

# Technical Data Sheet

## AlphaLISA® SureFire® Ultra™ Human p-c-Met (Tyr1234/1235) Detection Kit

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

### Kit-Specificity

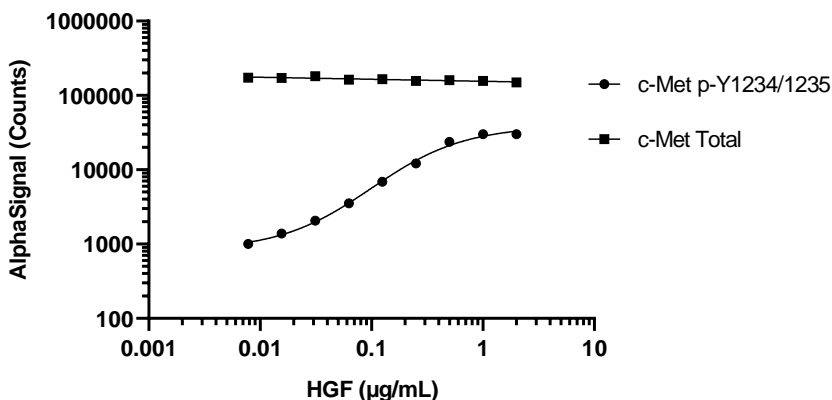
This assay kit contains antibodies which recognize the phospho-Tyr1234/1235 epitope and a distal epitope on c-Met. The protein detected by this kit corresponds to UniProt ID P08581. c-Met is also known as Hepatocyte growth factor receptor. These antibodies recognize c-Met of human origin. Other species 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 in 10% FBS containing medium, then treated with 1 mM Pervanadate for 15 minutes and lysed in 12 mL *SureFire Ultra* Lysis Buffer.

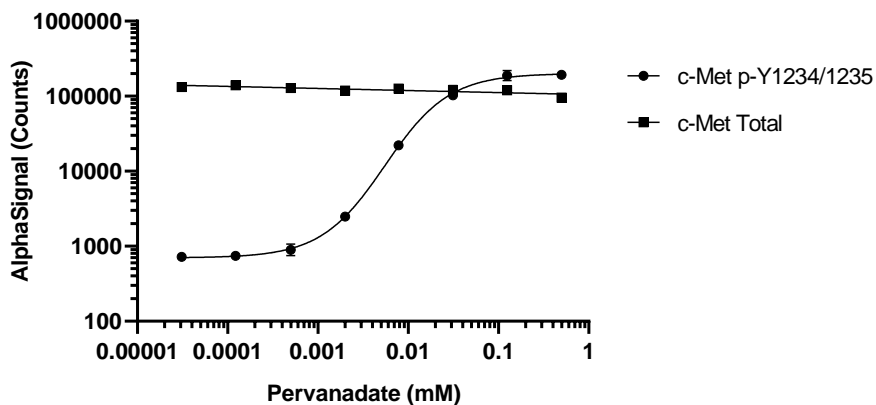
### Representative Data

Data obtained with a 2-plate, 2-incubation protocol A549 cells were plated at 40K cells/well in a 96 well plate and incubated overnight. Cells were treated with HGF at the indicated concentrations for 30 minutes. Cells were lysed with *SureFire Ultra* Lysis Buffer and assayed separately for Total and Phospho (Y1234/1235) c-Met, using respective *SureFire Ultra* kits. Equivalent to approximately 4,000 cells/datapoint.



## Technical Data Sheet

Data obtained with a 2-plate, 2-incubation protocol. A549 cells were plated at 40K cells/well in a 96 well plate and incubated overnight. Cells were treated with pervanadate at the indicated concentrations for 30 minutes. Cells were lysed with *SureFire Ultra* Lysis Buffer and assayed separately for Total and Phospho (Y1234/1235) c-Met, using respective *SureFire Ultra* kits. Equivalent to approximately 4,000 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>.