



**CERTIFICATE OF CONFORMANCE**

COC - 10294

**Product Description:** MSD<sup>®</sup> Diluent 2

**Catalog #/Size:** R51BB-2/200 mL; R51BB-3/40mL

**Lot Number:** M0020029

**Expiration Date:** 31 JAN 2012

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

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

**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>Preparer</b>	Meredith Yeager	Manufacturing Associate	<i>Meredith Yeager</i>	10/18/2010
<b>Review/Approval</b>	Staci Leffel	Manufacturing Manager	<i>Staci Leffel</i>	10/18/2010
<b>Review/Approval</b>	Richard Mann	Quality Assurance	<i>Richard Mann</i>	10/18/2010