

# Technical Data Sheet

Caution: Research use only.

## AlphaLISA® SureFire® Ultra™ p-SMAD1 (Ser463/465) Assay Kit

**Products: ALSU-PSM1-A500, ALSU-PSM1-A10K, ALSU-PSM1-A50K,  
ALSU-PSM1-A-HV**

### Kit-Specificity Information

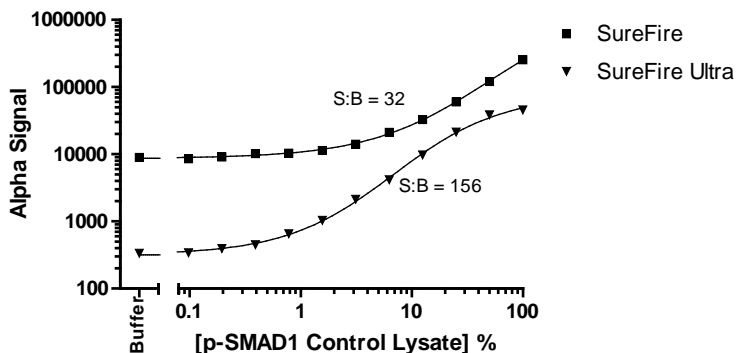
This assay kit contains antibodies which recognize the phospho-Ser463/465 epitope, and a distal epitope, on SMAD family member 1 (SMAD1). The protein detected by this kit corresponds to GenBank Accession NP\_001003688. Based on sequence similarity, cross reaction to SMAD5 and SMAD8 may occur. SMAD1 is also known as BSP1, JV41, JV4-1, MADH1 and MADR1. These antibodies recognize SMAD1 of human, mouse and rat origin. Other species should be tested on a case-by-case basis.

### Control Lysate Information

Positive Control Lysate: Prepared from C2C12 cells, cultured to confluence in T175 flasks in 10% FBS containing medium, then treated with 50ng/mL rhBMP4 for 30min and lysed in 12mL of 1X *SureFire Ultra* Lysis buffer.

### Representative data

Data obtained with a 2-plate, 2-incubation protocol. Data shows a comparison of AlphaScreen *SureFire* and AlphaLISA *SureFire Ultra* assay performance. Control lysate was diluted in the Lysis buffer specific for the assay type.



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>