



**CERTIFICATE OF CONFORMANCE**

Page 1 of 1

**Product Description:** Diluent 100

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


**Lot Number:** Y0270260

**Expiration Date:** 30 JUN 2021

**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>Review/Approval</b>	Dawn Stollar	Quality Assurance		14 MAR 2019