



CERTIFICATE OF CONFORMANCE

COC - 10504

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

**Catalog #/Size:** R51BB-3/40 mL; R51BB-4/8 mL

**Lot Number:** M0020079

**Expiration Date:** 30 NOV 2015

**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	Inventory Technician		15 Nov 2013
<b>Review/Approval</b>	Staci Leffel	Manufacturing Manager		15 NOV 2013