

Product Number:	103807
Lot Number:	502911
Paired Reagent Kit	502864 (Product No. 103806)
Expiration:	11-Oct-2021
Platform(s):	HD-1 / HD-X & SR-X

Component	Part Number	Lot Number
Sample Diluent	103800	110114
Calibrator Diluent	103802	110116
Calibrator Concentrate	103801	105602

Calibrator Concentrate Value Assignment (pg/mL)	3840
--	------

Review/Approval

Junpil Park
Name

QC Supervisor
Title

 06 May 2021
Signature/Date