



CERTIFICATE OF CONFORMANCE

COC - 10291

**Product Description:** MSD<sup>®</sup> Diluent 3

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

**Lot Number:** M0030034

**Expiration Date:** 31 OCT 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.**

	Name	Function	Signature	Date
Preparer	Meredith Yeager	Manufacturing Associate	<i>Meredith Yeager</i>	08/30/2010
Review/Approval	<i>Staci Leffel</i>	Kit Manufacturing Manager	<i>Staci Leffel</i>	08/25/2010
Review/Approval	<i>Richard Mann</i>	QA	<i>Richard Mann</i>	08/27/2010