

CERTIFICATE OF COMPLIANCE

COC-10027

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

Catalog #/Product Size: R51BB-4/8 mL, R51BB-3/40 mL, or R51BB-2/200 mL

Lot Number: M0020011

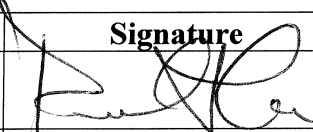

Expiration Date: November 2009

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:	July 30, 2009
---------------------	----------------	--------------	---------------

Reviewed and Approved By

Function	Name (Print)	Signature	Date
Operations	Paul Goodwin		7/31/09
Quality Assurance	Richard Mann		07/30/2009