



CERTIFICATE OF ANALYSIS

COA - 11326

Product Description: MSD[®] Diluent 17

Catalog #/Size: R50KA-3/30 mL

Lot Number: M0170044

Expiration Date: 31 MAY 2015

Statement: This product was manufactured and tested for interferents according to approved MSD documents.

Storage conditions: $\leq -10^{\circ}\text{C}$

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

	Name	Function	Signature	Date
Preparer	Meredith Yeager	Inventory Technician		03 Oct 2013
Review/Approval	Staci Leffel	Manufacturing Manager		03 Oct 2013