

Product Number:	101735
Lot Number:	504365
Expiration Date:	20-Dec-2025

Component	Part Number	Lot Number
SBG Dilution Buffer	100376	417210
SBG Concentrate	103397	420404

Laboratory Analysis

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

Result

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

Review/Approval

TAM LEE

DOC CONTROL II - QA

Tam Lee 03SEP2024

Name

Title

Signature/Date