



CERTIFICATE OF CONFORMANCE

COC - 00836

Product Description: MSD[®] Diluent 43
Catalog #/Size: R50AG-2/50 mL; R50AG-1/10 mL
Lot Number: M0430015
Expiration Date: 31 DEC 2017
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 Hutchinson	Production Control		27 SEP 2016
Review/Approval	Dawn Stollar	Quality Assurance		30 SEP 2016