

# Shp 1 & 2 substrate

This peptide sequence, EFpYAEVGRSPDPAK, was selected by phage display analysis for optimal Shp-1 and Shp-2 phosphatase activities.

Citations: 1

[View Online »](#)

## Ordering Information

[Order Online »](#)

BML-P157-0001	1mg
---------------	-----

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at -20°C for up to 1 year. Store solutions at -20°C for up to 3 months.
Long Term Storage	-20°C
Shipping	Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

Appearance	White to off-white solid.
Formulation	Lyophilized.
MW	1742.0
Purity	≥95% (HPLC)
Sequence	H-Glu-Phe- <b>pTyr</b> -Ala-Glu-Val-Gly-Arg-Ser-Pro-Pro-Asp-Pro-Ala-Lys
Solubility	Soluble in water
Source	Synthetic.



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)