

CERTIFICATE OF COMPLIANCE

COC-10104

Product Description: Diluent 1

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

Lot Number: M0010023

Expiration Date: 30 SEP 2010

Statement: This diluent was manufactured according to approved MSD requirements.

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

Prepared By:	Elise Siegel	Date:	September 21, 2009
---------------------	--------------	--------------	--------------------

Reviewed and Approved By

Function	Name (Print)	Signature	Date
Quality	Inderjit Kaushal	<i>[Signature]</i>	09/24/09
Kit Manufacturing	Staci Leffel	<i>[Signature]</i>	09/24/2009