[pThr¹⁸⁰/Tyr¹⁸²]p38 ELISA kit

The [pThr¹⁸⁰/Tyr¹⁸²]p38 ELISA kit is a colorimetric immunometric enzyme immunoassay kit

The [pThr¹⁸⁰/Tyr¹⁸²]p38 EIA kit is a colorimetric immunometric enzyme immunoassay kit with results in 3 hours. Absorbance is read at 450 nm. Use less precious sample per test compared to Western blot with higher throughput sample analysis; measure up to 40 samples in duplicate per kit.

Six proteins of the p38 subfamily of MAP kinases have been identified to date, including: p38α (SAPK2, CSBP) and its splice variant Mxi; p38β and β2; p38γ (ERK6, SAPK3); and p38δ (SAPK4). p38 MAP kinases contain a conserved TGY motif in their activation loop and are activated in response to cytokines, hormones, GPCRs, and osmotic and heat shock to mediate primarily pro-apoptotic and pro-inflammatory responses. MKK3, MKK4, and MKK6 are known to activate p38 family members with varying specificity. All four p38 isoforms are able to phosphorylate the activating transcription factor 2 (ATF2), and differentially activate numerous substrates including Mnk1/2, Msk, phospholipase A2, and the MAPKAP kinase-2 and -3. Activated MAPKAP kinase-2 subsequently phosphorylates the heat shock protein Hsp27, an F-actin binding protein involved in cytoskeletal maintenance.

Citations: 3

View Online »

Ordering Information

Order Online »

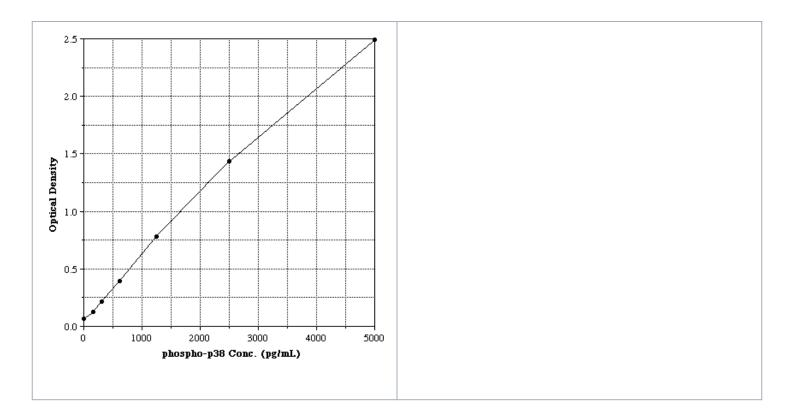
ADI-900-101

96 wells

Manuals, SDS & CofA

View Online »

- Sensitive use less precious sample per test compared to Western blot
- Higher throughput test up to 40 samples in duplicate in just 3 hours
- Quantitative obtain statistically signficant results and compare data between multiple experiments
- Ready-to-use save time and reduce errors with pre-coated plates



Handling & Storage

Use/Stability Store all components at 4°C, except standard at -20°C.

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name SAPK2, SAP kinase 2

Application Colorimetric detection, ELISA

Application Notes For the quantitative determination of [pThr¹⁸⁰/Tyr¹⁸²]p38

in cell lysates of human and mouse origin.

Assay Time 3 hours

Compatibility This product is compatible with the Absorbance 96 Plate

Reader.

Contents Microtiter plate, Conjugate, Antibody, Assay buffer 4

concentrate, Wash buffer concentrate, Standard, TMB

Substrate, RIPA Cell lysis buffer 2, Stop solution 2

Crossreactivity Phospho p38 (100%) and <0.01%: non-phosphorylated

p38, phospho Jnk, non-phosphorylated Jnk, AKT, non-

phosphorylated Erk, phospho Erk

Sensitivity 52.1 pg/ml (range 156 – 5,000 pg/ml)

Species Reactivity Human, Mouse

UniProt ID Q16539 (human p38α), Q15759 (human p38β), P53778

 $(p38\gamma)$, O15264 $(p38\delta)$

Wavelength 450 nm

