

# [pSer383]Elk1

## polyclonal antibody

This antibody is covered by our [Worry-Free Guarantee](#).

### Ordering Information

[Order Online »](#)

BML-SA381-0020	20µl
----------------	------

### Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage      -20°C

Shipping      Dry Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	ETS domain-containing protein Elk-1
Application	WB
Formulation	Liquid. In HEPES buffer containing BSA and glycerol.
Host	Rabbit
Immunogen	Synthetic peptide corresponding to a portion of human Elk1 phosphorylated at Ser383.
Species Reactivity	Human
Specificity	Recognizes Elk1 phosphorylated at Ser <sup>383</sup> .
UniProt ID	P19419
Worry-free Guarantee	This antibody is covered by our <a href="#">Worry-Free Guarantee</a> .

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)