

Product Number:	103397
Lot Number:	204912
Expiration:	8-Feb-2023
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
HD-1 Analyzer	PASS

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

Review/Approval

Richa Shrestha
Name

QC Supervisor
Title

 21A82021
Signature/Date