

Product Number:	101351
Lot Number:	504484
Expiration:	22-Feb-2025

Component	Part Number	Lot Number	Control Lot Number	Concentration
488-dyed Singleplex Beads ²	104007	403913	N/A	1.14 x 10 ⁹ beads/mL
Detector / Sample Diluent	101353	412313	N/A	
Bead Diluent	100458	414211	N/A	
SBG Dilution Buffer	100376	433720	N/A	
SBG Concentrate	103397	319213	319213A	50 nM
Bead Conjugation Buffer	100453	409213	N/A	
Bead Wash Buffer	101352	415103	N/A	
Bead Blocking Buffer	100457	411728	N/A	
Biotinylation Reaction Buffer	100460	405509	N/A	
EDC, 10mg	100022	RM2024070000010	N/A	
EZ-Link NHS-PEG4 Biotin	100348	RM2024100000053	N/A	
RGP ¹	103159	N/A ¹	N/A ¹	
PSA Capture Antibody	100463	414114	N/A	5.1 mg/mL
PSA Detector Antibody	100464	406102	N/A	1.1 mg/mL
Calibrator A-H	100758 - 100765	305311	N/A	
Control 1	100768	305312	N/A	
Control 2	100769	305312	N/A	

¹ Reagents are not Kit Lot Specific

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

Release Materials*			* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
	Lot	Spec. Range (pg/mL)	
Control 1	305312	2.05 - 3.07	
Control 2	305312	41.1 - 65.1	

Calibrator Levels

A	0.000
B	0.077
C	0.246
D	0.873
E	2.61
F	8.43
G	25.8
H	64.4

Review/Approval

TAM LEE DOC CONTROL II, QA *Tam Lee* 19 DEC 2024

Name

Title

Signature / Date