

CERTIFICATE OF COMPLIANCE

COC-10064

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: M0020021



Expiration Date: March 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:	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		08/26/2009