



CERTIFICATE OF CONFORMANCE

COC - 10459

**Product Description:** MSD<sup>®</sup> Diluent 3

**Catalog #/Size:** R51BA-4 / 5 mL; R51BA-5 / 25 mL; R51BA-6 / 125 mL

**Lot Number:** M0030051

**Expiration Date:** 31 OCT 2014

**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        | Rolanda Spratley | Biologics/Chemistry   |           | 13 Sep 12  |
| Review/Approval | Staci Leffel     | Manufacturing Manager |           | 09/13/2012 |