

Technical Data Sheet

Caution: Research use only.

AlphaLISA® SureFire® Ultra™ p-Btk (Tyr223) Assay Kit

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

Kit-Specificity Information

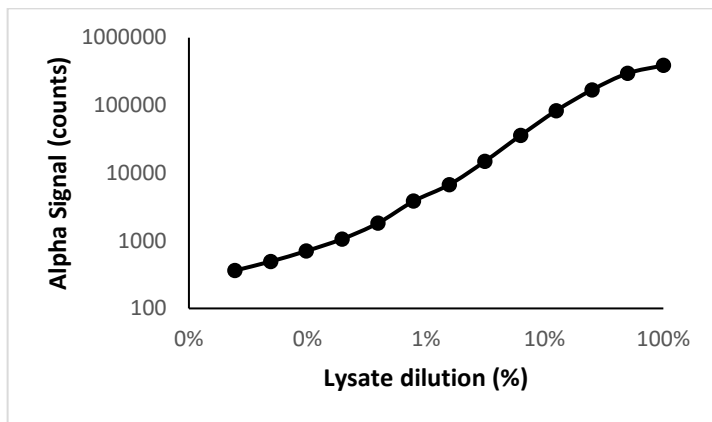
This assay kit contains antibodies which recognize the Tyr223 and an invariant epitope on Bruton's tyrosine kinase (Btk) . The protein detected by this kit corresponds to GenBank Accession NG_009616.1. These antibodies recognize Btk of human origin. Other species should be tested on a case-by-case basis.

Control Lysate Information

Positive Control Lysate: Prepared from U937 cells, cultured in a T175 flask to approximately 1×10^6 cells/mL in 10% FBS containing medium, then treated with 1% pervanadate for 30min and lysed with 5X *SureFire Ultra* Lysis buffer.

Representative data

Data obtained with a 2-plate, 2-incubation protocol. Control Lysate was diluted to the concentrations shown using 1X Lysis Buffer, and assayed for p-Btk (Tyr223) using the *SureFire Ultra* kit.



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>