

Product Number:	102040
Lot Number:	502499
Expiration:	17-Mar-2021

Component	Part Number	Lot Number
Detector/Sample Diluent	101359	019711
Biotinylation Reaction Buffer	101358	015108
Bead Conjugation Buffer	101357	015107
Bead Blocking Buffer	101356	019812
Bead Wash Buffer	101355	008007
Bead Diluent Buffer	100458	007710

Review/Approval

Shivani Goel  
Name

QC Manager  
Title

*Shivani Goel* 19 Aug 2020  
Signature/Date