

Product Number:	102907
Lot Number:	503662
Expiration:	17-Mar-2023
Storage:	2 - 8°C

Component	Part Number	Lot Number	Means of Analysis	Pass / Fail
Lysate Diluent A Reagent	102908	218909	pH meter	Pass
Lysate Diluent B Reagent	102910	218912		Pass
Lysate Diluent C Reagent	102912	207612		Pass
Lysate Diluent D Reagent	102914	219211		Pass

Review/Approval

Sonia Khillan

QC

Sonia Khillan / 27 JUL 2022

Name

Title

Signature/Date