

Alpha SureFire® Ultra™ Multiplex

Human p-Btk (Tyr223) + Total Btk Detection Kit

Product number: MPSU-PTBTK-K500, MPSU-PTBTK-K10K,
MPSU-PTBTK-K50K, MPSU-PTBTK-K-HV



Kit specificity:

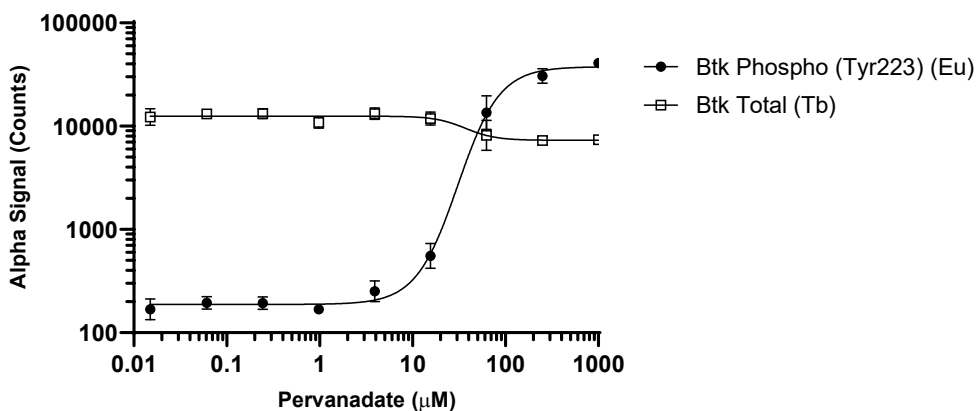
This assay kit contains antibodies which recognize the phospho-Tyr223 epitope and a distal epitope on BTK, plus antibodies to detect Total Btk. The proteins detected by this kit correspond to UniProt ID Q06187. These antibodies recognize BTK of human origin. Other species should be tested on a case-by-case basis.

Control lysate information:

Positive Control Lysate: Prepared from U-937 cells cultured in 10% FBS containing medium. Cells were harvested and washed with HBSS + 0.1% BSA, adjusted to 2×10^6 cells/mL and treated with 0.5 mM Pervanadate for 30 minutes. Cells were spun down at 300 RCF for 5 minutes and lysed with Lysis Buffer at 1×10^6 cells/mL.

Representative data

Data obtained with a 2-plate, 2-incubation protocol. U937 cells were harvested and washed in HBSS + 0.1% BSA and seeded at 200K cells/well in a 96 well plate. Cells were treated with Pervanadate at the indicated concentrations for 30 minutes. Cells were lysed with 5X Lysis Buffer and assayed for Phospho (Tyr223) and Total BTK in the same wells of an assay plate using the SureFire Ultra Multiplex kit. Equivalent to approximately 8,000 cells/datapoint.



The information provided in this document is for reference purposes only and may not be all-inclusive. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume liability for the accuracy or completeness of the information contained herein. Users should exercise caution when handling materials as they may present unknown hazards. Revvity shall not be liable for any damages or losses resulting from handling or contact with the product, as Revvity cannot control actual methods, volumes, or conditions of use. Users are responsible for ensuring the product's suitability for their specific application. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF"