

CERTIFICATE OF COMPLIANCE

COC-10063

Product Description: MSD[®] Human Serum Cytokine Antibody Diluent

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

Lot Number: M0030009

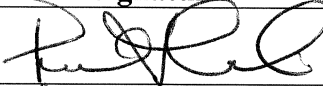

Expiration Date: December 2009

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:	Meagan McNamee	Date:	August 6, 2009
---------------------	-----------------------	--------------	-----------------------

Reviewed and Approved By

Function	Name (Print)	Signature	Date
Operations	Paul Goodwin		8/6/2009
Quality Assurance	Richard Mann		08/06/2009