



CERTIFICATE OF CONFORMANCE

COC - 10248

**Product Description:** MSD<sup>®</sup> Diluent 3  
**Catalog #/Size:** R51BA-5/25 mL; R51BA-6/125 mL  
**Lot Number:** M0030032  
**Expiration Date:** 31 AUG 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	Janice Kim	Manufacturing Associate	<i>Jan Kim</i>	04/23/2010
Review/Approval	<i>Staci Leffel</i>	Kit Mngg Manager	<i>Staci Leffel</i>	04/23/2010
Review/Approval	<i>Richard Mann</i>	Quality Assurance	<i>Richard Mann</i>	04/23/2010