

# AlphaLISA SIRT1 p53 Lysine 382 Deacetylase Assay

AlphaLISA #10

AlphaLISA®

## Authors

Mathieu Arcand  
Julie Blouin  
Mireille Caron  
Claire Normand  
Anne Labonté  
Lucille Beaudet  
Jaime Padrós

PerkinElmer, Inc.  
Montreal, QC  
Canada, H3J 1R4

This AlphaLISA immunodetection assay measures the deacetylation of a biotinylated p53 (368-393) peptide acetylated at lysine 382.

### Anti-acetyl-p53 Lysine 382 (p53K382ac) AlphaLISA® Acceptor Beads

- AL124C: 250 µg, 500 assay points\*
- AL124M: 5 mg, 10,000 assay points\*
- AL124R: 25 mg, 50,000 assay points\*

\*0.5 µg/assay point

### Peptidic Substrate Sequence:

(Biotin)K-GG-HLKSKKGQSTSRHKK(ac)LMFKTEGPDS-D-NH<sub>2</sub>

### AlphaLISA Assays

AlphaLISA technology is a powerful and versatile platform that offers highly sensitive, no-wash immunoassays using Alpha Streptavidin Donor and AlphaLISA Acceptor beads. In this technical note, we present the optimization of a signal decrease SIRT1 assay using as substrate a biotinylated p53-derived peptide acetylated at lysine 382. In the absence of enzyme or cofactor, the anti-p53K382ac Acceptor beads bind the acetylated residue on the peptide. Upon laser irradiation of the beads-target complexes at 680 nm, short-lived singlet oxygen molecules produced by the Donor beads can reach the Acceptor beads in proximity to generate maximal AlphaLISA signal at 615 nm (left panel). When enzyme and cofactor are added to the reaction, the peptide substrate is deacetylated and the anti-p53K382ac Acceptor beads do not recognize the biotinylated peptide anymore, leading to a signal decrease (right panel). This signal decrease is proportional to the deacetylation activity of the SIRT1 enzyme.

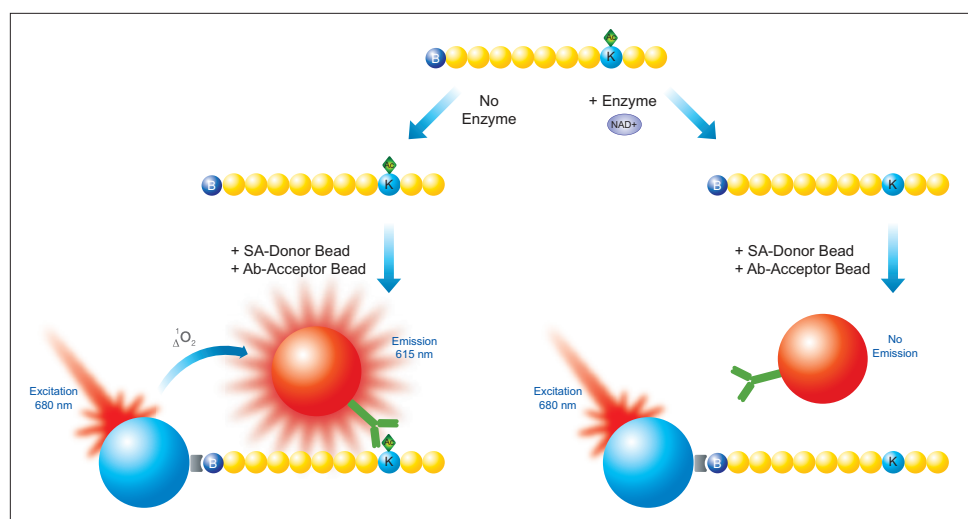


Figure 1. Schematic representation of the AlphaLISA detection of a modified p53-derived peptide.

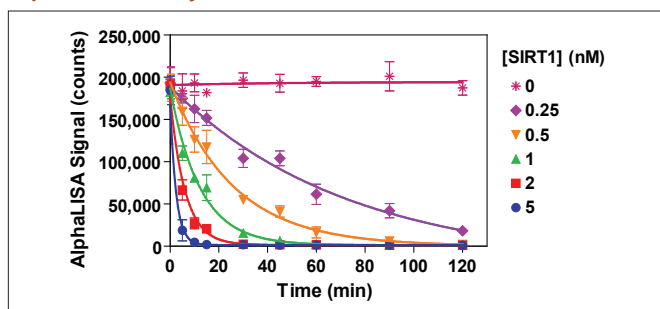
## Development of a SIRT1 p53 Lysine 382 Deacetylase Assay:

### Reagents needed for the assay:

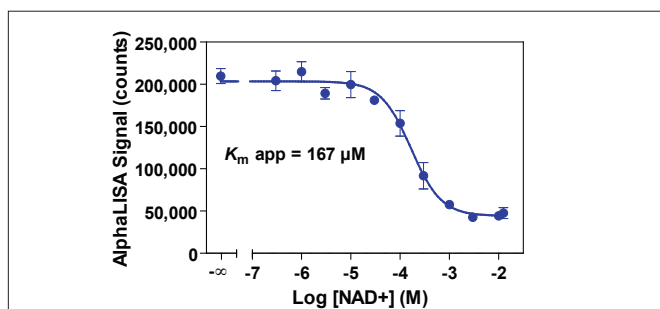
Anti-acetyl-p53 Lysine 382	
AlphaLISA Acceptor beads	PerkinElmer # AL124
p53 (368-393) acetyl-lysine 382 peptide (p53K382ac), biotinylated	AnaSpec # 64869
Alpha Streptavidin Donor beads	PerkinElmer # 6760002
AlphaLISA 5X Epigenetics Buffer 1 Kit	PerkinElmer # AL008
Sirtuin 1 (human), recombinant	BPS BioScience # 50012
EX-527	Tocris # 2780
Suramin	Calbiochem # 574625
SIRT1 inhibitor III	Calbiochem # 566322
Nicotinamide	Sigma # N3376
$\beta$ -Nicotinamide adenine dinucleotide hydrate (NAD <sup>+</sup> )	Sigma # N1636
White opaque OptiPlate™-384	PerkinElmer # 6007299
TopSeal™-A films	PerkinElmer # 6005185

Assay Buffer: 50 mM Tris-HCl pH 8.0, 150 mM NaCl, 1 mM DTT, 0.01% Tween-20 and 0.01% BSA.

### Experiment 1: Enzyme Titration and Time-Course



### Experiment 2: NAD<sup>+</sup> Titration

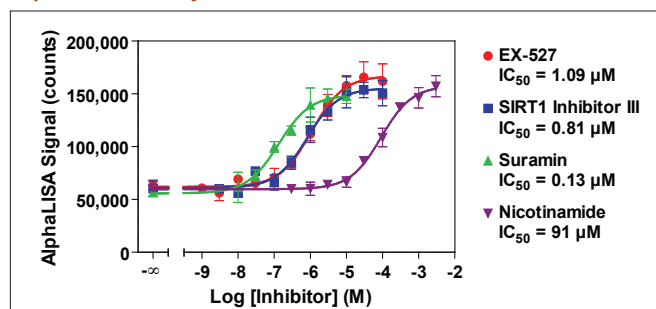


PerkinElmer, Inc.  
940 Winter Street  
Waltham, MA 02451 USA  
P: (800) 762-4000 or  
(+1) 203-925-4602  
[www.perkinelmer.com](http://www.perkinelmer.com)

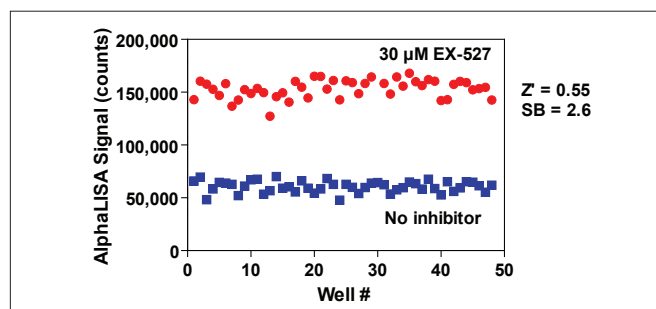
### Standard Protocol

- Dilute SIRT1 enzyme, inhibitors, biotinylated p53K382ac peptide substrate and NAD<sup>+</sup> in Assay Buffer just before use.
- Add to the wells of a white OptiPlate-384:
  - 2.5  $\mu$ L of enzyme (4X)
  - 2.5  $\mu$ L of inhibitor (4X) or Assay buffer
  - Incubate 5 min at RT
  - 2.5  $\mu$ L of biotinylated p53K382ac peptide substrate (4X)
  - 2.5  $\mu$ L of NAD<sup>+</sup> (4X)
- Cover the plate with TopSeal-A film and incubate at room temperature (RT).
- Prepare 1X Epigenetics Buffer 1 as recommended in the buffer technical data sheet.
- Prepare a 5X Stop Solution Mix containing 250  $\mu$ M of EX-527 and 100  $\mu$ g/mL of Acceptor Beads in 1X Epigenetics Buffer 1 (final concentration of 50  $\mu$ M EX-527 and 20  $\mu$ g/mL Acceptor Beads in 25  $\mu$ L total assay volume).
  - 5  $\mu$ L of Stop Solution Mix
- Cover with TopSeal-A film and incubate for 60 min at RT.
- Prepare a 2.5X Streptavidin Donor beads solution at 50  $\mu$ g/mL in 1X Epigenetics Buffer 1 (final concentration of 20  $\mu$ g/mL in 25  $\mu$ L total assay volume) in subdued light.
  - 10  $\mu$ L of Streptavidin Donor beads
- Cover with TopSeal-A film and incubate in subdued light for 30 min at RT.
- Read signal in Alpha mode with the EnVision® or EnSpire® reader.

### Experiment 3: Enzyme Inhibition



### Experiment 4: Z'-factor Determination



For a complete listing of our global offices, visit [www.perkinelmer.com/ContactUs](http://www.perkinelmer.com/ContactUs)

Copyright © 2010-2011, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.