

Product Number:	102907
Lot Number:	502481
Expiration:	22-Jul-2021
Storage:	2 - 8°C

Component	Part Number	Lot Number	Means of Analysis	Pass / Fail
Lysate Diluent A Reagent	102908	020415	pH meter	Pass
Lysate Diluent B Reagent	102910	020404		Pass
Lysate Diluent C Reagent	102912	020407		Pass
Lysate Diluent D Reagent	102914	020507		Pass

Review/Approval

Shivani Goel QC Manager
Name Title

Shivani Goel 28 Jul 2020
Signature/Date

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