

CERTIFICATE OF COMPLIANCE

COC-10151

Product Description: Diluent 1

Catalog #/Product Size: R50CK-4/50 mL

Lot Number: M0010022

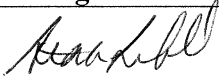

Expiration Date: 31 JUL 2010

Statement: This diluent was manufactured according to approved MSD Standard Operating Procedures.

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

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

Reviewed and Approved By

Function	Name (Print)	Signature	Date
Kit Manufacturing	Staci Leffel		10/24/2009
Quality Assurance	Richard Mann		10/26/2009