

AlphaLISA® SureFire® Ultra™

## Human and Mouse p- $\alpha$ -Synuclein (Ser129) Aggregate Detection Kit

**Product number:** ALSU-APASYN-A500, ALSU-APASYN-A10K,  
 ALSU-APASYN-A50K, ALSU-APASYN-A-HV



### Kit specificity:

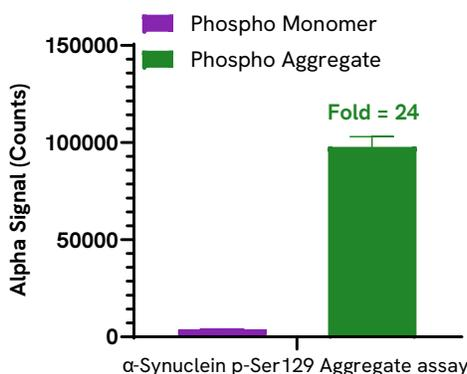
This assay kit contains antibodies which recognize the phospho-Ser129 epitope on aggregated  $\alpha$ -Synuclein. The protein detected by this kit corresponds to UniProt ID P37840.  $\alpha$ -Synuclein is also known as Non-A  $\beta$  component of AD amyloid. These antibodies recognize  $\alpha$ -Synuclein aggregate of human and mouse origin. Other species should be tested on a case-by-case basis.

### Control lysate information:

Positive Control: Prepared from recombinant human  $\alpha$ -Synuclein aggregate (active, phosphorylated) at a concentration of 1  $\mu$ g/mL (Abcam ab218819).

### Representative data:

Specificity of the  $\alpha$ -Synuclein Phospho (Ser129) Aggregate assay was assessed by assaying aggregated and monomeric proteins. Phosphorylated  $\alpha$ -Synuclein aggregate (Abcam ab218819) and phosphorylated  $\alpha$ -Synuclein monomer were prepared at approximately 250 ng/mL in Lysis Buffer and evaluated using the Phospho (Ser129) Aggregate AlphaLISA SureFire Ultra assay.  $\alpha$ -Synuclein Phospho (Ser129) Aggregate assay recognizes phosphorylated  $\alpha$ -Synuclein protein only when it is aggregated.



The information provided in this document is for reference purposes only and may not be all-inclusive. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume liability for the accuracy or completeness of the information contained herein. Users should exercise caution when handling materials as they may present unknown hazards. Revvity shall not be liable for any damages or losses resulting from handling or contact with the product, as Revvity cannot control actual methods, volumes, or conditions of use. Users are responsible for ensuring the product's suitability for their specific application. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF"