

Product Number:	101351
Lot Number:	536624
Expiration:	20-Jun-2026

Component	Part Number	Lot Number	Concentration
488-dyed Singleplex Beads ²	104007	520513	1.29 x 10 ⁹ beads/mL
Detector / Sample Diluent	101353	523826	50 nM
Bead Diluent	100458	530005	
SBG Dilution Buffer	100376	530913	
SBG Concentrate	103397	530003	
Bead Conjugation Buffer	100453	528907	
Bead Wash Buffer	101352	521021	
Bead Blocking Buffer	100457	521020	
Biotinylation Reaction Buffer	100460	523111	
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 fluorescent dye.

Review/Approval

Dhwani Patel *QA Specialist II* *SDP/20 Nov 2025*
 Name Title Signature/Date