

CERTIFICATE OF ANALYSIS

Monster Block™			
Catalog#	Size	Lot#	Expiry at 2-8°C
5149	Bulk	17R24	03/2019
6295	100 mL	NA	NA
6297	1 L	17R25	03/2019
6298	10 L	NA	NA

Specifications	Criteria	Lot Data	Results
pH	7.1-7.6	7.46	Pass
Appearance	Clear liquid	Conforms	Pass
Conductivity μ MHOs	13,000-20,000	16,190	Pass
A ₂₈₀ of 1:2 dilution	0.35-0.75	0.507	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

This is an assurance that the specified lot conforms to all requirements of the product specifications.
 This information is believed to be accurate and is offered in good faith.
 Product is for research use only. Not for use in diagnostic procedures.

Quality Assurance: 
 Kristi Strandberg

Date: 03/16/17
 03/16/2017