

# CERTIFICATE OF ANALYSIS

<b>Phosph-Free Block</b>			
<b>Catalog#</b>	<b>Size</b>	<b>Lot#</b>	<b>Expiry</b>
5141	Bulk	17A37	04/2019
6262	100 mL	17A38 18S19	04/2019 04/2019
6263	500 mL	18S20 18T1	04/2019 04/2019
6264	1 L	17A39	04/2019
6265	10 L	NA	NA

<b>Specifications</b>	<b>Criteria</b>	<b>Lot Data</b>	<b>Results</b>
pH	7.1-7.5	7.30	Pass
Appearance	Clear liquid	Conforms	Pass
Conductivity $\mu$ MHOs	Report measurement	14,010	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control   
 Kristi Strandberg

Date 10/01/18  
 10/1/2018