



RFE MEMO

Page 1 of 1

Product Description(s): Total Akt Kit
Catalog Number(s) / Size(s): K150MOD-Series
Lot Number: K0014232

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0014232 has been extended to 31 MAY 2021.

The expiration dates for the following kit components have also been extended:

Description	Lot Number	Original Expiration Date	New Expiration Date
Akt Plate	Z0056118	31 MAY 2020	31 MAY 2021
Akt Calibrator	A0010452	30 NOV 2020	31 MAY 2021
Anti-Total Akt Antibody	D0013861	31 MAY 2020	31 MAY 2021

Function	Name	Signature	Date
Quality	Karen Hamilla		25 MAR 2020

This memo supersedes all other statements regarding date of expiration for this product.

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: Total Akt Kit
Kit Catalog Numbers: K150MOD-Series
Kit Lot Number: K0014232
Expiration Date: 31 May 2020

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Akt Control 1		Akt Control 2	
		Recovery	CV	Recovery	CV	Recovery	CV
1	Akt	80% - 120%	< 20%	70% - 130%	< 20%	70% - 130%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	Akt	80X Calibrator	4104	ng/mL
		Akt Control 1	16.8	ng/mL
		Akt Control 2	0.985	ng/mL

Controls tested may not be the same lot as shipped.

Concentrations provided in this table refer to the stock concentration of the calibrators. For top of the standard curve concentrations, please refer to the Standard Curve Data

Standard Curve Data:

Akt - Spot 1		
Conc. ng/mL	ECL Signal	%CV
51.3	569,000	3.0
12.8	166,000	3.7
3.21	36,500	4.4
0.802	8,010	4.9
0.200	1,950	1.9
0.0501	533	3.2
0.0125	179	4.7
0.00	66	9.8

CERTIFICATE OF ANALYSIS

Page 2 of 2

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Akt Plate	Z0056118	2-8°C	31 May 2020
Anti-Total Akt Antibody	D0013861	2-8°C	31 May 2020
Akt Calibrator	A0010452	≤ -70°C	30 Nov 2020
Blocker A	Not Kit Specific	Not Kit Specific	N/A
Diluent 39	Not Kit Specific	Not Kit Specific	N/A
MSD Tris Wash Buffer	Not Kit Specific	Not Kit Specific	N/A
Phosphatase Inhibitor I	Not Kit Specific	Not Kit Specific	N/A
Phosphatase Inhibitor II	Not Kit Specific	Not Kit Specific	N/A
Protease Inhibitor Solution	Not Kit Specific	Not Kit Specific	N/A
MSD Read Buffer T	Not Kit Specific	Not Kit Specific	N/A

Statement:

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	28 AUG 2018