

Product Number:	101351
Lot Number:	503664
Expiration:	8-Mar-2023

Component	Part Number	Lot Number	Concentration
488-dyed Singleplex Beads <sup>2</sup>	104007	208312	1.28 x 10 <sup>9</sup> beads/mL
Detector/Sample Diluent	101353	217217	
Bead Diluent	100458	220106	
SBG Dilution Buffer	100376	217316	
SBG Concentrate	103397	215710	50nM
Bead Conjugation Buffer	100453	217503	
Bead Wash Buffer	101352	220310	
Bead Blocking Buffer	100457	210318	
Biotinylation Reaction Buffer	100460	215211	
EDC, 10mg	100022	RM2022070000003	
EZ-Link NHS-PEG4 Biotin	100348	RM2022030000015	
RGP <sup>1</sup>	103159	N/A	
PSA Capture Antibody	100463	204106	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	

<sup>1</sup>RGP is not Kit Lot specific.

<sup>2</sup>Assay 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

*Matt Smart*  
Name

QC  
Title

*MAT 03Aug2022*  
Signature/Date