

Int-H1-S6A F8 c-myc inhibitor

c-Myc inhibitor

H1 DNA-binding region of c-Myc containing Ser to Ala, and Phe to Ala substitutions (underlined) to confer an increase in its potency to inhibit c-Myc. The N-terminus of this peptide is the Int peptide sequence derived from the third Antennapedia homeodomain, to confer cell permeability. Control for this peptide is H1-S6A,F8A c-Myc inhibitor peptide (Prod. No. BML-P606). Inhibited cloning efficiency of MCF-7 human breast cancer cells by 90% at 10 μ M (IC_{50} =5.9 μ M). In MCF-7 cells, it inhibited cell growth and induced apoptosis. In addition, at 10 μ M it strongly inhibited transcription of the c-Myc regulated genes ODC and p53. The non-cell-permeable control for this peptide is H1-S6A, F8A c-Myc inhibitor peptide which lacks the N-terminal internalization sequence.

Citations: 3

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-P605-0500	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

Appearance White to off-white powder.

Formulation Lyophilized.

MW 3873.6

Purity ≥95% (HPLC)

Sequence Arg-Gln-Ile-Lys-Ile-Trp-Phe-Gln-Asn-Arg-Arg-Met-Lys-Trp-Lys-Lys-Asn-Glu-Leu-Lys-Arg-Ala-Phe-Ala-Ala-Leu-Arg-Asp-Gln-Ile

Solubility Soluble in water.

Source Synthetic.



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)