

Technical Data Sheet

Caution: Research use only.

Alpha SureFire® Ultra™ Multiplex

Terbium SureFire Ultra p-p38 MAPK (Thr180/Tyr182) Assay Kit

For multiplexing with AlphaLISA SureFire Ultra Assay Kits

**Products: TBSU-PP38-B500, TBSU-PP38-B10K, TBSU-PP38-B50K,
TBSU-PP38-B-HV**

Kit-Specificity Information

This assay kit contains antibodies which recognize the phospho-Thr180/Tyr182 epitope, and a distal epitope, on p38 MAPK. The protein detected by this kit corresponds to GenBank Accession NP_001306. Also known as p38; SAPK2A; p38alpha; MAPK14. These antibodies recognize p38 MAPK of human, mouse, 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-PP38-B kits, other than the CaptSure2 tag.

Control Lysate Information

Positive Control Lysate: Prepared from HeLa cells, cultured to confluence in T175 flasks in 10% FBS containing medium for 3 days, then treated with anisomycin (2 µg/mL) for 15 min and lysed in 8 mL of 1X SureFire Ultra Lysis buffer.

Representative data

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

