

Product Number:	103650
Lot Number:	205307
Expiration:	22-May-2022
Storage:	2-8°C



Data below represents results generated on the Simoa® HD-1 Analyzer.

Laboratory Analysis

Characteristic	Result
HD-1 Analyzer	PASS

SBG Concentrate Concentration (nM)	138
------------------------------------	-----

Review/Approval

Robert Sonogyi
Shivani Goel

QC Manager

Name

Title

Signature/Date