

Technical Data Sheet

Caution: Research use only.

AlphaLISA® SureFire® Ultra™ eIF4E Total Assay Kit

**Products: ALSU-TEIF4-A500, ALSU-TEIF4-A10K, ALSU-TEIF4-A50K,
ALSU-TEIF4-A-HV**

Kit-Specificity Information

This assay kit contains antibodies which recognize invariant epitopes on eukaryotic translation initiation factor 4E (eIF4E). The protein detected by this kit corresponds to Entrez Gene ID 1977. eIF4E is also known as CBP, EIF4F, EIF4E1, EIF4EL1 and MGC111573. These antibodies recognize eIF4E of human origin. Other species should be tested on a case-by-case basis.

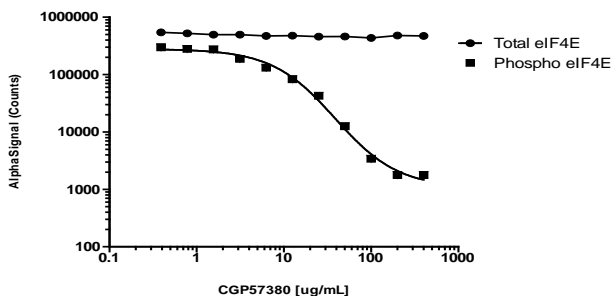
Control Lysate Information

Positive Control Lysate: Prepared from HEK293 cells, cultured to confluence in T175 flasks in 10% FBS containing medium, then treated with 200 ng/mL EGF for 30 min and lysed in 4 mL of 1X *SureFire Ultra* Lysis buffer.

Representative data

Data obtained with a 2-plate, 2-incubation protocol. HEK293 cells were plated overnight at 40 K cells/well, then inhibited with the indicated concentrations of CGP57380 for 2 hrs. Treated with 200 ng/mL of EGF for 30 min, then lysed with 100 µL of 1X Lysis Buffer. Samples were assayed separately for total and phospho (S209) eIF4E, using respective kits.

eIF4E Inhibition Dose Response - HEK293 cells



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>