

Product Number:	104074
Lot Number:	504226
Paired Reagent Kit: (104073)	504139
Expiration:	22-Jul-2025
Platform(s):	HD-X

Component	Part Number	Lot Number
NF-Light v2 Adv Sample Diluent Reagent	104071	407604
NF-Light v2 Adv Calibrator Dil Reagent	104072	402219
NF-light Calibrator Concentrate	102619	329214



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

Calibrator Concentrate Value Assignment (pg/mL)	9,983
--	-------

Review/Approval

Cody Carlson

QC Manager

Cody Carlson 24 MAY 2024

Name

Title

Signature/Date