

Product Number:	102040
Lot Number:	502170
Expiration:	11-Jul-2020

Component	Part Number	Lot Number
Detector/Sample Diluent	101359	928214
Biotinylation Reaction Buffer	101358	924104
Bead Conjugation Buffer	101357	920505
Bead Blocking Buffer	101356	935209
Bead Wash Buffer	101355	919213
Bead Diluent Buffer	100458	921319

Review/Approval

Patrizia Stadler

Assoc. Director, QC

p.stadler *Oct 14, 2020*

Name

Title

Signature/Date