

Product Number:	102040
Lot Number:	503039
Expiration:	29-Mar-2022

Component	Part Number	Lot Number
Detector/Sample Diluent	101359	117115
Biotinylation Reaction Buffer	101358	117222
Bead Conjugation Buffer	101357	108827
Bead Blocking Buffer	101356	110521
Bead Wash Buffer	101355	112103
Bead Diluent Buffer	100458	110117

Review/Approval

Shivani Goel
Name

QC
Title

Shivani Goel 12 Jul 2021
Signature/Date