



CERTIFICATE OF ANALYSIS

COA - 03927

Product Description: MSD® MA6000 Human Insulin Kit
MSD® Human Insulin Kit

Catalog #/Size: K111BZC-1 / 1 Plate; K111BZC-2 / 5 Plates; K111BZC-3 / 20 Plates
K151BZC-1 / 1 Plate; K151BZC-2 / 5 Plates; K151BZC-3 / 20 Plates

Lot Number: K0012988

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: Z0054086
Plate Lot Catalog Number: L451BZA
Plate Lot Expiration Date: 31 JUL 2013
Antibody blend Lot Number: D0012535
Antibody blend Catalog Number: D21BZ
Antibody blend Expiration Date: 31 JUL 2013
Calibrator blend Lot Number: A0010292
Calibrator blend Catalog Number: C00BZ
Calibrator blend Expiration Date: 15 Months From Date of Receipt

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.

Table with 5 columns: Name, Function, Signature, Date. Rows include Preparer (Rolanda Spratley) and Review/Approval (Dana Short).