



**CERTIFICATE OF CONFORMANCE**

COC - 10352

**Product Description:** MSD<sup>®</sup> Human Serum Cytokine Assay Diluent

**Catalog #/Size:** R51BB-2/200 mL; R51BB-3/40mL

**Lot Number:** M0020026

**Expiration Date:** 30 NOV 2011

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

Storage conditions: -20°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		04/20/2011
<b>Review/Approval</b>	Staci Leffel	Manufacturing Manager		04/20/2011
<b>Review/Approval</b>	Richard Mann	Quality Assurance		04/20/2011