

Technical Data Sheet

AlphaLISA® *SureFire® Ultra™*

Human and Mouse Androgen Receptor Total Detection Kit

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

Kit Specificity

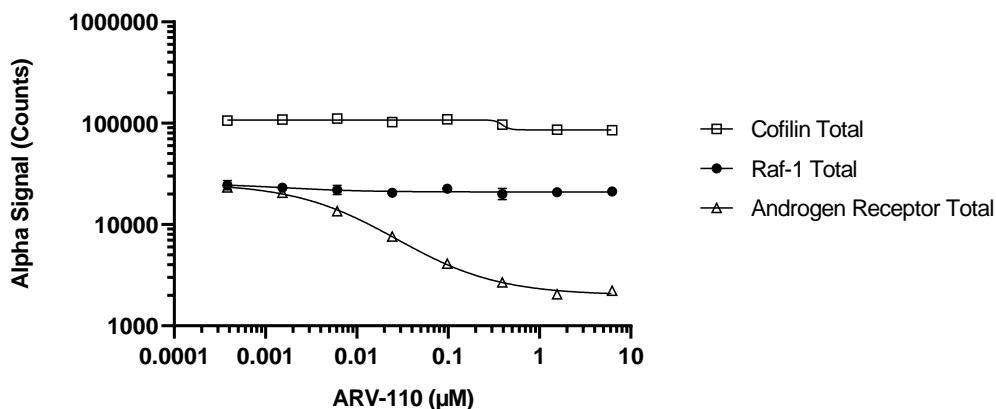
This assay kit contains antibodies which recognize invariant epitopes on Androgen Receptor. The protein detected by this kit corresponds to UniProt ID P10275. Androgen Receptor is also known as AR, NR3C4, AR8. These antibodies recognize Androgen Receptor of human and mouse origin. Other species should be tested on a case-by-case basis.

Control Lysate Information

Positive Control Lysate: Prepared from LNCaP cells, cultured to confluence in T175 flasks in 10% FBS containing medium and lysed in 4 mL of 1X *SureFire Ultra* Lysis Buffer.

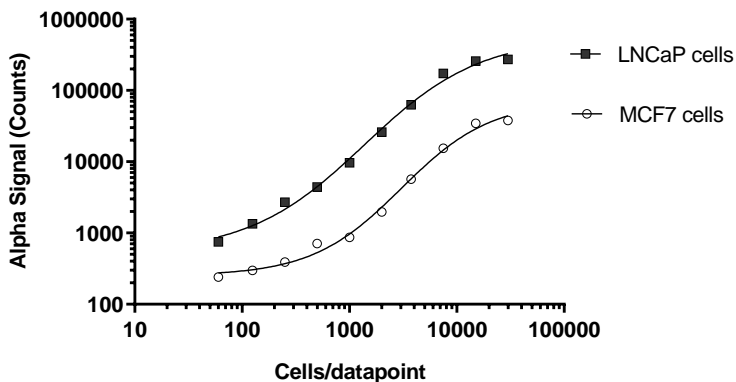
Representative Data

Data obtained with a 2-plate, 2-incubation protocol. MCF7 cells were plated overnight at 40K cells/well in a 96 well plate, then treated with the indicated concentrations of ARV-110 (Androgen Receptor PROTAC Degrader) for 24 h. Cells were lysed with *SureFire Ultra* Lysis Buffer and assayed separately for Androgen Receptor Total, Raf-1 Total and Cofilin Total using respective *SureFire Ultra* kits. Equivalent to approximately 4,000 cells/datapoint for Androgen Receptor Total and Raf-1 Total. Equivalent to approximately 400 cells/datapoint for Cofilin Total.

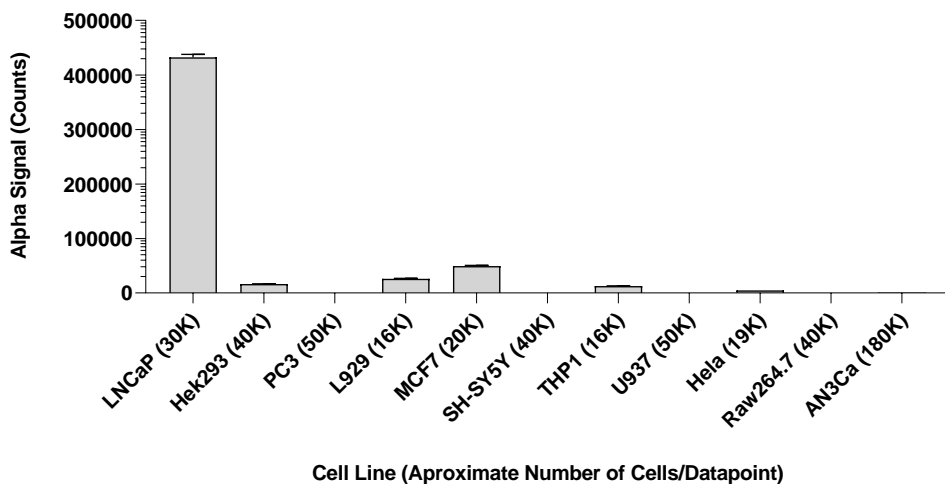


Technical Data Sheet

LNCaP and MCF7 cells were cultured to confluence in T175 flasks in 10% FBS containing medium and lysed in 4 mL of *SureFire Ultra* Lysis Buffer. Lysates were serially diluted in *SureFire Ultra* Lysis Buffer and assayed for Androgen Receptor Total. Approximate cells/datapoint are indicated.



Various lysates screened for Androgen Receptor Total. Lysates were prepared from cells lysed in *SureFire Ultra* Lysis Buffer and assayed for Androgen Receptor Total. Approximate cells/datapoint are indicated.



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>