



CERTIFICATE OF CONFORMANCE

COC - 10302

Product Description: MSD[®] Diluent 2
Catalog #/Size: 3-3027-020000/1700mL
Lot Number: M0020048
Expiration Date: 28 FEB 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		11/18/2010
Review/Approval	Staci Leffel	Manufacturing Manager		11/17/2010
Review/Approval	Richard Mann	Quality Assurance		11/18/2010