



**CERTIFICATE OF CONFORMANCE**

COC - 10311

**Product Description:** MSD® Diluent 2  
**Catalog #/Size:** R51BB-4/8 mL; R51BB-3/40 mL; R51BB-2/200 mL  
**Lot Number:** M0020046  
**Expiration Date:** 31 JAN 2013  
**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>	Meredith Yeager	Manufacturing Associate		12/09/2010
<b>Review/Approval</b>	Staci Leffel	Manufacturing Manager		12/09/2010
<b>Review/Approval</b>	Richard Mann	Quality Assurance		12/09/2010