



CERTIFICATE OF ANALYSIS

COA - 07686

Product Description: Diluent 100

Catalog #/Size: R50AA-4 / 50 mL; R50AA-2 / 200 mL; R50AA-3 / 1000 mL

Lot Number: Y0270246

Expiration Date: 31 MAR 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.

| | Name | Function | Signature | Date |
|------------------------|------------------|-------------------|---------------------|-------------|
| Preparer | Xiomara Talavera | Quality Control | <i>X Talavera</i> | 04 JAN 2017 |
| Review/Approval | Dawn Stollar | Quality Assurance | <i>Dawn Stollar</i> | 04 JAN 2017 |