

Product Number:	102907
Lot Number:	503780
Expiration:	22-Aug-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	235005		Pass
Lysate Diluent C Reagent	102912	235008		Pass
Lysate Diluent D Reagent	102914	232003		Pass

Review/Approval

Matt Koularis

QC Supervisor

[Signature] 09Jan2023

Name

Title

Signature/Date