



## CERTIFICATE OF ANALYSIS

COA - 06474

**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:** K00E0247

**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: Z00E0200

Plate Lot Catalog Number: N451TGA

Plate Lot Expiration Date: 30 APR 2018

Antibody blend Lot Number: D00E0228

Antibody blend Catalog Number: D21TG

Antibody blend Expiration Date: 30 APR 2018

Calibrator blend Lot Number: A00E0121

Calibrator blend Catalog Number: C01TG

Calibrator blend Expiration Date: 31 MAY 2018

### 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		25 FEB 2016
Review/Approval	Dawn Stollar	Quality Assurance		26 FEB 2016