

PRODUCT DATA SHEET



PEPTIDE NAME: Int-H1-S6A,F8A c-Myc inhibitor

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*

SEQUENCE DESCRIPTION: 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^{1,2}. The N-term of this peptide is the Int peptide sequence derived from the third antennapedia homeodomain (italics)³, to confer cell-permeability. Control for this peptide is H1-S6A,F8A c-Myc inhibitor peptide (BIOMOL cat.# P-606), which lacks the N-term internalization sequence.

CATALOG NO.: P-605

LOT NO.: P4876

MOLECULAR WEIGHT: 3873.6

PURITY: >95% by HPLC analysis.

AMINO ACID ANALYSIS AND IDENTITY: Confirms.

PEPTIDE CONTENT: 63%

SUPPLIED AS: Trifluoroacetate salt; 0.5 mg net peptide/vial.

PHYSICAL APPEARANCE: Lyophilized solid.

SOLUBILITY: H₂O

STORAGE: -20°C

USAGE: Inhibited cloning efficiency of MCF-7 human breast cancer cells by >90% at 10 μ M (IC₅₀=5.9 μ M)¹. In MCF-7 cells, it inhibited cell growth and induced apoptosis, and strongly inhibited transcription of the c-Myc-regulated genes ODC and p53 in cells treated for 48 hr with 10 μ M Int-H1-S6A,F8A peptide¹.

OTHER: See attached Guide to Handling and Storing Peptides.

LITERATURE REFERENCES:

1. L Giorello *et al.* *Cancer Res.* 1998 **58** 3654
2. L.J. Draeger *et al.* *J. Biol. Chem.* 1994 **269** 1785
3. D. Derossi *et al.* *J. Biol. Chem.* 1994 **269** 10444