

Product Number:	102040
Lot Number:	502918
Expiration:	8-Jan-2022

Component	Part Number	Lot Number
Detector/Sample Diluent	101359	108821
Biotinylation Reaction Buffer	101358	100825
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 11 May 2021
Signature/Date