

# CERTIFICATE OF ANALYSIS

| <b>5X HRP Conjugate Stock Stabilizer Diluent (5X CS6)</b> |             |                |                        |
|-----------------------------------------------------------|-------------|----------------|------------------------|
| <b>Catalog#</b>                                           | <b>Size</b> | <b>Lot#</b>    | <b>Expiry at 2-8°C</b> |
| 52                                                        | Bulk        | 18B46          | 11/2020                |
| 6173                                                      | 25 mL       | 19Y18          | 11/2020                |
| 667                                                       | 100 mL      | 18B47          | 11/2020                |
| 668                                                       | 500 mL      | 18B48          | 11/2020                |
| 669                                                       | 1 L         | 18B49<br>19D19 | 11/2020<br>11/2020     |
| 670                                                       | 10 L        | NA             | NA                     |

| <b>Specifications</b> | <b>Criteria</b>                                   | <b>Lot Data</b> | <b>Results</b> |
|-----------------------|---------------------------------------------------|-----------------|----------------|
| pH at 5X              | 7.2-7.6                                           | 7.41            | Pass           |
| Appearance            | Clear to slightly cloudy,<br>yellow tinted liquid | Conforms        | Pass           |
| Conductivity uMH0s    | 60,000-80,000 uMHO                                | 63,400          | Pass           |
| Storage               | Refrigerated                                      | Refrigerated    | Pass           |
| Documentation         | Approved                                          | Approved        | Pass           |

Quality Control

*Tracy Murphy*  
 Tracy Murphy

Date

*12/6/19*  
 12/06/2019