

# Technical Data Sheet

Caution: Research use only.

## Alpha SureFire® Ultra™ Multiplex

### Terbium SureFire Ultra p-SMAD3 (Ser423/425) Assay Kit

For multiplexing with AlphaLISA SureFire Ultra Assay Kits

**Products: TBSU-PSM3-A500, TBSU-PSM3-A10K, TBSU-PSM3-A50K,  
TBSU-PSM3-A-HV**

#### **Kit-Specificity Information**

This assay kit contains antibodies which recognize the phospho-Ser423/425 epitope, and a distal epitope, on SMAD3. The protein detected by this kit corresponds to GenBank Accession NP\_001138574. Based on sequence identity, cross reaction to SMAD2 is predicted to be likely. SMAD3 is also known as MADH3, JV15-2, HSPC193, HsT17436, MGC60396, DKFZp586N0721 and DKFZp686J10186. These antibodies recognize SMAD3 of human and mouse 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-PSM3-A kits, other than the CaptSure2 tag.

#### **Control Lysate Information**

Positive Control Lysate: Prepared from RAW 264.7 cells, cultured to confluence in T175 flasks in 10% FBS containing medium for 3 days, then treated with rhTGFβ (10ng/mL) for 90min and lysed in 6mL of 1X SureFire Ultra Lysis buffer.

#### **Representative data**

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

