

# Technical Data Sheet

## AlphaLISA® SureFire® Ultra™ β-Catenin Total Assay Kit

ALSU-TBCAT-A500, ALSU-TBCAT-A10K, ALSU-TBCAT-A50K, ALSU-TBCAT-A-HV

### Kit Specificity

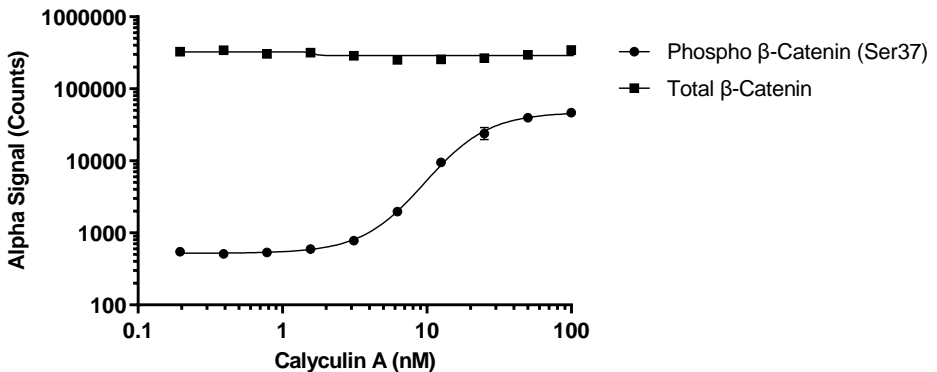
This assay kit contains antibodies which recognize invariant epitopes on β-Catenin. The protein detected by this kit corresponds to UniProt ID P35222. β-Catenin is also known as CTNNB1 and Catenin Beta 1. These antibodies recognize β-Catenin of human and mouse origin. Other species should be tested on a case-by-case basis.

### Control Lysate Information

Positive Control Lysate: Prepared from A549 cells, cultured to confluence in T175 flasks in 10% FBS containing medium. Cells were treated with 1 μg/mL PMA and 50 nM Calyculin A for 4 hours at 37°C. Cells were lysed in 5 mL of 1X *SureFire Ultra* Lysis Buffer.

### Representative Data

Data obtained with a 2-plate, 2-incubation protocol. A549 cells were seeded at 40K cells per well in a 96 well plate and incubated overnight. Cells were treated with 1 μg/mL PMA and Calyculin A at the indicated concentrations for 4 hours at 37°C. Cells were lysed with 1X *SureFire Ultra* Lysis Buffer, and assayed for Total or Phospho (Ser37) β-Catenin in separate wells of an assay plate with respective *SureFire Ultra* kits. Equivalent to approximately 4,000 cells/datapoint.



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>