

Int-H1-S6A F8 control peptide

Negative control for Int-H1-S6A F8

H1 DNA-binding region of c-Myc containing Ser to Ala and Phe to Ala substitutions (underlined) to confer an increase in its ability to inhibit c-Myc. It inhibits the CM1 consensus sequence binding of c-Myc-92 (C-terminus of c-Myc containing the bHLH region) with $IC_{50} = 1 \mu M$. It is a non-cell-permeable control for Int-H1-S6A,F8A.

Ordering Information

[Order Online »](#)

BML-P606-0500	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

MW 1644.9

Purity ≥95%

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

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)