

AlphaLISA® SureFire® Ultra™

Human and Mouse GSPT1 Total Detection Kit

Product number: ALSU-TGSPT1-A500, ALSU-TGSPT1-A10K,
ALSU-TGSPT1-A50K, ALSU-TGSPT1-A-HV



Kit specificity:

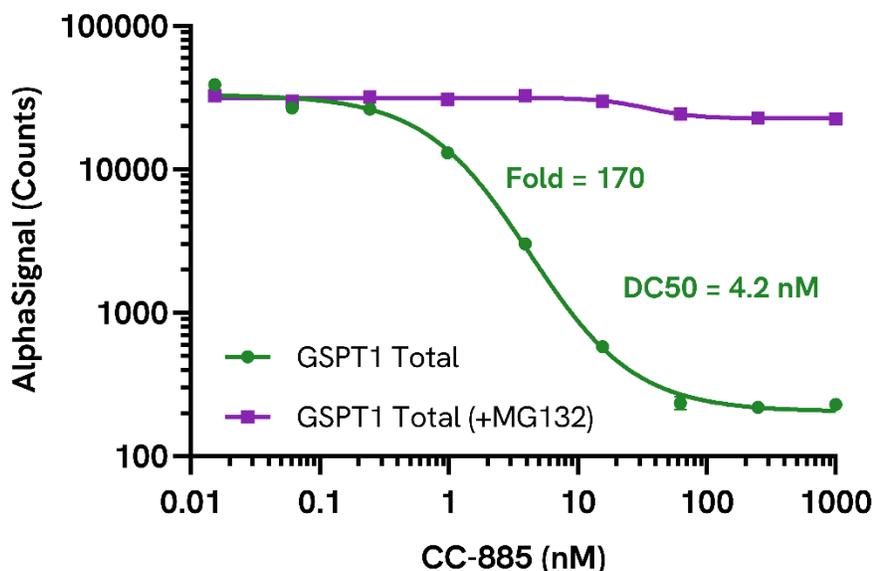
This assay kit contains antibodies which recognize distinct epitopes on GSPT1. The protein detected by this kit corresponds to UniProt ID P15170. GSPT1 is also known as eRF3a and G1 to S phase transition protein 1 homolog. These antibodies recognize GSPT1 of human and mouse origin. Other species should be tested on a case-by case basis.

Control lysate information:

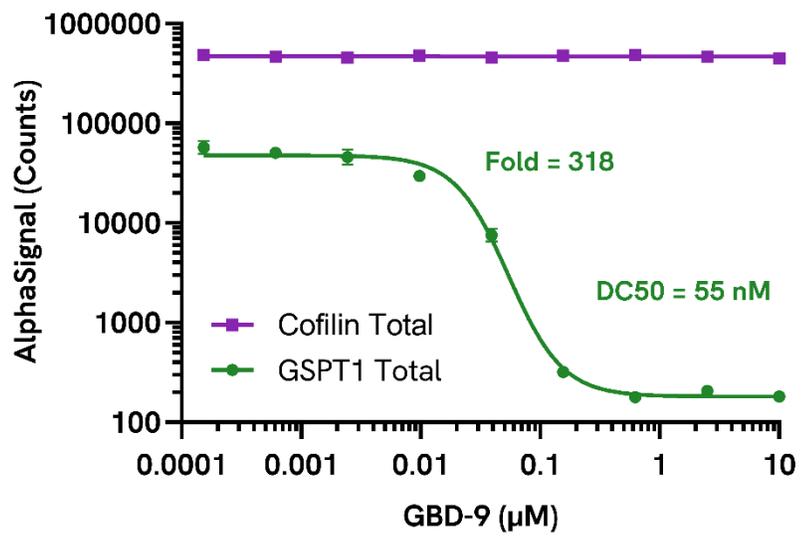
Positive Control Lysate: Prepared from HeLa cells, cultured to confluence in T175 flasks in 10% FBS containing medium and lysed with 10 mL of Lysis Buffer.

Representative data:

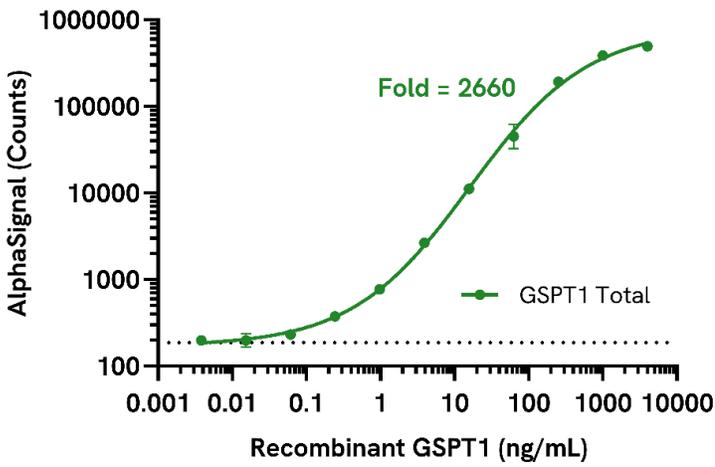
Data obtained with a 2-plate, 2-incubation protocol. MOLT-4 cells were harvested, washed with HBSS + 0.1% BSA and seeded at 250K cells/well in a 96 well plate. Cells were pretreated with 5 μM MG132 for 1 hour and treated with increasing concentrations of CC-885 for 2 hours. Cells were lysed with 5X Lysis Buffer and assayed for Total GSPT1 using the *SureFire Ultra* kit. Equivalent to approximately 10,000 cells/datapoint.



Data obtained with a 2-plate, 2-incubation protocol. Ramos cells were seeded at 100K cells/well in a 96 well plate in complete RPMI and treated with increasing concentrations of GBD-9 for 16 hours. The cells were washed with HBSS and then lysed with Lysis Buffer. Lysates were assayed separately for Total GSPT1 and Total Cofilin using respective *SureFire Ultra* kits. Equivalent to approximately 10,000 cells/datapoint.



Dilutions of recombinant GSPT1 protein (Abcam, ab12287) were prepared in Lysis Buffer and assayed using the *SureFire Ultra* GSPT1 Total kit.



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"

