



# CERTIFICATE OF ANALYSIS

COA - 07256

**Product Description:** Human (6E10) Abeta 3-Plex Kit

**Catalog #/Size:** K15148E-1 / 1 Plate; K15148E-2 / 5 Plates; K15148E-3 / 20 Plates

**Lot Number:** K0013738

**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: Z0055187
- Plate Lot Catalog Number: N45148A
- Plate Lot Expiration Date: 31 MAR 2017
- Antibody blend Lot Number: D0013305
- Antibody blend Catalog Number: N/A
- Antibody blend Expiration Date: 30 APR 2017
- Calibrator blend Lot Number: A0010373
- Calibrator blend Catalog Number: C01XX
- Calibrator blend Expiration Date: 31 MAY 2017
- Calibrator blend Lot Number: A0010374
- Calibrator blend Catalog Number: C01XX
- Calibrator blend Expiration Date: 31 MAY 2017
- Calibrator blend Lot Number: A0010375
- Calibrator blend Catalog Number: C01XX
- Calibrator blend Expiration Date: 31 MAY 2017

### 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
<b>Preparer</b>	Xiomara Talavera	Quality Control	<i>Talavera</i>	30 AUG 2016
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	01 SEP 2016