

CERTIFICATE OF COMPLIANCE

COC-10210

Product Description: Diluent 1

Catalog #/Product Size: R50CK-2/200 mL, R50CK-4/50 mL

Lot Number: M0010020

Expiration Date: 31 MAY 2010

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 8, 2010
---------------------	------------	--------------	-----------------

Reviewed and Approved By

Function	Name (Print)	Signature	Date
Operations	Staci Leffler	<i>Staci Leffler</i>	1/08/2010
Quality Assurance	Richard Mann	<i>Richard Mann</i>	01/11/2010