

Product Number:	101351
Lot Number:	504532
Expiration:	25-Oct-2025

Component	Part Number	Lot Number	Control Lot Number	Concentration
488-dyed Singleplex Beads ²	104007	505210	N/A	1.33 x 10 ⁹ beads/mL
Detector / Sample Diluent	101353	502213	N/A	
Bead Diluent	100458	502411	502411A	
SBG Dilution Buffer	100376	506219	506219A	50 nM
SBG Concentrate	103397	506201	506201A	
Bead Conjugation Buffer	100453	505807	505807A	
Bead Wash Buffer	101352	415103	N/A	
Bead Blocking Buffer	100457	411728	N/A	
Biotinylation Reaction Buffer	100460	416316	N/A	
EDC, 10mg	100022	RM2024110000028	N/A	
EZ-Link NHS-PEG4 Biotin	100348	RM2024100000053	N/A	
RGP ¹	103159	N/A ¹	N/A ¹	

¹ Reagents are not Kit Lot Specific

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

Review/Approval

Dipesh Patel
Name

Sr. Document
Title

AT / 12 MAR 2025
Signature/Date