

Product Number:	102040
Lot Number:	503532
Expiration:	12-Dec-2022

Component	Item Number	Lot Number
Detector/Sample Diluent	101359	217216
Biotinylation Reaction Buffer	101358	214005
Bead Conjugation Buffer	101357	216704
Bead Blocking Buffer	101356	206118
Bead Wash Buffer	101355	206115
Bead Diluent Buffer	100458	134609

Review/Approval

Matt Smart
Name

QC

Title

MA 29 Jun 2022
Signature/Date