



CERTIFICATE OF CONFORMANCE

COC - 10263

Product Description: MSD® Diluent 2

Catalog #/Size: R51BB-46/8 mL; R51BB-3/40 mL; R51BB-2/200 mL

Lot Number: M0020042

Expiration Date: 31 OCT 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		05/19/2010
Review/Approval		OPS		5/19/2010
Review/Approval	Richard Mann	QA		05/19/2010