

Product Number:	101351
Lot Number:	503677
Expiration:	16-Feb-2023

Component	Part Number	Lot Number	Concentration
488-dyed Singleplex Beads ²	104007	220210	1.34 x 10 ⁹ beads/mL
Detector/Sample Diluent	101353	218619	
Bead Diluent	100458	220106	
SBG Dilution Buffer	100376	222007	
SBG Concentrate	103397	209718	50nM
Bead Conjugation Buffer	100453	217505	
Bead Wash Buffer	101352	220310	
Bead Blocking Buffer	100457	220606	
Biotinylation Reaction Buffer	100460	215211	
EDC, 10mg	100022	RM2022070000003	
EZ-Link NHS-PEG4 Biotin	100348	RM2022030000015	
RGP ¹	103159	N/A	
PSA Capture Antibody	100463	220604	5.1 mg/ml
PSA Detector Antibody	100464	204107	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>



	Spec. Range (pg/mL)	* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
Control 1	1.70 - 3.55	
Control 2	43.1 - 64.6	

Calibrator Levels (pg/mL)	
A	0
B	0.043
C	0.237
D	0.937
E	2.62
F	8.76
G	26.2
H	61.6

Review/Approval: Sonia Khillan Name, QC Title, [Signature] 18 AUG 2022 Signature/Date