

Product Number:	104112
Lot Number:	503681
Paired Reagent Kit:	503680
Expiration:	13-Sep-2023
Platform(s):	HD-1 / HD-X & SR-X

Component	Part Number	Lot Number
Sample Diluent	104101	225605
Calibrator Diluent	104103	226215
Calibrator Concentrate	104102	226601

Calibrator Concentrate Value Assignment (pg/mL)	38819
---	-------

Review/Approval

Tony Lam

Name

QC

Title

[Signature]

Signature/Date

23 Nov 2022