

Product Number:	102040
Lot Number:	502107
Expiration:	6-Mar-2020

Component	Part Number	Lot Number
Detector/Sample Diluent	101359	926016
Biotinylation Reaction Buffer	101358	917620
Bead Conjugation Buffer	101357	916503
Bead Blocking Buffer	101356	906502
Bead Wash Buffer	101355	919213
Bead Diluent Buffer	100458	917022

Review/Approval

Patrizia Stadler

Assoc. Director, QC

P. Stadler 08 NOV 2019

Name

Title

Signature/Date