

Product Number:	102040	
Lot Number:	503210	
Expiration:	15-Apr-2022	

Component	Part Number	Lot Number
Detector/Sample Diluent	101359	126809
Biotinylation Reaction Buffer	101358	120909
Bead Conjugation Buffer	101357	127410
Bead Blocking Buffer	101356	110523
Bead Wash Buffer	101355	122510
Bead Diluent Buffer	100458	118710

Review/Approval

Shivani Goel
Name

QC
Title

Shivani Goel 08 NOV 2021
Signature/Date