

CERTIFICATE OF ANALYSIS

10X Assay Buffer

Catalog#	Size	Lot#	Expiry
529	Bulk	20R30	03/2022
686	15 mL	NA	NA
6161	30 mL	NA	NA
685	60 mL	20R31	03/2022
6259	125 mL	NA	NA

Specifications	Criteria	Lot Data	Results
pH at 10X	6.6-7.0	6.91	Pass
pH at 1X	7.2-7.6	7.52	Pass
Appearance	Clear liquid	Conforms	Pass
Conductivity μ MHOs at 10X	100,000-120,000	118,800	Pass
Storage	Refrigerated or room temperature	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control

Sally Hed Dahlquist
 Sally Hed Dahlquist

Date

3/20/2020
 3/20/2020