

Technical Data Sheet

Caution: Research use only.

AlphaLISA® SureFire® Ultra™ p-ACC (Ser79) Assay Kit

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

Kit-Specificity Information

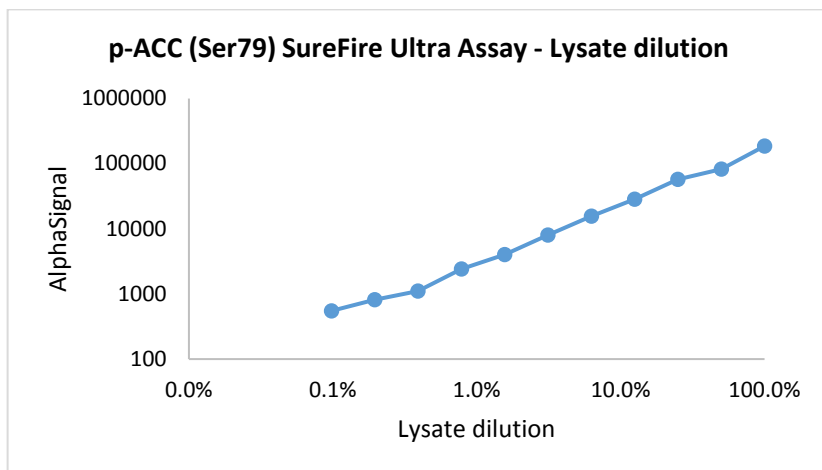
This assay kit contains antibodies which recognize the phospho-Ser79 epitope, and a distal invariant epitope, on p-ACC. The assay recognizes ACC corresponding to Entrez-Gene ID 31. These antibodies recognize ACC of human origin and is predicted to also recognize mouse and rat. Other species should be 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, then treated with 11mM H2O2 for 10min and lysed in 5ml 1X *SureFire Ultra* Lysis Buffer.

Representative data

Data obtained with a 2-plate, 2-incubation protocol.



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>.