

Technical Data Sheet

AlphaLISA® SureFire® Ultra™ IRAK4 Total Assay Kit

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

Kit Specificity

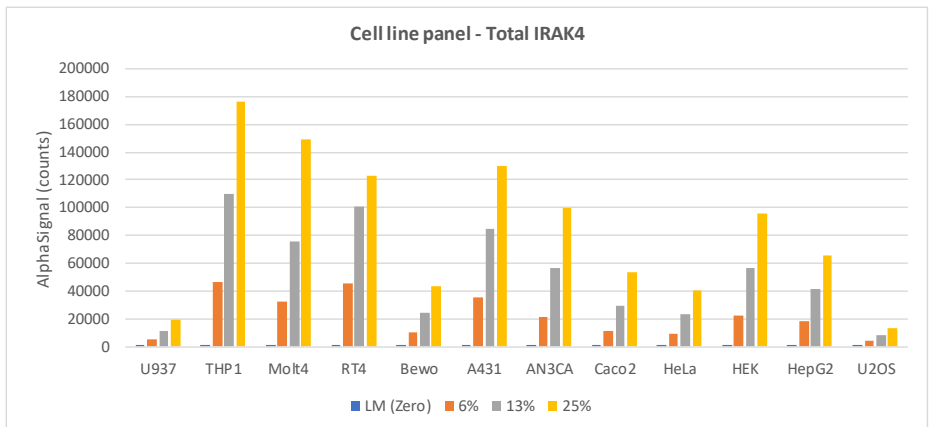
This assay kit contains antibodies which recognize two distinct invariant epitopes on IRAK4. The protein detected by this kit corresponds to UniProt ID Q9NWX3. IRAK4 is also known as Interleukin-1 receptor-associated kinase. These antibodies recognize human and mouse origin. Other species should be tested on a case-by-case basis.

Control Lysate Information

Positive Control Lysate: Prepared from THP-1 cells cultured in T175 flasks to 10^6 cells/mL in medium containing 10% FBS and lysed with 1x *SureFire Ultra* Lysis buffer at a final cell concentration of 2×10^6 cells/mL.

Representative Data

Data obtained with a 2-plate, 2-incubation protocol. A dilution series of various cell lysates from a panel of cell lines were prepared in 1X *SureFire Ultra* Lysis buffer at the approximate cell numbers indicated below. Lysates were assayed for Total IRAK4 by *SureFire Ultra*.



Cell Type	U937	THP1	Molt4	RT4	Bewo	A431	AN3CA	Caco2	HeLa	HEK	HEPG2	U2OS
Cells/datapoint at "25%"	3,000	8,000	20,000	9,000	2,750	7,500	13,750	2,250	8,000	10,000	7,000	5,000

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>.