

Product Number:	103397
Lot Number:	129210
Expiration:	19-Oct-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
HD-1 Analyzer	PASS

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

Review/Approval Shivani Goel QC Shivani Goel 15 Feb 2022
Name Title Signature/Date