

AlphaLISA® SureFire® Ultra™

Human p-TrkA (Tyr674/675) / p-TrkB (Tyr706/707) Detection Kit

Product number: ALSU-PTRKAB-A500, ALSU-PTRKAB-A10K,
ALSU-PTRKAB-A50K, ALSU-PTRKAB-A-HV



Kit specificity:

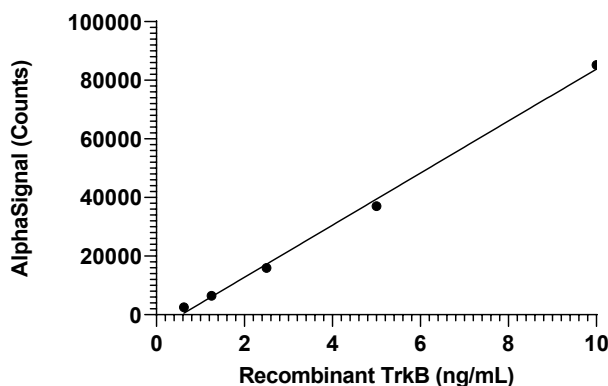
This assay kit contains antibodies which recognize the phospho-Tyr674/675 (TrkA) and/or phospho-Tyr706/707 (TrkB) epitopes, and a distal epitope common to TrkA and TrkB. The TrkA protein detected by this kit corresponds to UniProt ID P04629 and TrkB protein corresponds to UniProt ID Q16620. TrkA is also known as High affinity nerve growth factor receptor and Tropomyosin-related kinase A. TrkB is also known as BDNF/NT-3 growth factor receptor and Tropomyosin-related kinase B. These antibodies recognize p-Trk A and B of human origin. Other species should be tested on a case-by-case basis.

Control lysate information:

Positive Control Lysate: Prepared from purified, kinase active, recombinant TrkB protein, which has been autophosphorylated in vitro by incubation with ATP-Mg²⁺. Prepared at a concentration of 10 ng/mL.

Representative data

Data obtained with a 2-plate, 2-incubation protocol. Control lysate was serially diluted in Lysis Buffer and assayed for Phospho TrkA (Tyr674/675)/TrkB (Tyr706/707).



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"