

Product Number:	101735
Lot Number:	504398
Expiration Date:	3-Mar-2026

Component	Part Number	Lot Number	Control Lot Number
SBG Dilution Buffer	100376	424714	N/A
SBG Concentrate	103397	319213	319213A

Laboratory Analysis

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

Result

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

Review/Approval

Dilisha Patel
Name

Doc control II QA
Title

ENT / 18 SEP 2024
Signature/Date