

Product Number:	101351
Lot Number:	503373
Expiration:	13-Aug-2022

Component	Part Number	Lot Number	Concentration
488-dyed Singleplex Beads ²	104007	126705	1.53 x 10 ⁹ beads/mL
Detector/Sample Diluent	101353	131516	
Bead Diluent	100458	134609	
SBG Dilution Buffer	100376	203113	
SBG Concentrate	103397	129210	50nM
Bead Conjugation Buffer	100453	127413	
Bead Wash Buffer	101352	122511	
Bead Blocking Buffer	100457	127514	
Biotinylation Reaction Buffer	100460	203013	
EDC, 10mg	100022	RM2021100000035	
EZ-Link NHS-PEG4 Biotin	100348	RM2021120000042	
RGP ¹	103159	N/A	
PSA Capture Antibody	100463	122803	5.1 mg/ml
PSA Detector Antibody	100464	122804	1.0 mg/ml
Calibrator A-H	100758-100765	121603	
Control 1	100768	121604	
Control 2	100769	121604	

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>



	Mean Range (pg/mL)	* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
Control 1	1.69-3.70	
Control 2	42.7-64.0	

Calibrator Levels (pg/mL)	
A	0.000
B	0.043
C	0.237
D	0.937
E	2.62
F	8.76
G	26.2
H	61.6

Review/Approval

Shivani Goel
Name

QC
Title

Shivani Goel 28Feb2022
Signature/Date