

Product Number:	102040
Lot Number:	503433
Expiration:	28-Nov-2022

Component	Item Number	Lot Number
Detector/Sample Diluent	101359	134925
Biotinylation Reaction Buffer	101358	203014
Bead Conjugation Buffer	101357	134611
Bead Blocking Buffer	101356	206118
Bead Wash Buffer	101355	133220
Bead Diluent Buffer	100458	134609

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 06Apr2022
Signature/Date

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