



CERTIFICATE OF ANALYSIS

HRP Conjugate Stock Stabilizer, 5X			
Catalog#	Size	Lot#	Expiry at 2-8°C
52	Bulk	20R39	03/2022
6173	25 mL	20D2	03/2022
667	100 mL	20R40 20Y5	03/2022 03/2022
668	500 mL	20R41	03/2022
669	1 L	20R42 20D13	03/2022 03/2022
670	10 L	NA	NA

Specifications	Criteria	Lot Data	Results
pH at 5X	7.2-7.6	7.39	Pass
Appearance	Clear to light yellow liquid	Conforms	Pass
Conductivity μMHOs	55,000-70,000	63,000	Pass
A₂₈₀ of 1:100 dilution	0.5-0.85	0.691	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.

Quality Assurance Tracy Murphy
 Tracy Murphy
 Date 12/10/2020
 12/10/2020