

Technical Data Sheet

AlphaLISA® SureFire® Ultra™ p-GCN2 (Thr899) Assay Kit

ALSU-PGCN2-A500, ALSU-PGCN2-A10K, ALSU-PGCN2-A50K, ALSU-PGCN2-A-HV

Kit Specificity

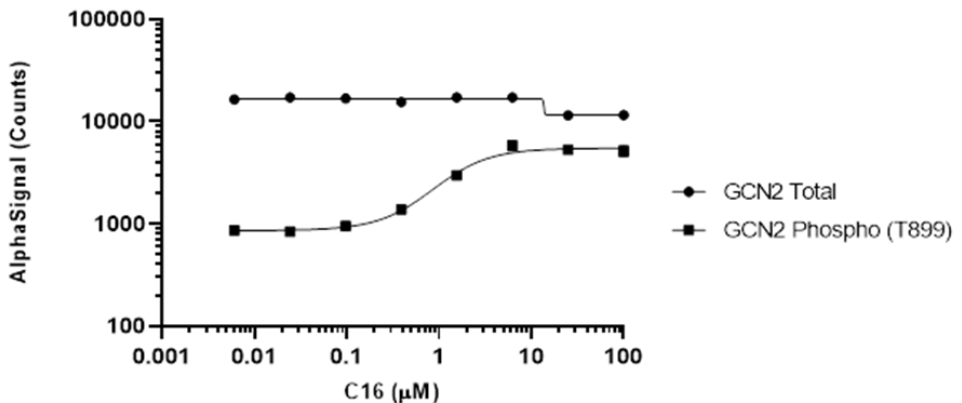
This assay kit contains antibodies which recognize the phospho-Thr899 epitope and a distal epitope on GCN2. The protein detected by this kit corresponds to UniProt ID Q9P2K8. GCN2 is also known as eIF-2-alpha kinase GCN2 and EIK2AK4. These antibodies recognize GCN2 of human origin, other species should 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, then treated with 20 μ M C16 for 15 minutes and lysed in 5 mL of 1X *SureFire Ultra* Lysis Buffer.

Representative Data

Data obtained with a 2-plate, 2-incubation protocol. AN3CA cells were plated at 40K cells per well in a 96 well plate and incubated overnight. Cells were then treated with C16 at the indicated concentrations for 15 minutes. Cells were lysed with 1X *SureFire Ultra* Lysis Buffer and assayed separately for Total or Phospho (Thr899) GCN2, using respective *SureFire Ultra* assay kits. Equivalent to approximately 4000 cells/datapoint.



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>