

Technical Data Sheet

Caution: Research use only.

Alpha SureFire® Ultra™ Multiplex

Terbium SureFire Ultra p-AKT (Thr308) Assay Kit For multiplexing with AlphaLISA SureFire Ultra Assay Kits

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

Kit-Specificity Information

This assay kit contains antibodies which recognize the phospho-Thr308 epitope, and a distal epitope, common to Akt 1, Akt 2 and Akt 3. The proteins detected by this kit correspond to GenBank Accessions NP_001014431, NP_001617, and NP_859029. Akt proteins are also known as PKB or RAC. These antibodies recognize Akt 1/2/3 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-PAKT-A kits, other than the CaptSure2 tag.

Control Lysate Information

Positive Control Lysate: Prepared from HEK293 cells, cultured to confluence in T175 flasks in 10% FBS containing medium for 3 days, then treated with 20% FBS for 15min and lysed in 10mL of 1X SureFire Ultra Lysis buffer.

Representative data

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

