

CERTIFICATE OF COMPLIANCE

COC-10223

Product Description: MSD[®] Diluent 3

Catalog #/Product Size: R51BA-4/5 mL, R51BA-5/25 mL, or R51BA-6/125 mL

Lot Number: M0030028

Expiration Date: 31 MAY 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 Lettel	<i>Staci Lettel</i>	02/02/2010
Quality Assurance	Richard Mann	<i>Richard Mann</i>	02/02/2010