



CERTIFICATE OF CONFORMANCE

COC - 10286

Product Description: Human Serum Cytokine Assay Diluent

Catalog #/Size: R51BB-3/40 mL; R51BB-2/200 mL

Lot Number: M0020025

Expiration Date: 30 APR 2010

Statement: This product was manufactured according to approved MSD documents.

Storage conditions: -20°C

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

	Name	Function	Signature	Date
Preparer	Meredith Yeager	Manufacturing Associate		07/30/2010
Review/Approval	Staci Leffel	Kit Manufacturing Manager		07/28/2010
Review/Approval	Richard Mann	QA		07/30/2010