

# Technical Data Sheet

Caution: Research use only.

## Alpha SureFire® Ultra™ Multiplex

### Terbium SureFire Ultra p-4E-BP1 (Thr37/46) Assay Kit

For multiplexing with AlphaLISA SureFire Ultra Assay Kits

**Products: TBSU-P4EBP-A500, TBSU-P4EBP-A10K, TBSU-P4EBP-A50K,  
TBSU-P4EBP-A-HV**

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

This assay kit contains antibodies which recognize the phospho-Thr37/46 epitope, and a distal epitope, on 4E-BP1. The protein detected by this kit corresponds to GenBank Accession NP\_004086. 4E-BP1 is also known as BP-1, 4EBP1; PHAS-I, MGC4316 and EIF4EBP1. These antibodies recognize 4E-BP1 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-P4EBP-A kits, other than the CaptSure2 tag.

#### **Control Lysate Information**

Positive Control Lysate: Prepared from MCF7 cells, cultured to confluence in T175 flasks in 10% FBS containing medium for 3 days, then treated with insulin (500ng/mL) for 30min and lysed in 30mL of 1X SureFire Ultra Lysis buffer.

#### **Representative data**

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

