



CERTIFICATE OF CONFORMANCE

COC - 10301

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

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

**Lot Number:** M0020045

**Expiration Date:** 31 DEC 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       |
|------------------------|-----------------|-------------------------|------------------------|------------|
| <b>Preparer</b>        | Meredith Yeager | Manufacturing Associate | <i>Meredith Yeager</i> | 11/05/2010 |
| <b>Review/Approval</b> | Staci Leffel    | Manufacturing Manager   | <i>Staci Leffel</i>    | 11/03/2010 |
| <b>Review/Approval</b> | Richard Mann    | Quality Assurance       | <i>Richard Mann</i>    | 11/05/2010 |