

Product Number:	101351
Lot Number:	502951
Expiration:	17-Dec-2021

Component	Part Number	Lot Number	Concentration
Beads	103611	108822	2.49X10 ⁹ beads/mL
Detector/Sample Diluent	101353	111312	
Bead Diluent	100458	110117	
SBG Dilution Buffer	100376	108323	
SBG Concentrate	103397	107422	50nM
Bead Conjugation Buffer	100453	100709	
Bead Wash Buffer	101352	111608	
Bead Blocking Buffer	100457	109014	
Biotinylation Reaction Buffer	100460	100824	
EDC, 10mg	100022	RM2021030000096	
EZ-Link NHS-PEG4 Biotin	100348	RM2020120000031	
RGP	103159	N/A ¹	
PSA Capture Antibody	100463	105804	5.1 mg/ml
PSA Detector Antibody	100464	108106	1.0 mg/ml
Calibrator A-H	100758-100765	027208	
Control 1	100768	027209	
Control 2	100769	027209	

¹ 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	2.25-3.38	
Control 2	47.1-70.6	

Calibrator Levels (pg/mL)

A	0.000
B	0.095
C	0.281
D	0.943
E	2.95
F	9.81
G	28.5
H	73.6

Review/Approval

Jungil Park
Name

QC Supervisor
Title

[Signature] 27 May 2021
Signature/Date