

PDKtide (biotinylated)

Protein kinase substrate

This peptide, Biotin-Ahx-KTFCGTPEYLAPEVRREPRILSEEEQEMFRDFDYIADWC, is derived from AKT1 and PKN2/PRK2. It is a substrate for PDK1, CDK9/Cyclin T1 and JAK2. The biotin allows peptide to be used in kinase assays with streptavidin-bound membranes.

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	5109.6
Purity	≥95%
Sequence	Biotin-Ahx-Lys-Thr-Phe-Cys-Gly-Thr-Pro-Glu-Tyr-Leu-Ala-Pro-Glu-Val-Arg-Arg-Glu-Pro-Arg-Ile-Leu-Ser-Glu-Glu-Glu-Gln-Glu-Mez-Phe-Arg-Asp-Phe-Asp-Tyr-Ile-Ala-Asp-Trp-Cys

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com