

Product Number:	102907
Lot Number:	503464
Expiration:	17-Mar-2023
Storage:	2 - 8°C

Component	Part Number	Lot Number	Means of Analysis	Pass / Fail
Lysate Diluent A Reagent	102908	210904	pH meter	Pass
Lysate Diluent B Reagent	102910	211307		Pass
Lysate Diluent C Reagent	102912	207609		Pass
Lysate Diluent D Reagent	102914	210907		Pass


Review/Approval

Richa Shrestha

QC Supervisor

Name

Title

 28Apr2022
Signature/Date

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