

CERTIFICATE OF ANALYSIS

10X Assay Buffer			
Catalog#	Size	Lot#	Expiry
529	Bulk	19N29	06/2021
686	15 mL	19L39	06/2021
6161	30 mL	20N40	06/2021
685	60 mL	19T3	06/2021
6259	125 mL	NA	NA

Specifications	Criteria	Lot Data	Results
pH at 10X	6.6-7.0	6.88	Pass
pH at 1X	7.2-7.6	7.45	Pass
Appearance	Clear liquid	Conforms	Pass
Conductivity μ MHOs at 10X	100,000-120,000	117,400	Pass
Storage	Refrigerated or room temperature	Room temperature	Pass
Documentation	Approved	Approved	Pass

Quality Control Tracy Murphy
 Tracy Murphy

Date 6/24/2020
 6/24/2020