

Product Number:	102040
Lot Number:	502181
Expiration:	1-Aug-2020

Component	Part Number	Lot Number
Detector/Sample Diluent	101359	932502
Biotinylation Reaction Buffer	101358	924108
Bead Conjugation Buffer	101357	926906
Bead Blocking Buffer	101356	935209
Bead Wash Buffer	101355	932505
Bead Diluent Buffer	100458	921319

Review/Approval

Patrizia Stadler

Assoc. Director, QC

*P. Stadler* 30 Dec 2019

Name

Title

Signature/Date