

Product Number:	103650
Lot Number:	207404
Expiration:	15-Jun-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)	129.8
------------------------------------	-------

Review/Approval	Shivani Goel	QC Manager	<i>Shivani Goel</i> 05Apr2022
	Name	Title	Signature/Date