



CERTIFICATE OF ANALYSIS

COA - 06416

Product Description: MSD[®] Human α -Synuclein Kit

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

Lot Number: K00E0182

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: Z00E0160

Plate Lot Catalog Number: N451TGA

Plate Lot Expiration Date: 30 JUN 2017

Antibody blend Lot Number: D00E0172

Antibody blend Catalog Number: D21TG

Antibody blend Expiration Date: 30 JUN 2017

Calibrator blend Lot Number: A00E0091

Calibrator blend Catalog Number: C01TG

Calibrator blend Expiration Date: 30 JUN 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
Preparer	Rolanda Spratley	Biologics/Chemistry		22 JAN 2016
Review/Approval	Dawn Stollar	Quality Assurance		27 JAN 2016