

Product Number:	103397
Lot Number:	209717
Expiration:	22-Feb-2023
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
HD-1 Analyzer	PASS

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

Review/Approval

Matt Smart
Name

QC
Title

[Signature] 26/01/2022
Signature/Date