

Technical Data Sheet

Caution: Research use only.

AlphaLISA® SureFire® Ultra™ α-Synuclein Aggregate-Specific Assay Kit

**Products: ALSU-AASYN-A500, ALSU-AASYN-A10K, ALSU-AASYN-A50K,
ALSU-AASYN-A-HV**

Kit-Specificity Information

This assay kit contains antibodies, recognising distinct epitopes specific to the Aggregated form of α-Synuclein. The human protein detected by this kit corresponds to Entre Gene ID 6622.

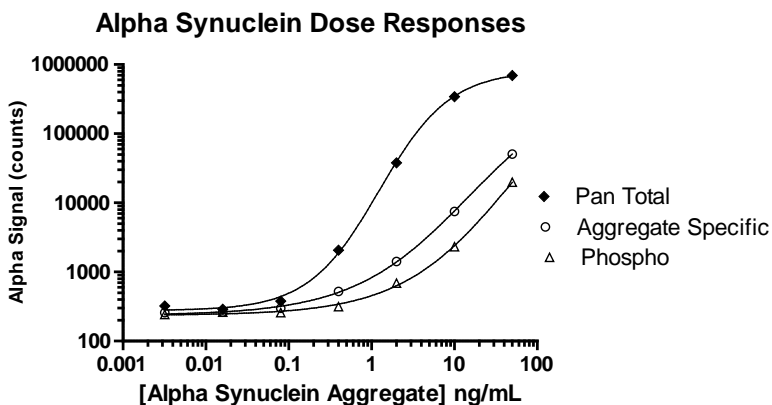
These antibodies recognize α-Synuclein of human, mouse and rat origin. Other species should be tested on a case-by-case basis.

Positive Control Information

Positive Control: Prepared from recombinant human Alpha-synuclein protein aggregate (Active) at a concentration of 1μg/ml (Abcam cat no. ab218819).

Representative data

Data obtained with a 2-plate, 2-incubation protocol. Recombinant α-Synuclein aggregate was diluted to the concentrations shown with 1X Lysis Buffer, and assayed for α-Synuclein Total, α-Synuclein Aggregate-Specific and p-α-Synuclein (Ser129) using respective *SureFire Ultra* kits.



For more information and detailed protocols, please refer to the generic AlphaLISA *SureFire Ultra* Assay Kits Manual, available from <http://www.perkinelmer.com/category/alpha-surefire-kits>.