



RFE MEMO

Page 1 of 1

Product Description(s): Phospho-Akt (Ser473) Kit

Catalog Number(s) / Size(s): K150MND - Series

Lot Number: K0014328

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0014328 has been extended to 30 NOV 2021.

The expiration date for the following kit component has been extended:

Description	Lot Number	Original Expiration Date	New Expiration Date
Akt Plate	Z0056281	30 NOV 2020	30 NOV 2021
Akt Calibrator	A0010452	30 NOV 2020	30 NOV 2021

Function	Name	Signature	Date
Quality	Karen Hamilla		04 NOV 2020

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

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: Phospho-Akt(Ser473) Kit
Kit Catalog Numbers: K150MND-Series
Kit Lot Number: K0014328
Expiration Date: 30 Nov 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	20X Calibrator	316	UP473/well
		Akt Control 1	3.77	UP473/well
		Akt Control 2	0.204	UP473/well

Control data was generated utilizing internal reference controls

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. UP473/well	ECL Signal	%CV
15.8	72,100	5.5
3.95	24,100	3.9
0.988	4,520	3.4
0.247	942	2.8
0.0617	301	2.4
0.0154	149	4.2
0.00386	113	7.4
0.00	101	4.5

CERTIFICATE OF ANALYSIS

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Akt Plate	Z0056281	2-8°C	30 Nov 2020
Anti-Phospho-Akt(Ser473) Antibody	D0013958	2-8°C	30 Nov 2021
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
Tris Wash Buffer (10X)	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	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		14 SEP 2018