

CERTIFICATE OF COMPLIANCE

COC-10214

Product Description: MSD[®] Diluent 2

Catalog #/Product Size: R51BB-4/8 mL or R51BB-3/40 mL

Lot Number: M0020039

Expiration Date: 30 JUN 2012

Statement: This diluent was manufactured at MSD[®] according to MSD[®] SOPs and specifications.

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

Prepared By:	Janice Kim	Date:	January 28, 2010
---------------------	------------	--------------	------------------

Reviewed and Approved By

Function	Name (Print)	Signature	Date
Operations	Staci Leffel	<i>Staci Leffel</i>	02/02/2010
Quality Assurance	Richard Mann	<i>Richard Mann</i>	02/02/2010