

Product Number:	102040
Lot Number:	502458
Expiration:	17-Dec-2020

Component	Part Number	Lot Number
Detector/Sample Diluent	101359	003515
Biotinylation Reaction Buffer	101358	015108
Bead Conjugation Buffer	101357	935108
Bead Blocking Buffer	101356	935209
Bead Wash Buffer	101355	008007
Bead Diluent Buffer	100458	001517

Review/Approval

Shivani Goel

QC Manager

Shivani Goel 08/01/2020

Name

Title

Signature/Date