

# HER2 substrate (biotinylated)

This peptide, Biotin-Ahx-GGMEDIYFEFMGGKKK, identified by screening a degenerate peptide library, is a substrate for several kinases, including HER-2/c-neu/erb-B-2, Tie-2, cFMS, and cKit. The biotin allows peptide to be used in kinase assays with streptavidin-bound membranes.

## Ordering Information

[Order Online »](#)

<b>BML-P243-0001</b>	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

Alternative Name	erbB-2 substrate (biotinylated)
Appearance	White to off-white solid.
Formulation	Lyophilized.
Identity	Determined by MS.
MW	2174.5
Purity	≥95% (HPLC)
Sequence	Biotin-Ahx-Gly-Gly-Met-Glu-Asp-Ile-Tyr-Phe-Glu-Phe-Met-Gly-Gly-Lys-Lys-Lys
Solubility	Soluble in Water.
Source	Synthetic.

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)