



CERTIFICATE OF CONFORMANCE

COC - 00689

Product Description: MSD[®] Diluent 43
Catalog #/Size: R50AG-2/50 mL; R50AG-1/10 mL
Lot Number: M0430007
Expiration Date: 30 APR 2016
Statement: This product was manufactured according to approved MSD documents.

Storage conditions: ≤-10°C

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

	Name	Function	Signature	Date
Preparer	Meredith Yeager	Inventory Technician		05 MAR 2014
Review/Approval	Staci Leffel	Manufacturing Manager		21 APR 2014