



CERTIFICATE OF ANALYSIS

COA - 06728

Product Description: MSD® Phospho/Total ERK1/2 Whole Cell Lysate Kit
Catalog #/Size: K15107D-1 / 1 Plate; K15107D-2 / 5 Plates; K15107D-3 / 20 Plates
Lot Number: K0013810
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: Z0055336
Plate Lot Catalog Number: N45107B
Plate Lot Expiration Date: 30 SEP 2017
Antibody blend Lot Number: D0013397
Antibody blend Catalog Number: D21CM
Antibody blend Expiration Date: 30 SEP 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.

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