

Product Number:	102907
Lot Number:	503182
Expiration:	1-Oct-2022
Storage:	2 - 8°C

Component	Part Number	Lot Number	Means of Analysis	Pass / Fail
Lysate Diluent A Reagent	102908	127407	pH meter	Pass
Lysate Diluent B Reagent	102910	127408		Pass
Lysate Diluent C Reagent	102912	127501		Pass
Lysate Diluent D Reagent	102914	127502		Pass

Review/Approval

Junil Park
Name

QC Supervisor
Title

[Signature] 13 Oct 2021
Signature/Date

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