

AlphaLISA® SureFire® Ultra™

Human and Mouse p- α -Synuclein (Ser129) Detection Kit

Product number: ALSU-PASYN-B500, ALSU-PASYN-B10K,
 ALSU-PASYN-B50K, ALSU-PASYN-B-HV



Kit specificity:

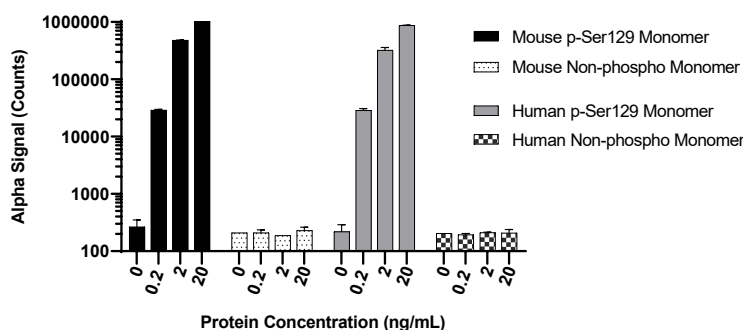
This assay kit contains antibodies which recognize the phospho-Ser129 epitope and a distal epitope on α -Synuclein. The protein detected by this kit corresponds to UniProt ID P37840. α -Synuclein is also known as Non-A β component of AD amyloid. These antibodies recognize α -Synuclein of human and mouse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

Control lysate information:

Positive Control: Prepared from recombinant human α -Synuclein aggregate (Active) at a concentration of 1 μ g/mL (Abcam catalogue number: ab218819).

Representative data

Data obtained with a 2-plate, 2-incubation protocol. Recombinant mouse and human Phospho (Ser129) and non-phosphorylated α -Synuclein monomers* were diluted with Lysis Buffer at the indicated concentrations and assayed for Phospho (Ser129) α -Synuclein using the SureFire Ultra kit. This assay has been validated in wildtype mouse brain tissue lysate, as well as lysate from PLK3 transfected HEK293 cells to promote significant phosphorylation of human α -Synuclein at Ser129 (unpublished data*). No assay signal was obtained using corresponding α -Synuclein knockout mouse brain tissue or HEK293 cell lysate (unpublished data*). *Recombinant proteins and unpublished data provided by Prof. Deniz Kirik and Dr. Benjamin Trist from The University of Sydney, Australia.



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"