

Product Number:	103397
Lot Number:	120035
Expiration:	19-Jul-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
HD-1 Analyzer	PASS

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

Review/Approval

Shivani Goel

Name

QC

Title

Shivani Goel 23 Sep 2021

Signature/Date