



CERTIFICATE OF ANALYSIS

COA - 06999

Product Description: MSD[®] Human (6E10) Abeta 42 Ultra-Sensitive Kit

Catalog #/Size: K151FUE-1 / 1 Plate; K151FUE-2 / 5 Plates; K151FUE-3 / 20 Plates

Lot Number: K0013415

Expiration Date: N/A (See expiration date on specific components below)

Statement: This product was manufactured and tested according to approved MSD documents. This lot meets MSD's specifications.

Kit Contents:

Plate Lot Number: Z0054683

Plate Lot Catalog Number: N451FSA

Plate Lot Expiration Date: 31 OCT 2016

Antibody blend Lot Number: D0012967

Antibody blend Catalog Number: D21FU

Antibody blend Expiration Date: 31 OCT 2016

Calibrator blend Lot Number: A0040072

Calibrator blend Catalog Number: C01LB

Calibrator blend Expiration Date: 31 OCT 2016

Specifications:

Release method: Coating uniformity



Mean Intraplate CV: $\leq 15\%$

Mean Interplate CV: $\leq 18\%$

No more than one plate with CV $> 18\%$

No plate with CV $> 20\%$

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

	Name	Function	Signature	Date
Preparer	Rolanda Spratley	Biologics/Chemistry		08 JUN 2016
Review/Approval	Vanessa Castleberry	Quality Assurance		08 JUN 2016