

Technical Data Sheet

Caution: Research use only.

AlphaLISA® SureFire® Ultra™ SMAD2 Total Assay Kit

**Products: ALSU-TSM2-A500, ALSU-TSM2-A10K, ALSU-TSM2-A50K,
ALSU-TSM2-A-HV**

Kit-Specificity Information

This assay kit contains antibodies which recognize two invariant epitopes, on SMAD2. The protein detected by this kit corresponds to Entre-Gene Id 4087. This kit will not cross react with SMAD3. These antibodies recognize SMAD2 of human, rat and mouse origin. Other species should be tested on a case-by-case basis.

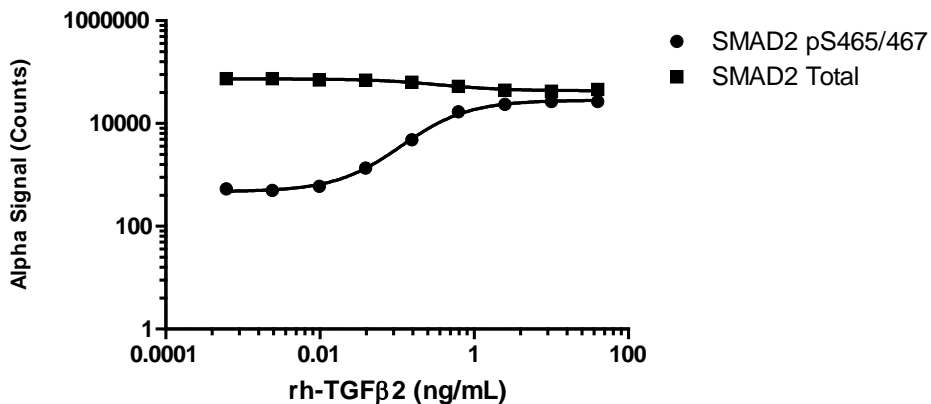
Control Lysate Information

Positive Control Lysate: Prepared from RAW 264.7 cells, cultured to confluence in T175 flasks in 10% FBS containing medium, then treated with 10ng/mL rhTGFβ for 90min and lysed in 6mL of 1X *SureFire Ultra* Lysis buffer.

Representative data

Data obtained with a 2-plate, 2-incubation protocol. Phosho and Total SMAD2 were measured in separate wells using respective *SureFire Ultra* assay kits.

rh-TGFβ2 Dose Reponse on RAW264.7 Cells



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>