

Technical Data Sheet

Caution: Research use only.

AlphaLISA® SureFire® Ultra™ p-c-Met (Tyr1234/1235) Assay Kit

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

Kit-Specificity Information

This assay kit contains antibodies which recognize the phospho-Tyr1234/1235 epitope, and a distal invariant epitope, on c-Met. The assay recognizes c-Met corresponding to Entrez-Gene ID 4233.

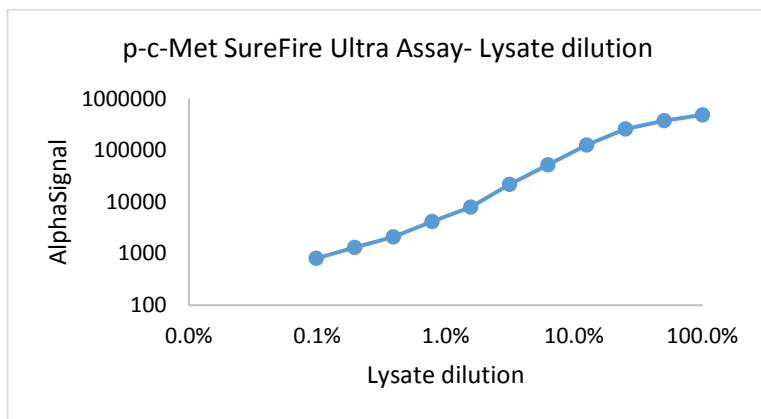
These antibodies recognize c-Met of human origin. Other species should be tested on a case-by-case basis.

Control Lysate Information

Positive Control Lysate: Prepared from HeLa cells, cultured to confluence in T175 flasks 10% FBS containing medium, then treated with 2% Pervanadate for 15min and lysed in 12ml 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>