

Product Number:	103397
Lot Number:	420404
Expiration:	22-Jan-2026
Storage:	2-8°C

**Laboratory Analysis**

Characteristic	Means of Analysis	PASS / FAIL
Functionality	HD(X/1) Analyzer	PASS

**RESULT**

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

Review/Approval

Robert Somogyi

QC Manager

 24 JUL 2024

Name

Title

Signature/Date