

Product Number:	101354
Lot Number:	502053
Expiration:	6-Mar-2020

Component	Item Number	Lot Number
Beads	103207	912113
Detector/Sample Diluent	101359	924905
Biotinylation Reaction Buffer	101358	924104
Bead Conjugation Buffer	101357	920505
Bead Blocking Buffer	101356	906502
Bead Wash Buffer	101355	919221
Bead Diluent Buffer	100458	921315

Bead Count (beads/mL)	3.46x10 ⁹
--------------------------	----------------------

Review/Approval

Patrizia Stadler QC Assoc. Director
Name Title

P. Stadler 10 OCT 2019
Signature/Date

Statement: The above product was tested using Quanterix procedures and conforms to Quanterix's specifications.