



CERTIFICATE OF ANALYSIS

COA - 07414

Product Description: MSD[®] PSD-95 Kit

Catalog #/Size: K150QND-1 / 1 Plate; K150QND-2 / 5 Plates; K150QND-4 / 25 Plates

Lot Number: K00E0177

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: Z00E0155

Plate Lot Catalog Number: N45QNA

Plate Lot Expiration Date: 30 APR 2017

Antibody blend Lot Number: D00E0165

Antibody blend Catalog Number: D20QN

Antibody blend Expiration Date: 30 APR 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	Xiomara Talavera	Quality Control		14 OCT 2016
Review/Approval	Dawn Stollar	Quality Assurance		17 OCT 2016