

Product Number:	101351
Lot Number:	502454
Expiration:	15-Jan-2021

Component	Part Number	Lot Number	Concentration
Beads	103611	005517	2.82X10 ⁹ beads/mL
Detector/Sample Diluent	101353	015116	
Bead Diluent	100458	007710	
SBG Dilution Buffer	100376	012309	
SBG Concentrate	103397	012511	50nM
Bead Conjugation Buffer	100453	011707	
Bead Wash Buffer	101352	014004	
Bead Blocking Buffer	100457	014011	
Biotinylation Reaction Buffer	100460	001504	
EDC, 10mg	100022	RM2020030000046	
EZ-Link NHS-PEG4 Biotin	100348	RM2020040000009	
RGP	103159	N/A ¹	
PSA Capture Antibody	100463	006901	5.0 mg/ml
PSA Detector Antibody	100464	008307	1.1 mg/mL
Calibrator A-H	100758-100765	921201	
Controls 1 and 2	100768-100769	921202	

¹ RGP is not Kit Lot specific.

	Mean Range (pg/mL)	
Control 1	2.16-3.24	* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
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

Shivani Goel
Name

QC Manager
Title

Shivani Goel (08/1/2020)
Signature/Date