



CERTIFICATE OF ANALYSIS

COA - 06616

**Product Description:** MSD® Human a-Synuclein Kit

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

**Lot Number:** K00E0191

**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: Z00E0170
- Plate Lot Catalog Number: N451TGA
- Plate Lot Expiration Date: 31 AUG 2017
- Antibody blend Lot Number: D00E0182
- Antibody blend Catalog Number: D21TG
- Antibody blend Expiration Date: 30 SEP 2017
- Calibrator blend Lot Number: A00E0098
- Calibrator blend Catalog Number: C01TG
- Calibrator blend Expiration Date: 31 AUG 2017

Specifications:

- Release method: Coating uniformity
- 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
<b>Preparer</b>	Rolanda Spratley	Biologics/Chemistry	<i>Rolanda Spratley</i>	09 MAR 2016
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	10 MAR 2016