

Product Number:	101351
Lot Number:	502200
Expiration:	11-Jul-2020

Component	Part Number	Lot Number	Concentration
Beads	103206	823622	3.62X10 <sup>9</sup> beads/mL
Detector/Sample Diluent	101353	934511	
Bead Diluent	100458	928215	
SBG Dilution Buffer	100376	930220	
SBG Concentrate	103397	930316	50nM
Bead Conjugation Buffer	100453	932512	
Bead Wash Buffer	101352	919219	
Bead Blocking Buffer	100457	926611	
Biotinylation Reaction Buffer	100460	924107	
EDC, 10mg	100022	RM2019120000025	
EZ-Link NHS-PEG4 Biotin	100348	RM2019120000049	
RGP	103159	N/A <sup>1</sup>	
PSA Capture Antibody	100463	903614	5.0 mg/mL
PSA Detector Antibody	100464	935009	1.1 mg/mL
Calibrator A-H	100758-100765	921201	
Controls 1 and 2	100768-100769	921202	

<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	2.16-3.24	
Control 2	51.2-76.8	

**Calibrator Levels (pg/mL)**

A	0.000
B	0.084
C	0.252
D	0.899
E	2.82
F	9.94
G	30.0
H	76.3

Review/Approval

Patrizia Stadler Assoc. Director, QC  
Name Title

*P. Stadler* 17 Jan 2020  
Signature/Date