

HER2 substrate

This peptide, GGMEIYFEFMGGKKK, 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.

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

Alternative Name	erbB-2 substrate
MW	1206.33
Purity	≥95%
Sequence	Glu-Gly-Met-Glu-Asp-Ile-Tyr-Phe-Glu-Phe-Met-Gly-Gly-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