

Product Number:	102907
Lot Number:	503748
Expiration:	19-Apr-2023
Storage:	2 - 8°C

Component	Part Number	Lot Number	Means of Analysis	Pass / Fail
Lysate Diluent A Reagent	102908	223408	pH meter	Pass
Lysate Diluent B Reagent	102910	218102		Pass
Lysate Diluent C Reagent	102912	210924		Pass
Lysate Diluent D Reagent	102914	232003		Pass

Review/Approval

Sonia Khillan QC Manager Sonia Khillan 01/DEC/2022

Name	Title	Signature/Date
------	-------	----------------