



**CERTIFICATE OF CONFORMANCE**

Page 1 of 1

**Product Description:** Diluent 2

**Catalog #/Size:** R51BB - Series

**Lot Number:** M0020094

**Expiration Date:** 30 SEP 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.**

|                        | <b>Name</b>  | <b>Function</b>   | <b>Signature</b>    | <b>Date</b> |
|------------------------|--------------|-------------------|---------------------|-------------|
| <b>Review/Approval</b> | Dawn Stollar | Quality Assurance | <i>Dawn Stollar</i> | 06 JUL 2017 |