



CERTIFICATE OF CONFORMANCE

COC - 10360

Product Description: MSD® Diluent 3
Catalog #/Size: R51BA-4/5 mL; R51BA-5/25 mL
Lot Number: M0030045
Expiration Date: 31 AUG 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		06/01/2011
Review/Approval	Staci Leffel	Manufacturing Manager		05/31/2011
Review/Approval	Richard Mann	Quality Assurance		06/01/2011