

# Fyn kinase substrate

This peptide, EFGTYGTLKSKK, is derived from a peptide (pF4) discovered, through phage display, to be a very good substrate for Fyn kinase. Its  $K_m$  is estimated to be 70  $\mu\text{M}$ . The addition of three lysines on the C-terminus allows this peptide to be used in phosphocellulose kinase assays.

Citations: 3

[View Online »](#)

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	1358.5
Purity	≥95%
Sequence	Glu-Phe-Gly-Thr-Tyr-Gly-Thr-Leu-Ser-Lys-Lys-Lys

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)