

CERTIFICATE OF ANALYSIS

Phosph-Free Block			
Catalog#	Size	Lot#	Expiry
5141	Bulk	19N40	06/2021
6262	100 mL	19N41	06/2021
6263	500 mL	19G33	06/2021
6264	1 L	19N42	06/2021
6265	10 L	NA	NA

Specifications	Criteria	Lot Data	Results
pH	7.1-7.5	7.32	Pass
Appearance	Clear liquid	Conforms	Pass
Conductivity μ MHOs	Report measurement	14,270	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control Tracy Murphy
 Tracy Murphy

Date 8/13/19
 8/13/2019