

Product Number:	103397
Lot Number:	330603
Expiration:	2-Nov-2024
Storage:	2-8°C

Laboratory Analysis

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

RESULT

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

Review/Approval

*Sonia Khillan*  
Name

*Senior QC Analyst*  
Title

*Sonia Khillan 14 NOV 2023*  
Signature/Date