

Product Number:	103397
Lot Number:	215712
Expiration:	6-Jun-2023
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
HD-1 Analyzer	PASS

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

Review/Approval

Sonia Khillan

Name

QC

Title

Sonia Khillan / 17 AUG 2022

Signature/Date