

Technical Data Sheet

Caution: Research use only.

Alpha SureFire® Ultra™ Multiplex p-Btk (Tyr223) + Total Btk Assay Kit

**Products: MPSU-PTBTK-K500, MPSU-PTBTK-K10K, MPSU-PTBTK-K50K,
MPSU-PTBTK-K-HV**

Kit-Specificity Information

This assay kit contains antibodies which recognize the phospho-Tyr223 epitope and an invariant epitope on Bruton's tyrosine kinase (Btk), plus antibodies to measure the total levels of Btk. The proteins detected by this kit correspond to Entre Gene ID 695. These antibodies recognize BTK of human origin. Other species should be tested on a case-by-case basis.

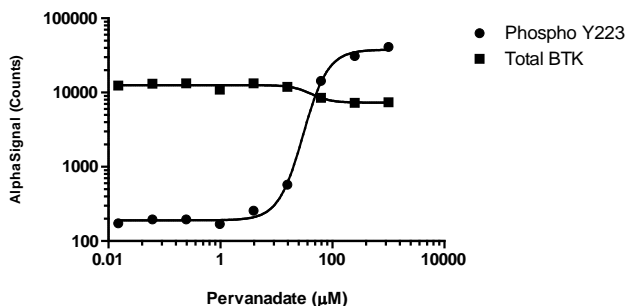
Control Lysate Information

Positive Control Lysate: Prepared from U937 cells, cultured in a T175 flask to approximately 1×10^6 cells/mL in 10% FBS containing medium, the 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. U937 cells were plated at 80K cells per well, and treated with Pervanadate at the indicated concentrations for 30min, then lysed with 5X Lysis buffer. Samples were assayed for both phospho and total BTK as detailed in the MPSU manual.

Pervanadate DR in U937 cells - Total-Phospho BTK Multiplex



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>