

Technical Data Sheet

Caution: Research use only.

Alpha SureFire® Ultra™ Multiplex

Terbium SureFire Ultra p-SMAD1 (Ser463/465) Assay Kit

For multiplexing with AlphaLISA SureFire Ultra Assay Kits

**Products: TBSU-PSM1-A500, TBSU-PSM1-A10K, TBSU-PSM1-A50K,
TBSU-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.

The antibodies in this kit are identical to those in the corresponding ALSU-PSM1-A kits, other than the CaptSure2 tag.

Control Lysate Information

Positive Control Lysate: Prepared from C2C12 cells, cultured to confluence in T175 flasks in 10% FBS containing medium for 3 days, then treated with rhBMP4 (50ng/mL) for 30min and lysed in 13mL of 1X SureFire Ultra Lysis buffer.

Representative data

Control lysate samples measured in the terbium channel using an Envision plate reader.

