

Product Number:	101735
Lot Number:	504199
Expiration Date:	12-Aug-25

Component	Part Number	Lot Number	Control Number
SBG Dilution Buffer	100376	404318	N/A
SBG Concentrate	103397	319211	319211A

Laboratory Analysis

Characteristic	Means of Analysis	PASS / FAIL
Functionality	HD(X/1) Analyzer	PASS

Result

SBG Concentrate Concentration (nM)	50
------------------------------------	----

Review/Approval

ROBERT SOMOGYI

Name

QC MANAGER

Title

[Handwritten Signature] 29 MAR 2024

Signature/Date