



**CERTIFICATE OF CONFORMANCE**

COC - 10230

**Product Description:** MSD® Diluent 3

**Catalog #/Size:** R51BA-4/5 mL; R51BA-5 25 mL

**Lot Number:** M0030030

**Expiration Date:** 30 JUN 2012

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

Storage conditions: ≤ -10°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>	Janice Kim	Manufacturing Associate	<i>Janice Kim</i>	03/10/2010
<b>Review/Approval</b>	<i>Stacy Leffel</i>	Kit Manufacturing Mgr	<i>Stacy Leffel</i>	03/09/2010
<b>Review/Approval</b>	<i>Richard Mann</i>	Quality Assurance	<i>Richard Mann</i>	03/09/2010