

CERTIFICATE OF ANALYSIS

Phosph-Free Block			
Catalog#	Size	Lot#	Expiry
5141	Bulk	17A37	04/2019
6262	100 mL	17A38 18S19	04/2019 04/2019
6263	500 mL	18S20	04/2019
6264	1 L	17A39	04/2019
6265	10 L	NA	NA

Specifications	Criteria	Lot Data	Results
pH	7.1-7.5	7.30	Pass
Appearance	Clear liquid	Conforms	Pass
Conductivity μ MHOs	Report measurement	14,010	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control 
 Kristi Strandberg

Date 09/10/18
 09/10/2018