

Product Number:	101351
Lot Number:	502741
Expiration:	18-Jul-2021

Component	Part Number	Lot Number	Concentration
Beads	103611	031522	3.20X10 ⁹ beads/mL
Detector/Sample Diluent	101353	031415	
Bead Diluent	100458	035807	
SBG Dilution Buffer	100376	100805	
SBG Concentrate	103397	101116	50nM
Bead Conjugation Buffer	100453	100706	
Bead Wash Buffer	101352	028314	
Bead Blocking Buffer	100457	024110	
Biotinylation Reaction Buffer	100460	100821	
EDC, 10mg	100022	RM2020120000032	
EZ-Link NHS-PEG4 Biotin	100348	RM2020100000049	
RGP	103159	N/A ¹	
PSA Capture Antibody	100463	034006	5.0 mg/ml
PSA Detector Antibody	100464	034005	1.1 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)	
Control 1	2.25-3.38	* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
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

Junpil Park
Name

QC, Supervisor
Title

 15 Feb 2021
Signature/Date