

# 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  $\mu$ M ( $IC_{50}$ =5.9  $\mu$ M). In MCF-7 cells, it inhibited cell growth and induced apoptosis. In addition, at 10  $\mu$ 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.

**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)