

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 His²⁷³ 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.

Ordering Information

[Order Online »](#)

BML-P603-0500	0.5mg
----------------------	-------

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

Alternative Name Fusion peptide 46

Appearance Lyophilized solid.

MW 4434.1

Purity ≥97% (HPLC)

Sequence 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

Solubility Soluble in Water.



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)