

Product Number:	102907
Lot Number:	502724
Expiration:	25-Jan-2022
Storage:	2 - 8°C

Component	Part Number	Lot Number	Means of Analysis	Pass / Fail
Lysate Diluent A Reagent	102908	102507	pH meter	Pass
Lysate Diluent B Reagent	102910	102508		Pass
Lysate Diluent C Reagent	102912	102509		Pass
Lysate Diluent D Reagent	102914	102510		Pass

Review/Approval

Shivani Goel QC Manager
Name Title

Shivani Goel 08 Feb 2021
Signature/Date

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