

<b>Product Number:</b>	101735
<b>Lot Number:</b>	504502
<b>Expiration Date:</b>	2-Jun-2026

Component	Part Number	Lot Number	Control Lot Number
SBG Dilution Buffer	100376	433720	N/A
SBG Concentrate	103397	502104	502104A

**Laboratory Analysis**

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

**Result**

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

Review/Approval

*Dhruvi Patel*

Name

*DOC Control I, QA*

Title

*DP 30 Jan 2025*

Signature/Date