

Product Number:	102040
Lot Number:	501926
Expiration:	12-Feb-2020

Component	Part Number	Lot Number
Detector/Sample Diluent	101359	917613
Bead Conjugation Buffer	101357	919109
Bead Diluent Buffer	100458	917020
Biotinylation Reaction Buffer	101358	904309
Bead Blocking Buffer	101356	906514
Bead Wash Buffer	101355	919213

Review/Approval

Patrizia Stadler

Assoc. Director, QC

P. Stadler 26 Jul 2019

Name

Title

Signature/Date