



CERTIFICATE OF CONFORMANCE

COC - 10231

Product Description: MSD® Diluent 3

Catalog #/Size: R51BA-6/125 mL

Lot Number: M0030031

Expiration Date: 31 JUL 2012

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	Janice Kim	Manufacturing Associate	<i>Jan Kim</i>	03/14/2010
Review/Approval	<i>Staci Leffel</i>	Kit Manufacturing mg	<i>Staci Leffel</i>	03/09/2010
Review/Approval	<i>Richard Mann</i>	Quality Assurance	<i>Richard Mann</i>	03/09/2010