

Alpha SureFire® Ultra™ Multiplex

Human and Mouse p-SMAD2 (Ser465/467) + Total SMAD2 Detection Kit

Product number: MPSU-PTSM2-K500, MPSU-PTSM2-K10K,
MPSU-PTSM2-K50K, MPSU-PTSM2-K-HV



Kit specificity:

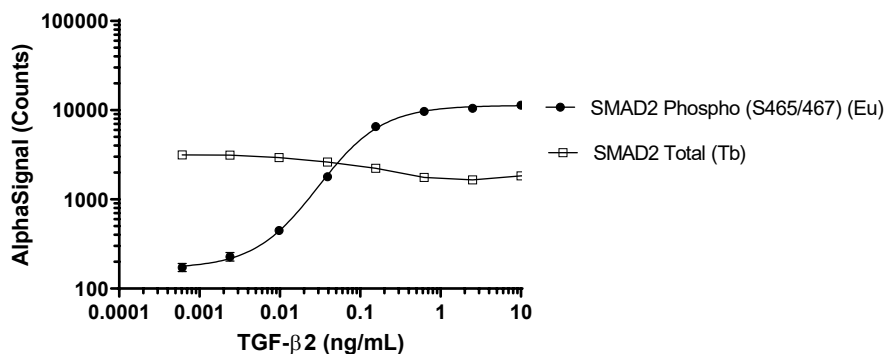
This assay kit contains antibodies which recognize the phospho-Ser465/467 epitope and a distal epitope on SMAD2, plus antibodies to detect Total SMAD2. The proteins detected by this kit correspond to UniProt ID Q15796. This kit will not cross react with SMAD3. These antibodies recognize SMAD2 of human and mouse origin. Other species should be tested on a case-by-case basis.

Control lysate information:

Positive Control Lysate: Prepared from RAW 264.7 cells cultured to confluence in T175 flasks in 10% FBS containing medium, treated with 50 ng/mL rhTGFβ for 90 minutes and lysed with 6 mL of Lysis buffer.

Representative data

Data obtained with a 2-plate, 2-incubation protocol. RAW 264.7 cells were seeded at 40K cells/well in a 96 well plate and incubated for 48 hours. Cells were serum starved for 1 hour and then treated with rhTGFβ at the indicated concentrations for 90 minutes. Cells were with lysed with Lysis Buffer and assayed for Phospho (Ser465/467) and Total SMAD2 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"