

Product Number:	101351
Lot Number:	502147
Expiration:	23-Sep-2020

Component	Part Number	Lot Number	Concentration
Beads	103206	823622	3.62X10 <sup>9</sup> beads/mL
Detector/Sample Diluent	101353	924008	
Bead Diluent	100458	921315	
SBG Dilution Buffer	100376	930219	
SBG Concentrate	103397	927507	50nM
Bead Conjugation Buffer	100453	930407	
Bead Wash Buffer	101352	928910	
Bead Blocking Buffer	100457	926611	
Biotinylation Reaction Buffer	100460	924107	
EDC, 10mg	100022	RM2019100000039	
EZ-Link NHS-PEG4 Biotin	100348	RM2019080000025	
RGP	103159	N/A <sup>1</sup>	
PSA Capture Antibody	100463	925412	5.0 mg/mL
PSA Detector Antibody	100464	931607	1.0 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)	
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

Patrizia Stadler Assoc. Director, QC

*P. Stadler* 04 Dec 2019

Name

Title

Signature/Date