

CERTIFICATE OF ANALYSIS

10X PBS (PB1)

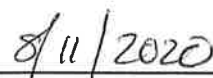
Catalog#	Size	Lot#	Expiry
529	Bulk	20A6	04/2022
6252	25 mL	NA	NA
6157	100 mL	20G6	04/2022
6158	500 mL	NA	NA
6159	1 L	NA	NA
6160	10 L	NA	NA

Specifications	Criteria	Lot Data	Results
pH at 10X	6.6-7.0	6.86	Pass
pH at 1X	7.2-7.6	7.44	Pass
Appearance	Clear liquid	Conforms	Pass
Conductivity μ MHOs at 10X	100,000-120,000	118,500	Pass
Storage	Refrigerated or room temperature	Room temperature	Pass
Documentation	Approved	Approved	Pass

Quality Control


Sally Hed Dahlquist

Date



8/11/2020