



CERTIFICATE OF CONFORMANCE

COC - 10346

**Product Description:** MSD® Diluent 3  
**Catalog #/Size:** R51BA-5/25 mL; R51BA-6/125 mL  
**Lot Number:** M0030044  
**Expiration Date:** 31 JUL 2013  
**Statement:** This product was manufactured according to approved MSD documents.

Storage conditions: ≤ -10°C

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Preparer	Meredith Yeager	Manufacturing Associate		04/19/2011
Review/Approval	Staci Leffel	Manufacturing Manager		04/12/2011
Review/Approval	Richard Mann	Quality Assurance		04/19/2011