

Product Number:	102040
Lot Number:	503487
Expiration:	2-Mar-2023

Component	Item Number	Lot Number
Detector/Sample Diluent	101359	212705
Biotinylation Reaction Buffer	101358	209003
Bead Conjugation Buffer	101357	210809
Bead Blocking Buffer	101356	206118
Bead Wash Buffer	101355	206115
Bead Diluent Buffer	100458	210310

Review/Approval

Matt Smart
Name

QC
Title

MSA 19 May 2022
Signature/Date