



**CERTIFICATE OF CONFORMANCE**

Page 1 of 1

**Product Description:** Diluent 2

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

**Lot Number:** M0020098

**Expiration Date:** 28 FEB 2020

**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>	19 APR 2018