

Product Number:	101351
Lot Number:	501851
Expiration:	9-Jan-2020

Component	Part Number	Lot Number	Concentration
Beads	103206	823620	3.41X10 <sup>9</sup> beads/mL
Detector/Sample Diluent	101353	913111	
Bead Diluent	100458	908916	
SBG Dilution Buffer	100376	916108	
SBG Concentrate	103397	910511	50nM
Bead Conjugation Buffer	100453	909308	
Bead Wash Buffer	101352	903602	
Bead Blocking Buffer	100457	900921	
Biotinylation Reaction Buffer	100460	900919	
EDC, 10mg	100022	UB281110	
EZ-Link NHS-PEG4 Biotin	100348	UA272225	
RGP	103159	N/A <sup>1</sup>	
PSA Capture Antibody	100463	903614	5.0 mg/mL
PSA Detector Antibody	100464	831601	1.0 mg/mL
Calibrator A-H	100758-100765	911301/912101	
Controls 1 and 2	100768-100769	912102	

<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	2.29-3.60	
Control 2	54.3-81.5	

Calibrator Levels (pg/mL)	
A	0.000
B	0.086
C	0.288
D	0.980
E	3.23
F	10.6
G	34.6
H	82.9

Review/Approval

Patrizia Stadler

Assoc. Director, QC

*P. Stadler* 24 JUN 2019

Name

Title

Signature/Date

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