

Product Number:	103397
Lot Number:	515406
Expiration:	3-Dec-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

Kathleen Cornell

Name

*Quality Control
Supervisor*

Title

[Signature] 10 JUN 2025

Signature/Date