

Product Number:	102907
Lot Number:	503508
Expiration:	17-Mar-2023
Storage:	2 - 8°C

Component	Part Number	Lot Number	Means of Analysis	Pass / Fail
Lysate Diluent A Reagent	102908	210921	pH meter	Pass
Lysate Diluent B Reagent	102910	214703		Pass
Lysate Diluent C Reagent	102912	207609		Pass
Lysate Diluent D Reagent	102914	210922		Pass

Review/Approval

Matt Smart
Name

QC

Title

MAT 03 Jun 2022
Signature/Date