



**CERTIFICATE OF ANALYSIS**

COA - 07873

**Product Description:** Diluent 100

**Catalog #/Size:** R50AA - Series

**Lot Number:** Y0270247

**Expiration Date:** 30 JUN 2019

**Statement:** This product was manufactured and tested according to approved MSD documents. This lot meets MSD's specifications.

pH: 7.4 ± 0.1

**The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.**

|                        | <b>Name</b>         | <b>Function</b>   | <b>Signature</b>           | <b>Date</b> |
|------------------------|---------------------|-------------------|----------------------------|-------------|
| <b>Preparer</b>        | Xiomara Talavera    | Quality Control   | <i>X Talavera</i>          | 08 FEB 2017 |
| <b>Review/Approval</b> | Vanessa Castleberry | Quality Assurance | <i>Vanessa Castleberry</i> | 10 FEB 2017 |