

## Certificate of Analysis

### sirtuin 2

#### NAD-dependent protein deacetylase sirtuin-2 isoform 1 (Human)

(Recombinant protein expressed in *E.coli*)

Item # EPI060

Lot # 139684

**Product Description:** N-terminal 6His-tagged, recombinant, amino acids 50-356, human sirtuin 2, expressed in *E.coli*. Purified using immobilised metal affinity chromatography.  
MW = 37.3kDa.

**Formulation:** 1mg/ml of protein in 25mM Hepes/NaOH pH7.4, 50mM NaCl, 50% glycerol. Frozen solution.

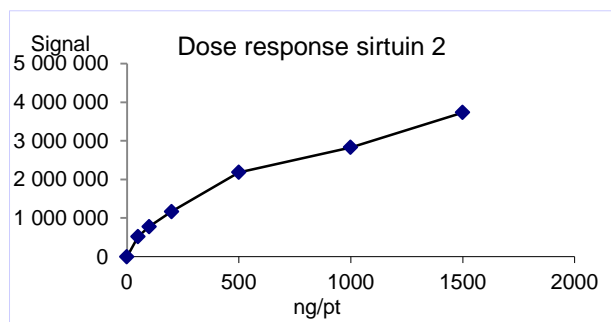
**Storage and Stability:** Stable for 1 year at -70°C from date of shipment. For maximum recovery of product, centrifuge original vial prior to removing the cap.

**Handling Recommendations:** Rapidly thaw the vial under cold water and immediately place on ice. Aliquot unused material into pre-chilled microcentrifuge tubes and store at -70°C.

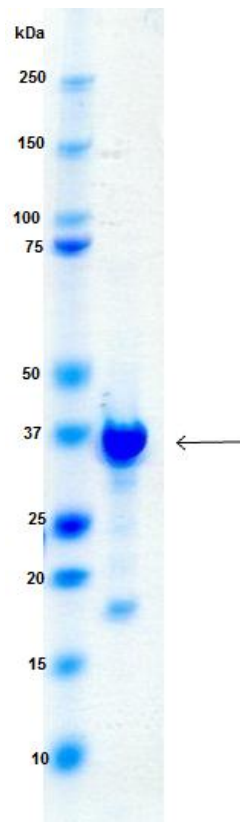
**FOR IN VITRO RESEARCH USE ONLY  
NOT FOR USE IN HUMANS OR ANIMALS**

### Quality Control Testing

**Enzymatic Assay:** 50-1500ng of this lot of enzyme deacetylated 150µM Fluor de Lys-SIRT2, deacetylase substrate in the assay described on page two. Assay background was subtracted from the actual counts to yield the results shown below.



**MS:** Size was confirmed by mass spectrometry using a Q-TOF.



**SDS-PAGE and Coomassie Stain:** Purity was assessed by SDS-PAGE and Coomassie blue staining using 4µg of sirtuin 2

## Certificate of Analysis

### sirtuin 2 Assay Protocol

#### Stock Solutions:

1. **Reaction buffer:** 50mM Tris/HCl pH8, 137mM NaCl, 2.7mM KCl, 1mM MgCl<sub>2</sub>, 0.2% BSA.
2. **sirtuin 2:** Dilute with reaction buffer. Use 50-1500ng per assay point.
3. **β-NAD:** Dilute with reaction buffer to 800μM.
4. **Fluor de Lys-SIRT2, Deacetylase Substrate:** Dilute with β-NAD solution to 300μM.

#### Assay Procedure (384 well plate format):

1. Add 2μl of H<sub>2</sub>O per assay to each well.
2. Add 8μl (**50-1500ng**) **sirtuin 2**.
3. Add 10μl of Fluor de Lys-SIRT2, deacetylase substrate/β-NAD.
4. Incubate for 60 minutes at 22°C.
5. Add 20μl of HDAC Developer.
6. Incubate for 15 minutes at 22°C.
7. Read at 460nm after an excitation flash at 355nm. Compare signal of enzyme samples with value of Back Ground that contains all assay components except sirtuin 2.

## Certificate of Analysis

### sirtuin 2 Sequence Information

**Protein** Human sirtuin 2  
**Tags** N-terminal 6His  
**Accession number** GenBank NP\_036369.2

***Recombinant sirtuin 2 amino acid sequence:***

1 MHHHHHSSG VDLGTENLYF QSMSLGSQKE RLLDELTLEG VARYMQSERC  
51 RRVICLVGAG ISTSAGIPDF RSPSTGLYDN LEKYHLPYPE AIFEISYFKK  
101 HPEPFFALAK ELYPGQFKPT ICHYFMRLK DKGLLLRCYT QNIDTLERIA  
151 GLEQEDLVEA HGTFYTSHCV SASCRHEYPL SWMKEKIFSE VTPKCEDCQS  
201 LVKPDIVFFG ESLPARFFSC MQSDFLKVDL LLVMGTSLQV QPFASLISKA  
251 PLSTPRLIN KEKAGQSDPF LGMIMGLGGG MDFDSKKAYR DVAWLGEDQ  
301 GCLALAELLG WKKELEDLVR REHASIDAQS

Reviewed and approved by site quality representative.

Unless otherwise stated in our catalogue or other company documentation accompanying the product(s), our products are intended for research use only and are not to be used for any other purpose, which includes but is not limited to, unauthorized commercial uses, in vitro diagnostic uses, ex vivo or in vivo therapeutic uses or any type of consumption or application to humans or animals.

© 2014 Eurofins Pharma Discovery Services UK Limited is an independent member of Eurofins Discovery Services.