

Chk 1 & 2 substrate

This peptide sequence, KKKVSRSGLYRSPSPENLNRPR, corresponds to amino acids 206-225 of the phosphatase CDC25C, which is a phosphorylation target of Chk1 on Ser²¹⁶. The addition of three N-terminal lysines allows the peptide to be used in phosphocellulose kinase assays. Both Chk1 and Chk2 will phosphorylate this peptide.

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

Alternative Name	CDC25C (206-225)
MW	2702.0
Sequence	Lys-Lys-Lys-Val-Ser-Arg-Ser-Gly-Leu-Tyr-Arg-Ser-Pro-Ser-Met-Pro-Glu-Asn-Leu-Asn-Arg-Pro-Arg

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