

Technical Data Sheet

Caution: Research use only.

AlphaLISA® SureFire® Ultra™ p-Src (Tyr419) Assay Kit

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

Kit-Specificity Information

This assay kit contains antibodies that specifically recognize the phospho-Tyr419-epitope, or an invariant epitope on Src. The human protein detected by this kit corresponds to Entrez Gene ID 6714.

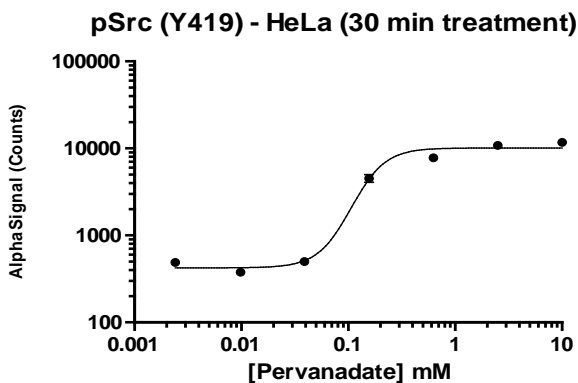
These antibodies recognize Src of human origin. Other species should be tested on a case-by-case basis.

Control Lysate Information

Positive Control Lysate: Prepared from HeLa cells, cultured to confluence in T175 flask in 10% FBS medium, then treated with 2.5% Pervanadate for 30 mins, and lysed in 4mL of 1X *SureFire Ultra* Lysis buffer.

Representative data

Data obtained with a 2-plate, 2-incubation protocol. HeLa cells were grown to confluence and stimulated with Pervanadate for 30 minutes.



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>