

Lyn & Syk kinase substrate

This peptide, EDPIYEFLPAKKK, was identified through phage display as a high affinity substrate for both Lyn kinase ($K_m=24\text{ }\mu\text{M}$) and Syk kinase ($K_m=17\text{ }\mu\text{M}$). The three lysines on the C-terminus allow this peptide to be used in phosphocellulose kinase assays.

Ordering Information

[Order Online »](#)

BML-P217-0001	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

MW	1577.8
Purity	≥95%
Sequence	Glu-Asp-Pro-Ile-Tyr-Glu-Phe-Leu-Pro-Ala-Lys-Lys-Lys

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