

CERTIFICATE OF ANALYSIS

Monster Block™			
Catalog#	Size	Lot#	Expiry at 2-8°C
5149	Bulk	19F20	02/2021
6295	100 mL	19F21 19D21	02/2021 02/2021
6296	500 mL	20S28	02/2021
6297	1 L	19F22	02/2021
6298	10 L	NA	NA

Specifications	Criteria	Lot Data	Results
pH	7.1-7.6	7.37	Pass
Appearance	Clear liquid	Conforms	Pass
Conductivity μ MHOs	13,000-20,000	16,290	Pass
A ₂₈₀ of 1:2 dilution	0.35-0.75	0.549	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: Sally Hed Dahlquist
 Sally Hed Dahlquist

Date: 9/21/2020
 9/21/2020