

Product Number:	102040
Lot Number:	502785
Expiration:	28-Aug-2021

Component	Part Number	Lot Number
Detector/Sample Diluent	101359	102708
Biotinylation Reaction Buffer	101358	100818
Bead Conjugation Buffer	101357	100703
Bead Blocking Buffer	101356	024111
Bead Wash Buffer	101355	031127
Bead Diluent Buffer	100458	101606

Review/Approval

Tumpil Park

Name

QC Supervisor

Title

[Signature] 09 Mar 2021

Signature/Date