

Product Number:	101351
Lot Number:	501577
Expiration:	16-Mar-2019

Component	Part Number	Lot Number	Concentration
Beads	103206	823619	3.29X10 ⁹ beads/mL
Detector/Sample Diluent	101353	MF-052318-14	
Bead Diluent	100458	830608	
SBG Dilution Buffer	100376	829916	
SBG Concentrate	100439	824115	155 nM
Bead Conjugation Buffer	100453	830908	
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 ¹	
PSA Capture Antibody	100463	81440	5.1 mg/mL
PSA Detector Antibody	100464	831601	1.0 mg/mL
Calibrator A-H	100758-100765	833801	
Controls 1 and 2	100768 & 100769	833802	

¹ RGP is not Kit Lot specific.

	Mean Range (pg/mL)	
Control 1	2.28-3.42	* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
Control 2	49.8-74.6	

Calibrator Levels (pg/mL)	
A	0.000
B	0.093
C	0.278
D	0.948
E	2.95
F	10.6
G	32.1
H	75.1

Review/Approval

Patrizia Stadler

QC Associate
Director

P. Stadler 1/10/2019

Name

Title

Signature/Date

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