

Product Number:	101351
Lot Number:	503774
Expiration:	1-Jun-2023

Component	Part Number	Lot Number	Concentration
488-dyed Singleplex Beads ²	104007	231417	1.11 x 10 ⁹ beads/mL
Detector/Sample Diluent	101353	234015	
Bead Diluent	100458	234304	
SBG Dilution Buffer	100376	231814	
SBG Concentrate	103397	222904	50nM
Bead Conjugation Buffer	100453	233906	
Bead Wash Buffer	101352	220310	
Bead Blocking Buffer	100457	227911	
Biotinylation Reaction Buffer	100460	215211	
EDC, 10mg	100022	RM2022100000009	
EZ-Link NHS-PEG4 Biotin	100348	XH359829	
RGP ¹	103159	N/A	
PSA Capture Antibody	100463	220604	5.1 mg/ml
PSA Detector Antibody	100464	225612	1.0 mg/ml
Calibrator A-H	100758-100765	213502	
Control 1	100768	213503	
Control 2	100769	213503	

¹RGP is not Kit Lot specific.

²Assay Beads provided are conjugated with 488 nm fluorescent dye.

For more information, please visit

<https://go.quanterix.com/HomebrewBeadUpdate>



	Spec. Range (pg/mL)	* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
Control 1	2.09 - 3.14	
Control 2	45.9 - 68.9	

Calibrator Levels (pg/mL)	
A	0.000
B	0.090
C	0.260
D	0.878
E	2.71
F	9.56
G	26.9
H	70

Review/Approval Sonia Khillen AC Manager Sonia Koll / 22 DEC 2022
 Name Title Signature/Date