

CERTIFICATE OF ANALYSIS

COA - 06463

Product Description: MSD® Mouse/Rat Insulin Kit

Catalog #/Size: K152BZC-1 / 1 Plate; K152BZC-2 / 5 Plates; K152BZC-3 / 20 Plates

Lot Number: K0013856

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: Z0055465

Plate Lot Catalog Number: L452BZA
Plate Lot Expiration Date: 31 MAR 2018
Antibody blend Lot Number: D0013444
Antibody blend Catalog Number: D22BZ
Antibody blend Expiration Date: 28 FEB 2018
Calibrator blend Lot Number: A0010380
Calibrator blend Catalog Number: C00BZ
Calibrator blend Expiration Date: 30 NOV 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	Bo-Jodon Falley	08 FEB 2016
Review/Approval	Dawn Stollar	Quality Assurance	Dawn Stollar	08 FEB 2016