

Product Number:	101351
Lot Number:	536661
Expiration:	20-Jun-2026

Component	Part Number	Lot Number	Concentration
488-dyed Singleplex Beads ²	104007	513911	1.20 x 10 ⁹ beads/mL
Detector / Sample Diluent	101353	523824	
Bead Diluent	100458	530005	
SBG Dilution Buffer	100376	530913	
SBG Concentrate	103397	515406	
Bead Conjugation Buffer	100453	528907	50 nM
Bead Wash Buffer	101352	521010	
Bead Blocking Buffer	100457	521011	
Biotinylation Reaction Buffer	100460	510810	
EDC, 10mg	100022	RM2025060000020	
EZ-Link NHS-PEG4 Biotin	100348	RM2025110000001	
RGP ¹	103159	N/A ¹	

¹ Reagents are not Kit Lot Specific

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