

Product Number:	101351
Lot Number:	501938
Expiration:	5-Feb-2020

Component	Part Number	Lot Number	Concentration
Beads	103206	823620	3.41X10 <sup>9</sup> beads/mL
Detector/Sample Diluent	101353	913109	
Bead Diluent	100458	917020	
SBG Dilution Buffer	100376	919202	
SBG Concentrate	103397	911210	50nM
Bead Conjugation Buffer	100453	919112	
Bead Wash Buffer	101352	903602	
Bead Blocking Buffer	100457	920006	
Biotinylation Reaction Buffer	100460	917619	
EDC, 10mg	100022	RM2019020000034	
EZ-Link NHS-PEG4 Biotin	100348	RM2019020000036	
RGP	103159	N/A <sup>1</sup>	
PSA Capture Antibody	100463	903614	5.0 mg/mL
PSA Detector Antibody	100464	919203	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)	
Control 1	2.34-3.60	* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
Control 2	54.9-82.3	

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* 06 Aug 2019  
 Name Title Signature/Date