

Product Number:	101351
Lot Number:	501640
Expiration:	5-Nov-2019

Component	Part Number	Lot Number	Concentration
Beads	103206	823619	3.29X10 ⁹ beads/mL
Detector/Sample Diluent	101353	834714	
Bead Diluent	100458	833308	
SBG Dilution Buffer	100376	834903	
SBG Concentrate	103397	832316	50nM
Bead Conjugation Buffer	100453	830908	
Bead Wash Buffer	101352	903602	
Bead Blocking Buffer	100457	900921	
Biotinylation Reaction Buffer	100460	900919	
EDC, 10mg	100022	TI276003	
EZ-Link NHS-PEG4 Biotin	100348	TH271791	
RGP	103159	N/A ¹	
PSA Capture Antibody	100463	903614	5.0 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 07 Feb 2019

Name

Title

Signature/Date

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