

# CERTIFICATE OF ANALYSIS

| <b>10X ELISA Wash Buffer (WB1)</b> |             |             |                                   |
|------------------------------------|-------------|-------------|-----------------------------------|
| <b>Catalog#</b>                    | <b>Size</b> | <b>Lot#</b> | <b>Expiry at room temperature</b> |
| 527                                | Bulk        | 17G25       | 08/2019                           |
| 6251                               | 25 mL       | NA          | NA                                |
| 650                                | 100 mL      | NA          | NA                                |
| 651                                | 500 mL      | NA          | NA                                |
| 652                                | 1 L         | NA          | NA                                |
| 676                                | 10 L        | 17G26       | 08/2019                           |

| <b>Specifications</b> | <b>Criteria</b>                  | <b>Lot Data</b>  | <b>Results</b> |
|-----------------------|----------------------------------|------------------|----------------|
| pH at 10X             | 6.8 – 7.0                        | 6.84             | Pass           |
| pH at 1X              | 7.2 – 7.6                        | 7.41             | Pass           |
| Appearance            | Clear liquid                     | Conforms         | Pass           |
| Conductivity uMHOs    | 99,000 - 121,000 uMHO            | 115,733          | Pass           |
| Storage               | Room temperature or refrigerated | Room temperature | Pass           |

Quality Control

*Tracy Murphy*

Tracy Murphy

Date

*08/21/2017*

08/21/2017