

Product Number:	101351
Lot Number:	503855
Expiration:	26-Oct-2023

Component	Part Number	Lot Number	Concentration
488-dyed Singleplex Beads ²	104007	307212	1.32 x 10 ⁹ beads/mL
Detector/Sample Diluent	101353	308307	
Bead Diluent	100458	306809	
SBG Dilution Buffer	100376	308609	
SBG Concentrate	103397	301009	50nM
Bead Conjugation Buffer	100453	306010	
Bead Wash Buffer	101352	302410	
Bead Blocking Buffer	100457	304807	
Biotinylation Reaction Buffer	100460	306812	
EDC, 10mg	100022	RM2023010000055	
EZ-Link NHS-PEG4 Biotin	100348	RM2023020000034	
RGP ¹	103159	N/A	
PSA Capture Antibody	100463	303212	5.1 mg/ml
PSA Detector Antibody	100464	303213	1.1 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 Khillan Q.C. Manager Sonia Khillan 12/16/2023
 Name Title Signature/Date