



CERTIFICATE OF CONFORMANCE

COC - 10271

**Product Description:** MSD® Diluent 3

**Catalog #/Size:** N/A

**Lot Number:** M0030035

**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	<i>Janice Kim</i>	06/03/2010
Review/Approval	<i>Staci Leffel</i>	Kit Manufacturing Mgr	<i>Staci Leffel</i>	06/03/2010
Review/Approval	<i>Richard Mann</i>	QA	<i>Richard Mann</i>	06/03/2010