351
Lot Number:	501221
Expiration:	24-Nov-2018
Platform(s):	HD-1

Component	Part Number	Lot Number	Concentration
Beads	103206	MF-041018-10	3.21X10 ⁹ beads/mL
Detector/Sample Diluent	101353	MF-112417-06	
Bead Diluent	100458	MF-112817-23	
SBG Dilution Buffer	100376	MF-041518-04	
SBG Concentrate	100439	80924	165nM
Bead Conjugation Buffer	100453	MF-031418-09	
Bead Wash Buffer	101352	MF-112417-09	
Bead Blocking Buffer	100457	MF-112417-08	
Biotinylation Reaction Buffer	100460	MF-112817-14	
EDC, 10mg	100022	SK253612	
EZ-Link NHS-PEG4 Biotin	100348	SI254541	
RGP	103159	N/A ¹	
PSA Capture Antibody	100463	72371	5.3 mg/mL
PSA Detector Antibody	100464	72133	1.08 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	2.54-3.81	
Control 2	57.7-86.6	

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

 23 APR 2018
Signature/Date

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