

CSH103 substrate

CDK substrate

This sequence, ADAQHATPPKKRKVED, is SV40 L Tag (118-138) with Ser to Ala¹²⁰ and Ala¹²³ substitutions to increase specificity. Substrate is specific for p34^{cdc2} (CDK1) and is a poor substrate for CKI, II, PKA, GSK3, PKC, CaMKII, and phosphorylase b kinase. Can be used in cell lysates using TCA/phosphocellulose assay. $K_m = 74 \mu\text{M}$ for active complex with cyclin B.

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-P104-0500	500µg
---------------	-------

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

CAS 135546-44-0

MW 2406.7

Sequence Ala-Asp-Ala-Gln-His-Ala-Thr-Pro-Pro-Lys-Lys-Lys-Arg-Lys-Val-Glu-Asp-Pro-Lys-Asp-Phe

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)