

Product Number:	103770
Lot Number:	503002
Paired Reagent Kit	502935 (Product No. 103769)
Expiration:	15-Mar-2022
Platform(s):	HD-X

Component	Part Number	Lot Number
Sample Diluent	103763	112807
Calibrator Diluent	103765	107424
Calibrator Concentrate	103764	113406

Calibrator Concentrate Value Assignment (ng/mL)	5086
--	------

Review/Approval

Junpil Park
Name

QC Supervisor
Title

 24 Jun 2021
Signature/Date