

AlphaLISA® SureFire® Ultra™

## Human and Mouse DJ1 Total Detection Kit

**Product number:** ALSU-TDJ1-A500, ALSU-TDJ1-A10K,  
ALSU-TDJ1-A50K, ALSU-TDJ1-A-HV



### **Kit specificity:**

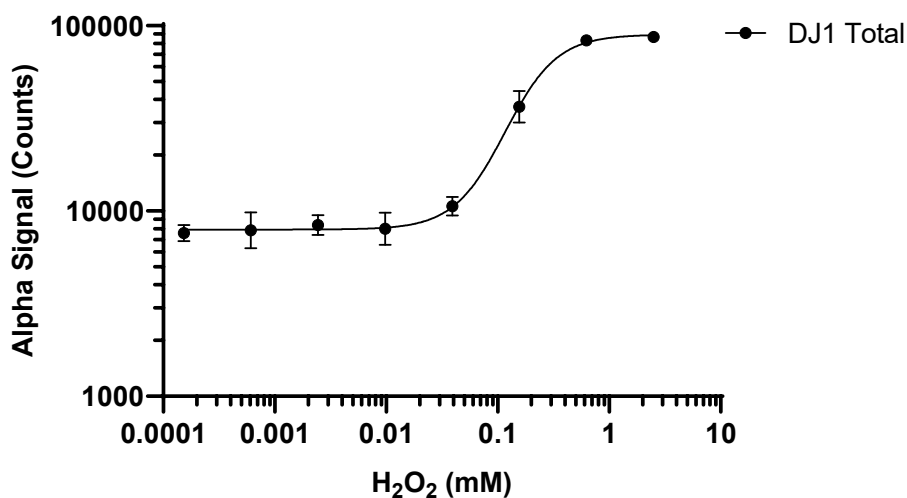
This assay kit contains antibodies which recognize distinct epitopes on DJ1. The protein detected by this kit corresponds to UniProt ID Q99497. DJ1 is also known as Parkinson disease protein 7 (PARK7) and Parkinsonism-associated deglycase. These antibodies recognize DJ1 of human and mouse origin. Other species should be tested on a case-by-case basis.

### **Control lysate information:**

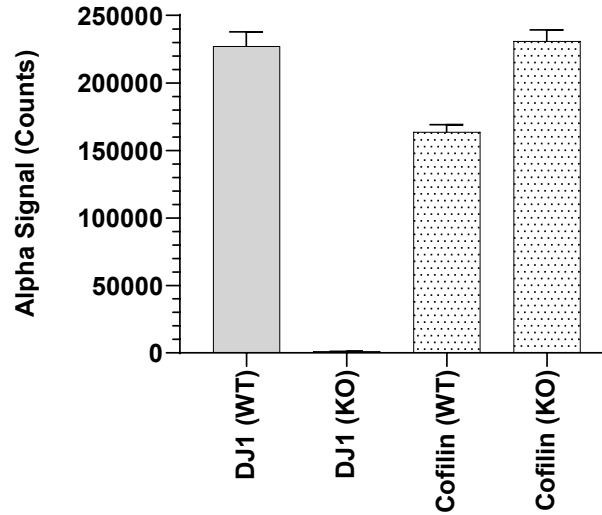
Positive Control Lysate: Prepared from HEK293 cells, cultured to confluence in T175 flasks in 10% FBS containing medium, treated with 22 mM H<sub>2</sub>O<sub>2</sub> for 10 minutes and lysed with 5 mL of Lysis Buffer.

### **Representative data**

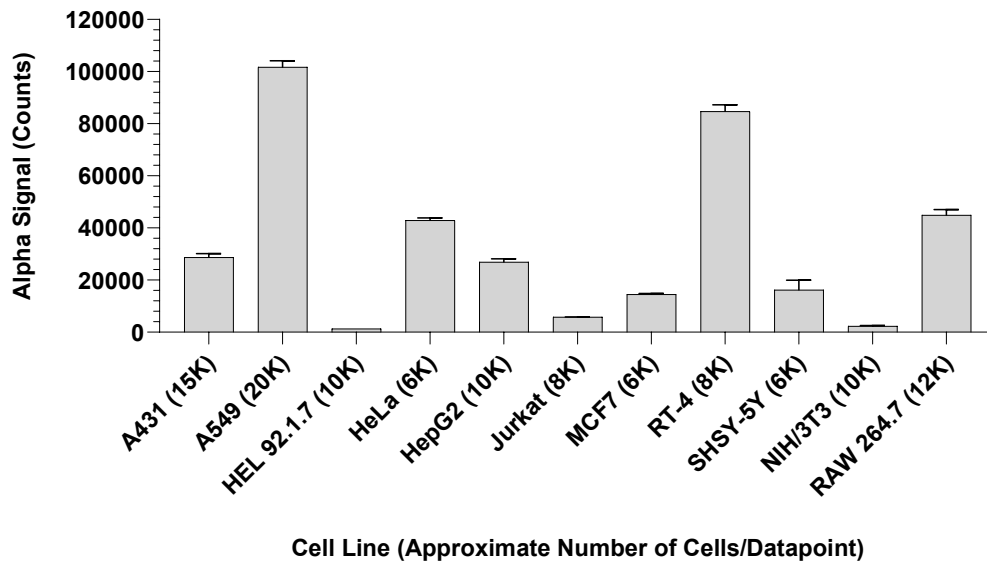
Data obtained with a 2-plate, 2-incubation protocol. MCF7 cells were seeded at 40K cells/well in a 96 well plate and incubated overnight. Cells were treated with H<sub>2</sub>O<sub>2</sub> at the indicated concentrations for 10 minutes. Cells were lysed with Lysis Buffer B and assayed for Total DJ1 using the *SureFire Ultra* kit. Equivalent to approximately 4,000 cells/datapoint.



Data obtained with a 2-plate, 2-incubation protocol. HEK293 DJ1-KO and HEK293 DJ1-WT cells were cultured to confluence in T75 flasks in 10% FBS containing medium. Cells were lysed with Lysis Buffer and assayed separately for Total DJ1 and Cofilin using respective *SureFire Ultra* kits. Equivalent to approximately 5,000 cells/datapoint for DJ1 and approximately 100 cells/datapoint for Cofilin.



Data obtained from measurement of CDK8 Total in various cell lines lysed with Lysis Buffer. Approximate number of cells/datapoint is indicated.



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"

