

[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: p38 α (SAPK2, CSBP) and its splice variant Mxi; p38 β and β 2; p38 γ (ERK6, SAPK3); and p38 δ (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

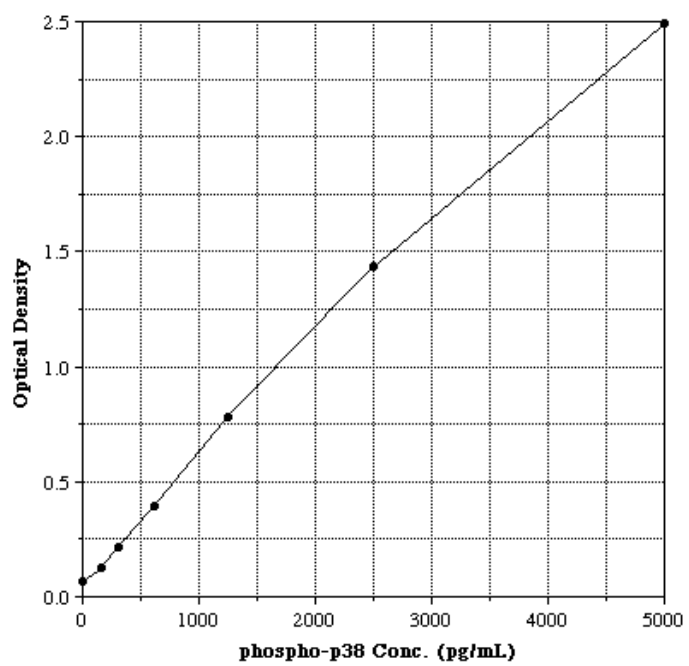
[Order Online »](#)

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 [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 p38 α), Q15759 (human p38 β), P53778 (p38 γ), O15264 (p38 δ)

Wavelength 450 nm



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-be@enzolifesciences.com)

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