

CDK5 substrate (biotinylated)

This peptide, Biotin-Ahx-PKTPKKAKKL, is derived from histone H1 and is a substrate for CDK5. The biotin allows the peptide to be used in kinase assays with streptavidine-bound membranes.

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

| | |
|---------------|-----|
| BML-P264-0001 | 1mg |
|---------------|-----|

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 1475.5

Purity ≥95% (HPLC)

Sequence Biotin-Ahx-Pro-Lys-Thr-Pro-Lys-Lys-Ala-Lys-Lys-Leu-OH

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)