

# Phosphothreonine detection set

This kit offers a set of three different monoclonal antibodies specific for threonine phosphorylation sites. Phosphorylation patterns in a given cell extract may differ when probed with different antibodies due to sequence specificity. Recognition therefore depends on two criteria: Phosphorylation and surrounding aa motif. If one of the two criteria is not fulfilled, the antibody will not detect the phosphorylation site.

## Ordering Information

[Order Online »](#)

ALX-850-024-KI01	1Set
------------------	------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Long Term Storage	-20°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Application	IP, WB
Contents	Contains a positive control for immunoblot applications as well as 25µg each of the following MAbs: 1E11 (Prod. No. ALX-804-168) 4D11 (Prod. No. ALX-804-169) 14B3 (Prod. No. ALX-804-170)
Recommendation Dilutions/Conditions	Western Blot (20µl/lane (mini gel) for HRPO/DAB detection, or 5µl/lane (mini gel) for HRPO/ECL detection. Do not use milk/casein for blocking and dilution! Replace by BSA! We recommend to use 1% BSA, 1% PVP-10 (polyvinylpyrrolidone) 1% PEG 3500, 0.2% Tween 20 in 2x PBS) Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

## Technical Info / Product Notes

Antibodies directed against phosphorylated epitopes recognize the phosphorylated amino acid in the context of the surrounding amino acid sequence. Recognition is therefore dependent on 2 criteria: 1) phosphorylation and 2) the surrounding amino acid motif. If one of the two criteria is not fulfilled, the antibody will not detect the phosphorylation site. Since the amino acid sequence varies between different phosphorylation sites, certain proteins – though phosphorylated – may not be detected by the antibody. Phosphorylation patterns in a given cell extract may differ when probed with different antibodies due to sequence specificity.

### Positive Control Cell Lysate

Formulation: The pSer/pThr molecular weight marker contains rabbit muscle phosphoproteins isolated by Fe<sup>3+</sup>/IDA – affinity chromatography. Proteins are lyophilized from PBS/NaF/PEG/Sucrose and Bromophenolblue.

Reconstitution: Reconstitute by addition of 200 µl H<sub>2</sub>O. After complete solubilization add 200 µl 2x SDS-PAGE sample buffer, mix and incubate at 90°C for 5 min. Aliquote and store frozen. Avoid repeated freeze/thaw cycles.

Application: The pSer/pThr molecular weight marker is recommended for immunoblot applications. Use 20µl molecular weight marker per lane. Note: Use BSA-based blot incubation buffers. Milk, Casein and Blotto might interfere with antibody – antigen interaction.



ENZO LIFE SCIENCES INC. Phone: 800.942.0430 <a href="mailto:info-usa@enzolifesciences.com">info-usa@enzolifesciences.com</a>	ENZO LIFE SCIENCES & Luxembourg Phone: +41 61 926 8989 <a href="mailto:info-eu@enzolifesciences.com">info-eu@enzolifesciences.com</a>	ENZO LIFE SCIENCES France Phone: +33 472 440 655 <a href="mailto:info-fr@enzolifesciences.com">info-fr@enzolifesciences.com</a>	ENZO LIFE SCIENCES Germany Phone: +49 7621 5500 526 <a href="mailto:info-de@enzolifesciences.com">info-de@enzolifesciences.com</a>	ENZO LIFE SCIENCES Phone (UK customers): +44 1392 825900 <a href="mailto:info-uk@enzolifesciences.com">info-uk@enzolifesciences.com</a>
---	--	--	---	--