



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: MSD® Human (6E10) Abeta 40 Ultra-Sensitive Kit

Catalog Number: K151FTE - Series

Lot Number: K0013858

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 #	Z0055443	Catalog #	N451FSA	Expiration Date:	28 FEB 2018

Antibody Blend	
Lot Number:	D0013446
Catalog Number:	D21FT
Expiration Date:	31 JAN 2018
Storage Temperature:	2-8°C

Calibrator Blend	
Lot Number:	A0040118
Catalog Number:	C000A
Expiration Date:	31 OCT 2017
Storage Temperature:	≤ -70°C

Specifications:

Release method:

Coating Uniformity	Specifications
Mean Intraplate CV	≤ 15%
Mean Interplate CV	≤ 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
Review/Approval	Vanessa Castleberry	Quality Assurance	<i>Vanessa Castleberry</i>	06JUN2017