

Product Number:	103397
Lot Number:	125015
Expiration:	7-Sep-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
HD-1 Analyzer	PASS

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

Review/Approval

Shivam Goel
Name

QC
Title

Shivam Goel 01 Nov 2021
Signature/Date