



Document #: **QA-FM-24**

| | | | |
|--|-----------------|-------------------------------------|-----------------|
| Title: Certificate of Conformance | | | |
| Revision Code: B | Level: 3 | Effective Date: JAN 21, 2010 | Pages: 2 |

| Reviewed & Approved By | | | |
|-----------------------------------|------------------|------------------|-------------|
| Department/Function | Name | Signature | Date |
| Operations | Alan Kishbaugh | | |
| Quality Control | Robert Clegg | | |
| Kit Manufacturing | Staci Leffel | | |
| Scientific Support | Jenny Ly | | |
| Scientific Support | Pankaj Oberoi | | |
| Corporate Operations | Hans Biebuyck | | |
| Corporate Operations | George Sigal | | |
| Quality Assurance | Inderjit Kaushal | | |
| Author | Bob Clegg | | |

REVISION SUMMARY

| Revision Code | CR/CO # | Description of Changes | Initiated By |
|----------------------|----------------|---|---------------------|
| A | 09132 | Initial Release | Inderjit Kaushal |
| B | 02618 | Used new form template. Added product size and a reference to QA-FM-124 for statements to be included based on each product type. | Inderjit Kaushal |



CERTIFICATE OF CONFORMANCE

COC - 10466

Product Description: MSD[®] Diluent 2

Catalog #/Size: R51BB-2/200 mL; R51BB-3/40 mL; R51BB-4/8 mL

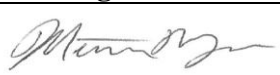

Lot Number: M0020076

Expiration Date: 28 FEB 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.

| | Name | Function | Signature | Date |
|------------------------|-----------------|-----------------------|---|------------|
| Preparer | Meredith Yeager | Inventory Technician |  | 12/13/2012 |
| Review/Approval | Staci Leffel | Manufacturing Manager |  | 12/13/2012 |