

# Technical Data Sheet

Caution: Research use only.

## AlphaLISA® SureFire® Ultra™ p-Btk (Tyr551) Assay Kit

**Products: ALSU-PBTK-B500, ALSU-PBTK-B10K, ALSU-PBTK-B50K,  
ALSU-PBTK-B-HV**

### **Kit-Specificity Information**

This assay kit contains antibodies which recognize the Tyr551 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, and are predicted to recognise the rat and mouse proteins. Other species should be tested on a case-by-case basis.

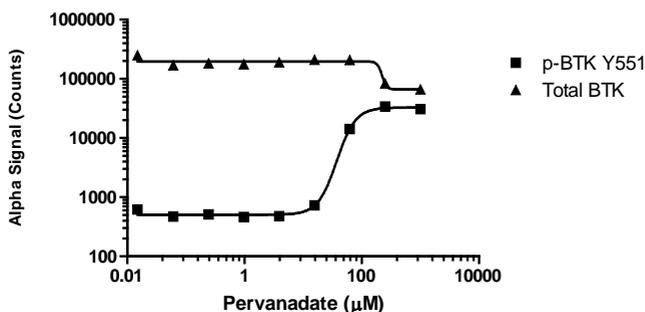
### **Control Lysate Information**

Positive Control Lysate: Prepared from U937 cells, cultured in a T175 flask to approximately  $1 \times 10^6$  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. U937 cells were plated at 200K cells per well, and treated with Pervanadate at the indicated concentrations for 30 minutes, then lysed in medium with 5X Lysis buffer. Levels of Btk (Tyr551) phosphorylation, and Total Btk were assayed with respective *SureFire Ultra* assay kits.

#### **Pervanadate Dose Response in U937 cells - BTK assays**



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>