

# [pThr<sup>180</sup>/Tyr<sup>182</sup>]p38

## ELISA kit

The [pThr<sup>180</sup>/Tyr<sup>182</sup>]p38 ELISA kit is a colorimetric immunometric enzyme immunoassay kit

The [pThr<sup>180</sup>/Tyr<sup>182</sup>]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: p38 $\alpha$  (SAPK2, CSBP) and its splice variant Mxi; p38 $\beta$  and  $\beta$ 2; p38 $\gamma$  (ERK6, SAPK3); and p38 $\delta$  (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

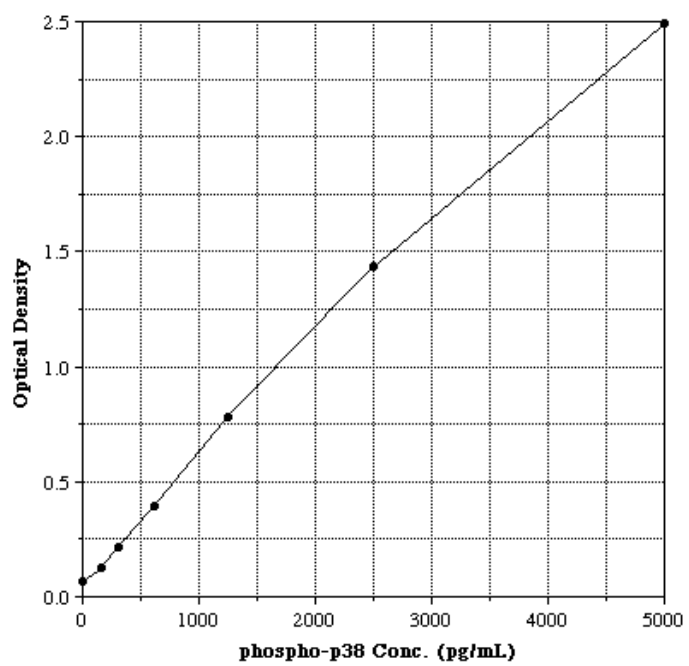
[Order Online »](#)

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 significant 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 [pThr<sup>180</sup>/Tyr<sup>182</sup>]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 p38 $\alpha$ ), Q15759 (human p38 $\beta$ ), P53778 (p38 $\gamma$ ), O15264 (p38 $\delta$ )

**Wavelength** 450 nm

Last modified: May 29, 2024



Enzo Life Sciences, INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)