

Product Number:	101735
Lot Number:	504117
Expiration Date:	12-Dec-2024

Component	Part Number	Lot Number
SBG Dilution Buffer	100376	401617
SBG Concentrate	103397	319210A

Laboratory Analysis

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

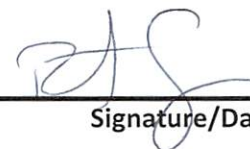
Result

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

Review/Approval

ROBERT SOMOGYI

QC MANAGER



25 JAN 2024

Name

Title

Signature/Date