

# Technical Data Sheet

Caution: Research use only.

## AlphaLISA® SureFire® Ultra™ MKK4 Total Assay Kit

**Products: ALSU-TMKK4-A500, ALSU-TMKK4-A10K, ALSU-TMKK4-A50K,  
ALSU-TMKK4-A-HV**

### **Kit-Specificity Information**

This assay kit contains antibodies which recognize two invariant epitopes on mitogen-activated protein kinase kinase 4 (MKK4). The protein detected by this kit correspond to GenBank Accession NP\_003001. MKK4 is also known as JNKK, MEK4, SEK1, JNKK1, SERK1, MAPKK4, PRKMK4 and MAP2K4.

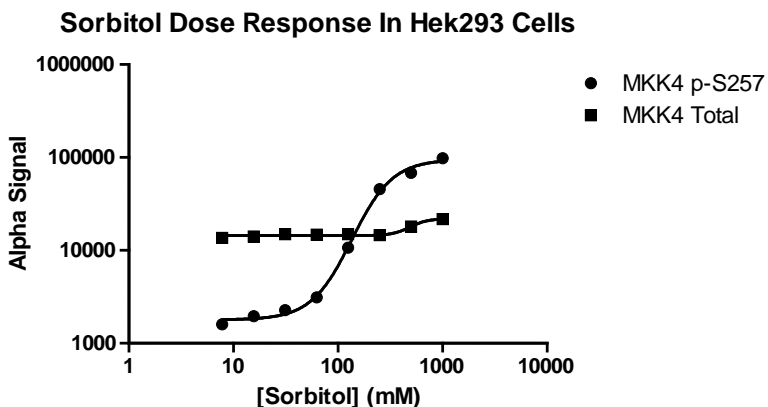
These antibodies recognize MKK4 of human origin. Other species should be tested on a case-by-case basis.

### **Control Lysate Information**

Positive Control Lysate: Prepared from HEK293 cells, cultured to confluence in T175 flasks in 10% FBS containing medium, then treated with 1M Sorbitol for 45min and lysed in 4mL of 1X *SureFire Ultra* Lysis buffer.

### **Representative data**

Data obtained with a 2-plate, 2-incubation protocol. HEK293 cells were grown to confluence and stimulated with Sorbitol for 45 minutes.



For more information and detailed protocols, please refer to the generic AlphaLISA *SureFire Ultra* Assay Kits Manual, available from <http://www.perkinelmer.com/category/alpha-surefire-kits>