

Technical Data Sheet

AlphaLISA® SureFire® Ultra™ Human and Mouse Ret Total Detection Kit

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

Kit Specificity

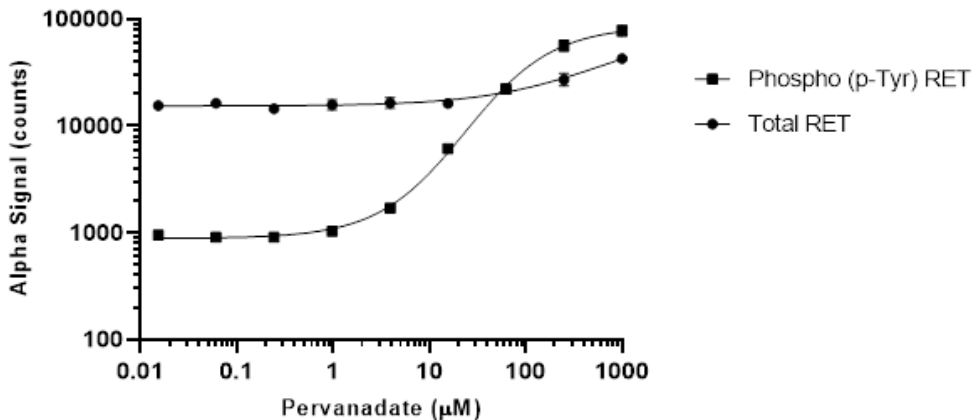
This assay kit contains antibodies which recognize invariant epitopes on Ret. The protein detected by this kit corresponds to UniProt ID P07949. Ret is also known as Cadherin family member 12 and Proto-oncogene c-Ret. These antibodies recognize Ret of human and mouse origin. Other species should be tested on a case-by-case basis.

Control Lysate Information

Positive Control Lysate: Prepared from SH-SY5Y cells, cultured to confluence in T175 flasks in 10% FBS containing medium, then treated with 12.5 μM retinoic acid for 16 hours and lysed in 5 mL of 1X *SureFire Ultra* Lysis Buffer.

Representative Data

Data obtained with a 2-plate, 2-incubation protocol. SH-SY5Y cells were plated at 40K cells/well for 6 h and treated with 12.5 μM retinoic acid for 16 h. Cells were then 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 (p-Tyr) Ret using respective *SureFire Ultra* detection kits. Equivalent to approximately 4000 cells/datapoint.



For more information and detailed protocols, please refer to the generic AlphaLISA *SureFire Ultra* Detection Kits Manual, available from <http://www.perkinelmer.com/category/alpha-surefire-kits>.