

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 »](#)

BML-P243-0001	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

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.



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)