

# p53 activator

## p53 activator

A peptide corresponding to 'peptide 46', (aa 361-382), which corresponds to the single-stranded DNA-end binding site of p53. This sequence is an activator of wild-type and certain p53 mutants. Fusion Peptide 46, a 35-mer peptide, consists of peptide 46 linked to a sequence from the third helix of *Drosophila* Antennapedia homeodomain protein in order to confer cell permeability. The p53 Activator (Fusion Peptide 46) has been demonstrated to have numerous effects including: rescue of p53 growth inhibition (5 µg/ml) in a human osteosarcoma Saos-2 subline transfected with His<sup>273</sup> p53 mutant; at 20 µM, it induced 56% apoptosis in Burkitt lymphoma cell line Ew36 carrying wt p53; and it induced apoptosis in mutant p53 cell lines BL41 (20%) and SW480 (25%). It has no effect in p53 null cell lines HL60 and Saos-21. The p53 activator peptide is a useful tool to study p53.

**Manuals, SDS & CofA**

[View Online »](#)

## Handling & Storage

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

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	Fusion peptide 46
<b>Appearance</b>	Lyophilized solid.
<b>MW</b>	4434.1
<b>Purity</b>	≥97% (HPLC)
<b>Sequence</b>	Gly-Ser-Arg-Ala-His-Ser-Ser-His-Leu-Lys-Ser-Lys-Lys-Gly-Gln-Ser-Thr-Ser-Arg-His-Lys-Lys-Trp-Lys-Met-Arg-Arg-Asn-Gln-Phe-Trp-Val-Lys-Val-Gln-Arg-Gly
<b>Solubility</b>	Soluble in Water.

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)