

Product Number:	102040
Lot Number:	501856
Expiration:	11-Feb-2020

Component	Part Number	Lot Number
Detector/Sample Diluent	101359	913110
Bead Conjugation Buffer	101357	909319
Bead Diluent Buffer	100458	904223
Biotinylation Reaction Buffer	101358	904308
Bead Blocking Buffer	101356	906502
Bead Wash Buffer	101355	910013

Review/Approval

Patrizia Stadler

Assoc. Director, QC

P. Stadler 18 JUN 2019

Name

Title

Signature/Date