



CERTIFICATE OF CONFORMANCE

COC - 10249

**Product Description:** MSD® Diluent 3

**Catalog #/Size:** R51BA-4/5 mL

**Lot Number:** M0030033

**Expiration Date:** 30 SEP 2012

**Statement:** This product was manufactured according to approved MSD documents.

Storage conditions:  $\leq -10^{\circ}\text{C}$

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	Paul Goodwin	OPS		5/19/2010
Review/Approval	Richard Mann	QA		05/19/2010