



## CERTIFICATE OF ANALYSIS

COA - 05084

**Product Description:** MSD<sup>®</sup> Human a-Synuclein Kit

**Catalog #/Size:** K151TGD-1 / 1 Plate; K151TGD-2 / 5 Plates; K151TGD-4 / 25 Plates

**Lot Number:** K00E0171

**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: Z00E0150

Plate Lot Catalog Number: N451TGA

Plate Lot Expiration Date: 28 FEB 2017

Antibody blend Lot Number: D00E0160

Antibody blend Catalog Number: D21TG

Antibody blend Expiration Date: 28 FEB 2017

Calibrator blend Lot Number: A00E0088

Calibrator blend Catalog Number: C01TG

Calibrator blend Expiration Date: 31 MAR 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>	Rolanda Spratley	Biologics/Chemistry		28 JAN 2015
<b>Review/Approval</b>	Anastasia Belilovets	Quality Assurance		28 JAN 2015