

Product Number:	102040
Lot Number:	503665
Expiration:	2-Mar-2023

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	220106

Review/Approval

Matt Sinsat
Name

QC
Title

[Signature] 03Apr2022
Signature/Date