



CERTIFICATE OF CONFORMANCE

Page 1 of 1

Product Description: Diluent 2

Catalog #/Size: R51BB - Series

Lot Number: M0020093

Expiration Date: 30 JUN 2019

Storage Condition: $\leq -10^{\circ}\text{C}$

Statement: This product was manufactured according to approved MSD documents.

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

| | Name | Function | Signature | Date |
|-----------------|--------------|-------------------|---------------------|-------------|
| Review/Approval | Dawn Stollar | Quality Assurance | <i>Dawn Stollar</i> | 07 NOV 2017 |