

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: Phospho-Akt (Ser473) Kit
Kit Catalog Numbers: K150MND-Series
Kit Lot Number: K0014302
Expiration Date: 31 Oct 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 | 332 | Up473/well |
| | | Akt Control 1 | 3.77 | Up473/well |
| | | Akt Control 2 | 0.204 | Up473/well |

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. Up473/well | ECL Signal | %CV |
| 16.6 | 88,400 | 6.5 |
| 4.15 | 27,100 | 2.2 |
| 1.04 | 5,060 | 4.3 |
| 0.259 | 1,090 | 3.4 |
| 0.0648 | 342 | 3.4 |
| 0.0162 | 164 | 5.0 |
| 0.00405 | 116 | 3.0 |
| 0.00 | 103 | 3.6 |

CERTIFICATE OF ANALYSIS


Page 2 of 2

Components:

| Description | Lot Number | Storage Temperature | Expiration Date |
|------------------------------------|------------------|---------------------|-----------------|
| Akt Plate | Z0056250 | 2-8°C | 31 Oct 2020 |
| Anti-Phospho-Akt (Ser473) Antibody | D0013930 | 2-8°C | 31 Oct 2021 |
| Akt Calibrator | A0010456 | ≤ -70°C | 31 Dec 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 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 |  | 13 AUG 2018 |