

Product Number:	102907
Lot Number:	503391
Expiration:	26-Feb-2023
Storage:	2 - 8°C

Component	Part Number	Lot Number	Means of Analysis	Pass / Fail
Lysate Diluent A Reagent	102908	205702	pH meter	Pass
Lysate Diluent B Reagent	102910	205705		Pass
Lysate Diluent C Reagent	102912	205708		Pass
Lysate Diluent D Reagent	102914	205711		Pass

Review/Approval

Shivani Goel QC Manager
Name Title

Shivani Goel 08 Mar 2022
Signature/Date

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