



CERTIFICATE OF CONFORMANCE

COC - 10235

Product Description: MSD® Diluent 2

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

Lot Number: M0020040

Expiration Date: 31 JUL 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 | | 03/07/2010 |
| Review/Approval | | K.I. Manufacturing | | 03/07/2010 |
| Review/Approval | | Quality Assurance | | 03/09/2010 |