[pThr¹⁸³/Tyr¹⁸⁵]Jnk1/2 ELISA kit

The [pThr¹⁸³/Tyr¹⁸⁵]Jnk1/2 ELISA kit is a colorimetric immunometric enzyme immunoassay kit

The [pThr¹⁸³/Tyr¹⁸⁵]Jnk1/2 EIA kit is a colorimetric immunometric enzyme immunoassay kit with results in 3 hours. Absorbance is read at 450 nm. Save time, money, and precious sample with fully quantitative results and increased sensitivity compared to Western blot analysis.

The Stress-activated protein kinases alpha and γ (SAPK α and γ), also know as c-Jun NH2-terminal kinases 1 and 2 (Jnk1/2) are serine/ threonine protein kinase belonging to the MAP kinase family. The extracellular stimuli leading to cell stress include inflammatory cytokines (via Rho and Rac), growth factors, protein synthesis inhibitors and osmotic stress amongst others. Both Jnk1 and Jnk2 must be dually phosphorylated in the activation loop, at threonine 183 and tyrosine 185, to achieve full enzymatic activity.

Citations: 1

View Online »

Ordering Information

Order Online »

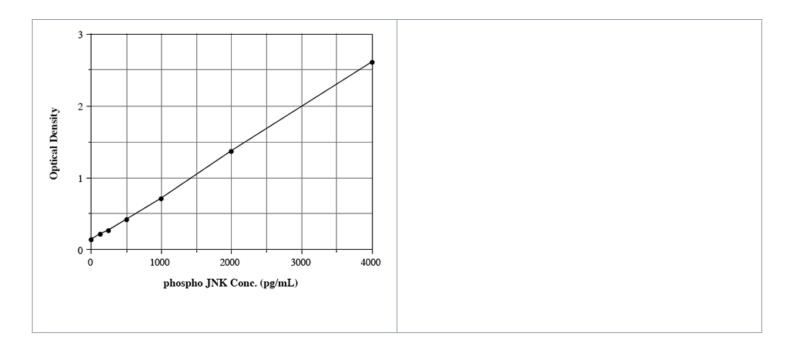
ADI-900-106

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 Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name c-Jun N-terminal kinase 1/2, SAP kinase 1a/c, MAP kinase

8/9, MAPK8/9

Application Colorimetric detection, ELISA

Application Notes For the quantitative determination of [pThr¹⁸³/Tyr¹⁸⁵

]Jnk1/2 in cell lysates of human, mouse, and rat 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 Jnk1 (100%), phospho Jnk2 (98.6%), phospho

p38 (0.42%), phospho Erk2 (0.16%) and <0.01%: non-

phosphorylated Erk2, non-phosphorylated Jnk1

Sensitivity 75.8 pg/ml (range 125 – 4,000 pg/ml)

Species Reactivity Human, Mouse, Rat

UniProt ID P45983 (Jnk1), P45984 (Jnk2)

Wavelength 450 nm

